

Agilent InfinityLab LC Series 1220 Infinity II LC System

User Manual



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2016

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Manual Part Number

G4280-90020 Rev. B

Edition

09/2016

Printed in Germany

Agilent Technologies
Hewlett-Packard-Strasse 8
76337 Waldbronn

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as "Commercial computer software" as defined in DFAR 252.227-7014 (June 1995), or as a "commercial item" as defined in FAR 2.101(a) or as "Restricted computer software" as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies' standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will

receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

Safety Notices

CAUTION

A CAUTION notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a CAUTION notice until the indicated conditions are fully understood and met.

WARNING

A WARNING notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a WARNING notice until the indicated conditions are fully understood and met.

In This Guide...

This manual covers the Agilent 1220 Infinity II LC System configurations:

- G4286B
- G4288B/C
- G4290B/C
- G4294B

1 Introduction

This chapter provides an overview of the Agilent 1220 Infinity II LC available configurations.

2 Site Requirements and Specifications

This chapter provides information on environmental requirements, physical and performance specifications.

3 Installation

This chapter provides an overview on shipment content and installation.

4 LAN Configuration

This chapter provides information on connecting the instrument to the Agilent ChemStation PC.

5 Solvent Delivery System Description

This chapter provides an overview on the operational principles of the solvent delivery system (pump and optional degasser).

6 Injection System Description

This chapter provides an overview of the operational principles of the injection systems: manual injector and autosampler.

[In This Guide...](#)

7 Column Oven Description

This chapter provides an overview of the operational principles of the column oven.

8 Detector Description

This chapter provides an overview of the operational principles of the detector.

9 Test Functions and Calibration

This chapter describes the tests, calibrations and tools that are available with the Instrument Utilities software or the Lab Advisor (Service and Diagnostics Section).

10 Error Information

This chapter provides information on the error messages that might be displayed, and gives the possible causes and suggestions on their solutions.

11 Maintenance and Repair

This chapter provides general information on maintenance and repair of the instrument.

12 Parts for Maintenance and Repair

This chapter provides information on parts for maintenance and repair.

13 Upgrading the 1220 Infinity II LC

This chapter provides information for upgrading the LC system.

14 Identifying Cables

This chapter provides information on cables used with the Agilent 1200 Infinity Series modules.

15 Appendix

This chapter provides addition information on safety, legal and web.

Contents

1 Introduction 9

- Agilent 1220 Infinity II LC Configurations 10
- Agilent 1220 Infinity II LC VL Configurations 11
- Early Maintenance Feedback 12

2 Site Requirements and Specifications 15

- Site Requirements 16
- Physical Specifications 19
- Performance Specifications 20

3 Installation 29

- Unpacking Your System 30
- Installing the Hardware 35
- Connecting and Configuring the Instrument to the Chromatographic Data System 44
- Connecting the Agilent 1220 Infinity II LC to the PC 45
- The Instrument Utility / Lab Advisor Software 47
- Configuration of the Instrument After an Upgrade Installation 48
- Priming the System and Performing the 'Installation Check' 49
- Performing a 'Checkout Run' 50
- Remove the Restriction Capillary 51
- Install a Column 53
- Remove a Column 55
- Install the Restriction Capillary 57

4 LAN Configuration 59

- To do first 60
- TCP/IP parameter configuration 62
- Configuration Switches 63
- Initialization mode selection 64
- Dynamic Host Configuration Protocol (DHCP) 66
- Link configuration selection 69

Automatic Configuration with BootP	70
Storing the settings permanently with Bootp	79
Manual Configuration	80

5 Solvent Delivery System Description 85

Overview	86
Degasser	87
Principles of Operation	88
Compressibility Compensation	92
Variable Stroke Volume	95
Using the Pump	96

6 Injection System Description 99

Manual Injector	100
Autosampler	103

7 Column Oven Description 115

Column Oven	116
-------------	-----

8 Detector Description 117

Detector Types	118
Agilent 1220 Infinity II LC Variable Wavelength Detector (VWD)	119
Agilent 1220 Infinity II LC Diode Array Detector (DAD)	120
Match the Flow Cell to the Column	136

9 Test Functions and Calibration 141

1220 Infinity II LC System	143
Solvent Delivery System	145
Autosampler	154
Column Oven	160
Variable Wavelength Detector (VWD)	162
Diode Array Detector (DAD)	171

10 Error Information 193

- What are Error Messages? 196
- General Error Messages 197
- Pump Error Messages 203
- Autosampler Error Messages 216
- General Detector Error Messages 226
- VWD Detector Error Messages 229
- DAD Detector Error Messages 235

11 Maintenance and Repair 241

- PM Scope of Work and Checklist 243
- Cautions and Warnings 244
- Solvent Delivery System 246
- Manual Injector 267
- Autosampler 271
- Variable Wavelength Detector (VWD) 290
- Diode Array Detector (DAD) 302
- Algae Growth in HPLC Systems 323
- Replacing the Module's Firmware 325

12 Parts for Maintenance and Repair 327

- 1220 Infinity II LC System 328
- Solvent Delivery System 331
- Injection System 340
- Column Oven 347
- Detector 348

13 Upgrading the 1220 Infinity II LC 355

- Oven Upgrade 356
- Install the (optional) External Tray 359
- Install the Active Inlet Valve 363

Contents

14 Identifying Cables 367

Cable Overview	368
Analog Cables	370
Remote Cables	372
BCD Cables	375
CAN/LAN Cables	377
Agilent 1200 module to PC	378

15 Appendix 379

General Safety Information	380
Solvent Information	383
Radio Interference	385
UV Radiation	386
Sound Emission	387
Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC)	388
Declaration of Conformity for HOX2 Filter	389
Agilent Technologies on Internet	390

1 Introduction

Agilent 1220 Infinity II LC Configurations	10
Agilent 1220 Infinity II LC VL Configurations	11
Early Maintenance Feedback	12
EMF counters for the pump	12
EMF counters for the autosampler	12
EMF counters for the variable wavelength detector	13
EMF counters for the diode array detector	14

This chapter provides an overview of the Agilent 1220 Infinity II LC available configurations.



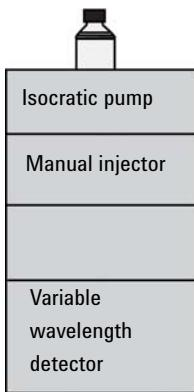
1 Introduction

Agilent 1220 Infinity II LC Configurations

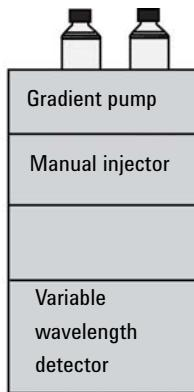
Agilent 1220 Infinity II LC Configurations

Available configurations of Agilent 1220 Infinity II LC

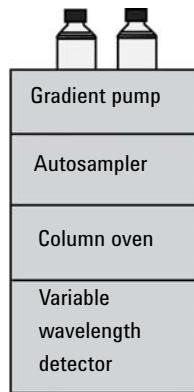
The Agilent 1220 Infinity II LC is available in four different configurations. Possible components include isocratic pump, dual-channel gradient pump (with degasser), manual injector, autosampler, column oven and detector. Each configuration comes with at least one pump, one injection system and one detector and includes Agilent Instrument Utilities Software.



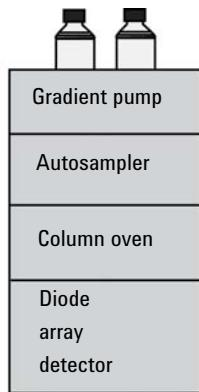
G4286B



G4288B



G4290B



G4294B

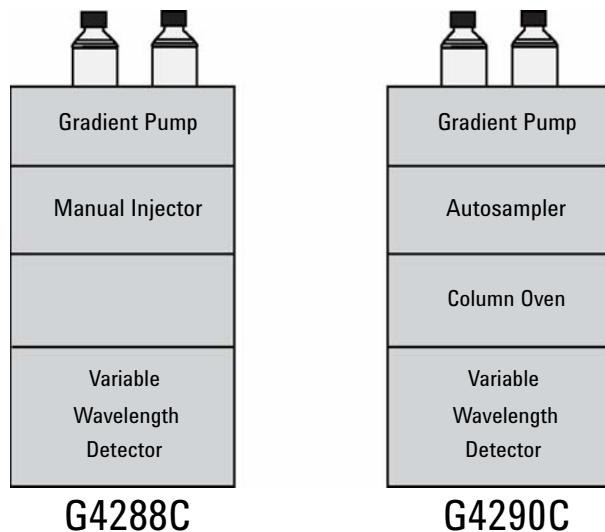
A Solvent Selection Valve (SSV) Upgrade Kit (G4280-68708) is available.

A 1220 Infinity II AIV Upgrade Kit (G5699B) is available to handle applications with high buffer concentration.

Agilent 1220 Infinity II LC VL Configurations

Available configurations of Agilent 1220 Infinity II LC VL

The Agilent 1220 Infinity II LC VL is available in two different configurations. Possible components include isocratic pump, dual-channel gradient pump (with degasser), manual injector, autosampler, column oven and detector. Each configuration comes with at least one pump, one injection system and one detector and includes Agilent Instrument Utilities Software.



A Solvent Selection Valve (SSV) Upgrade Kit (G4280-68708) is available.

A 1220 Infinity II AIV Upgrade Kit (G5699B) is available to handle applications with high buffer concentration.

1 Introduction

Early Maintenance Feedback

Early Maintenance Feedback

EMF counters for the pump

The user-settable EMF limits for the EMF counters enable the early maintenance feedback to be adapted to specific user requirements. The wear of pump components is dependent on the analytical conditions. Therefore, the definition of the maximum limits needs to be determined based on the specific operating conditions of the instrument.

The Agilent 1220 Infinity II LC pump provides a series of EMF counters for the pump head. Each counter increments with pump use, and can be assigned a maximum limit that provides visual feedback in the user interface when the limit is exceeded. Each counter can be reset to zero after maintenance has been done. The pump provides the following EMF counters:

Pump liquimeter

The pump liquimeter displays the total volume of solvent pumped by the pump head since the last reset of the counters. The pump liquimeter can be assigned an EMF (maximum) limit. When the limit is exceeded, the EMF flag in the user interface is displayed.

Seal wear counters

The seal wear counters display a value derived from pressure and flow (both contribute to seal wear). The values increment with pump usage until the counters are reset after seal maintenance. Both seal wear counters can be assigned an EMF (maximum) limit. When the limit is exceeded, the EMF flag in the user interface is displayed.

EMF counters for the autosampler

The user-settable EMF limits for the EMF counters enable the early maintenance feedback to be adapted to specific user requirements. The wear of autosampler components is dependent on the analytical conditions.

Therefore, the definition of the maximum limits need to be determined based on the specific operating conditions of the instrument.

The autosampler provides two EMF counters. Each counter increments with autosampler use, and can be assigned a maximum limit which provides visual feedback in the user interface when the limit is exceeded. Each counter can be reset to zero after maintenance has been done. The autosampler provides the following EMF counters:

Injection valve counter

This counter displays the total number of switches of the injection valve since the last reset of the counter.

Needle movements counter

This counter displays the total number of movements of the needle into the seat since the last reset of the counter.

EMF counters for the variable wavelength detector

The user-settable EMF limits for the EMF counter enables the early maintenance feedback to be adapted to specific user requirements. The useful lamp burn time is dependent on the requirements for the analysis (high or low sensitivity analysis, wavelength, and so on). Therefore, the definition of the maximum limits need to be determined based on the specific operating conditions of the instrument.

The detector module provides a EMF counter for the lamp. The counter increments with lamp use, and can be assigned a maximum limit which provides visual feedback in the user interface when the limit is exceeded. The counter can be reset to zero after the lamp is exchanged. The detector provides the following EMF counters:

Deuterium lamp on-time

This counter shows the total burn time of the deuterium lamp in hours.

1 Introduction

Early Maintenance Feedback

EMF counters for the diode array detector

Using the EMF Counters

The user-settable **EMF** limits for the **EMF Counters** enable the early maintenance feedback to be adapted to specific user requirements. The useful maintenance cycle is dependent on the requirements for use. Therefore, the definition of the maximum limits need to be determined based on the specific operating conditions of the instrument.

Setting the EMF Limits

The setting of the **EMF** limits must be optimized over one or two maintenance cycles. Initially the default **EMF** limits should be set. When instrument performance indicates maintenance is necessary, take note of the values displayed by the **EMF counters**. Enter these values (or values slightly less than the displayed values) as **EMF** limits, and then reset the **EMF counters** to zero. The next time the **EMF counters** exceed the new **EMF** limits, the **EMF** flag will be displayed, providing a reminder that maintenance needs to be scheduled.

2

Site Requirements and Specifications

Site Requirements	16
Power Considerations	16
Power Cord	17
Bench Space	17
Environment	18
Physical Specifications	19
Performance Specifications	20
Specification Conditions	27

This chapter provides information on environmental requirements, physical and performance specifications.



2 Site Requirements and Specifications

Site Requirements

Site Requirements

A suitable environment is important to ensure optimal performance of the instrument.

Power Considerations

The Agilent 1220 Infinity II LC power supply has wide-ranging capabilities. Consequently, there is no voltage selector at the instrument.

WARNING

Instrument is partially energized when switched off

The power supply still uses some power even when the power switch on the front panel is turned OFF. Repair work at the detector can lead to personal injuries, e. g. shock hazard, when the detector cover is opened and the instrument is connected to power.

- To disconnect the detector from the power line, unplug the power cord.
-

WARNING

Incorrect line voltage to the instrument

Shock hazard or damage to your instrumentation can result if the devices are connected to a line voltage higher than specified.

- Connect your instrument only to the specified line voltage.
-

CAUTION

In case of an emergency, it must be possible to disconnect the instrument from the power line at any time.

Make sure that there is easy access to the power cable of the instrument so that the instrument can quickly and easily be disconnected from the line voltage.

- Provide sufficient space next to the power socket of the instrument to allow the cable to be unplugged.
-

Power Cord

Different power cords are offered as options with the system. The female ends of all power cords are identical. The female end plugs into the power-input socket at the rear left side of the instrument. The male end of each power cord is different and designed to match the wall socket of a particular country or region.

WARNING

Absence of ground connection or use of unspecified power cord

The absence of ground connection or the use of unspecified power cord can lead to electric shock or short circuit.

- Never operate your instrument from a power outlet that has no ground connection.
 - Never use a power cord other than the Agilent Technologies power cord designed for your region.
-

WARNING

Use of cables not supplied by Agilent

Using cables that have not been supplied by Agilent Technologies can lead to damage of the electronic components or personal injury.

- Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.
-

Bench Space

The dimensions and weight of the Agilent 1220 Infinity II LC allow it to be placed on almost any desk or laboratory bench. It needs an additional 2.5 cm (1.0 in) of space on either side and approximately 8 cm (3.1 in) at the rear for air circulation and electric connections.

Make sure that the bench intended to carry the Agilent 1220 Infinity II LC is designed to bear the weight of the instrument.

The Agilent 1220 Infinity II LC should be operated upright.

2 Site Requirements and Specifications

Site Requirements

Environment

Your Agilent 1220 Infinity II LC will work within specifications at ambient temperatures and relative humidity as described in the following sections.

ASTM drift tests require a temperature change below 2 °C/hour (3.6 °F/hour) measured over one hour period. Our published drift specification is based on these conditions. Larger ambient temperature changes will result in larger drift.

Better drift performance depends on better control of the temperature fluctuations. To realize the highest performance, minimize the frequency and the amplitude of the temperature changes to below 1 °C/hour (1.8 °F/hour). Turbulences around one minute or less can be ignored.

CAUTION

Condensation within the module

Condensation can damage the system electronics.

- Do not store, ship or use your module under conditions where temperature fluctuations could cause condensation within the module.
 - If your module was shipped in cold weather, leave it in its box and allow it to warm slowly to room temperature to avoid condensation.
-

Physical Specifications

Table 1 Physical Specifications

Type	Specification	Comments
Weight	30 kg (66 lbs) G4294B: 43 kg (94 lbs)	
Dimensions (height × width × depth)	640 × 370 × 420 mm (25.2 × 14.6 × 16.5 inches) G4294B: 640×370×485 mm (25.2×14.6×19.1 inches)	
Line voltage	100 – 240 V~, ± 10 %	Wide-ranging capability
Line frequency	50 or 60 Hz, ± 5 %	
Power consumption	240 VA / 210 W / 717 BTU	Maximum
Ambient operating temperature	4–55 °C (39–131 °F)	
Ambient non-operating temperature	-40 – 70 °C (-40 – 158 °F)	
Humidity	< 95 % r.h. at 40 °C (104 °F)	Non-condensing
Operating altitude	Up to 2000 m (6562 ft)	
Non-operating altitude	Up to 4600 m (15092 ft)	For storing the module
Safety standards: IEC, CSA, UL	Installation category II, Pollution degree 2	For indoor use only.

2 Site Requirements and Specifications

Performance Specifications

Performance Specifications

Performance Specifications Agilent 1220 Infinity II LC

Table 2 Performance Specifications Agilent 1220 Infinity II LC

Type	Specification
Safety features	Extensive diagnostics, error detection and display, leak detection, safe leak handling, leak output signal for shutdown of pumping system. Low voltages in major maintenance areas.
Control and data evaluation	Agilent EZChrom Compact, Agilent ChemStation, Agilent Instrument Utilities, Agilent Lab Advisor
Communications	Controller-area network (CAN), RS-232C, APG Remote: ready, start, stop and shut-down signals, LAN
GLP features	Early maintenance feedback (EMF), electronic records of maintenance and errors

Performance Specifications Agilent 1220 Infinity II LC Pump

Table 3 Performance Specifications Agilent 1220 Infinity II LC Pump

Type	Specification	Comments
Hydraulic system	Dual plunger in series pump with proprietary servo-controlled variable stroke drive, floating plungers and passive inlet valve	
Flow range	Settable: 0.001 – 10 mL/min Recommended: 0.2 – 10.0 mL/min	Set points in 0.001 mL/min increments
Flow precision	$\leq 0.07\%$ RSD, or ≤ 0.02 min SD whichever is greater	Based on retention time at constant temperature
Flow accuracy	$\pm 1\%$ or 10 μ L/min whatever is greater	Pumping degassed H ₂ O at 10 MPa (100 bar)
Pressure operating range	Up to 60 MPa (600 bar, 8702 psi) up to 5 mL/min Up to 20 MPa (200 bar, 2901 psi) up to 10 mL/min	
Pressure pulsation	< 2 % amplitude (typically < 1.3 %), or < 0.3 MPa (3 bar, 44 psi), whichever is greater	
Compressibility compensation	User-selectable, based on mobile phase compressibility	
Recommended pH	1.0 – 12.5	Solvents with pH < 2.3 should not contain acids which attack stainless steel
Gradient formation (optional)	Low pressure binary mixing/gradient capability using proprietary high-speed proportioning valve	
Delay volume	600 – 900 μ L, dependent on back pressure	Measured with water at 1 mL/min (water/water with tracer)
Composition range	Settable 0 – 100 % Recommended 5 – 95 %	In 0.1 % increments
Composition precision	< 0.2 % RSD or < 0.04 min SD, whichever is greater	At 1 mL/min; based on retention time at constant temperature

2 Site Requirements and Specifications

Performance Specifications

Performance Specifications Agilent 1220 Infinity II LC Pump VL

Table 4 Performance Specifications Agilent 1220 Infinity II LC Pump VL

Type	Specification	Comments
Hydraulic system	Dual plunger in series pump with proprietary servo-controlled variable stroke drive, floating plungers and passive inlet valve	
Flow range	Settable 0.001 – 10 mL/min Recommended: 0.2 – 10.0 mL/min	Set points in 0.001 mL/min increments
Flow precision	$\leq 0.07\%$ RSD, or ≤ 0.02 min SD whichever is greater	Based on retention time at constant temperature
Flow accuracy	$\pm 1\%$ or 10 μ L/min whatever is greater	Pumping degassed H ₂ O at 10 MPa (100 bar)
Pressure operating range	Up to 40 MPa (400 bar, 5801 psi) up to 5 mL/min Up to 20 MPa (200 bar, 2901 psi) up to 10 mL/min	
Pressure pulsation	< 2 % amplitude (typically < 1.3 %), or < 0.3 MPa (3 bar, 44 psi), whichever is greater	
Compressibility compensation	User-selectable, based on mobile phase compressibility	
Recommended pH range	1.0 – 12.5	Solvents with pH < 2.3 should not contain acids which attack stainless steel
Gradient formation	Low pressure binary mixing/gradient capability using proprietary high-speed proportioning valve	
Delay volume	600 – 900 μ L, dependent on back pressure	Measured with water at 1 mL/min (water/water with tracer)
Composition range	Settable: 0 – 100 % Recommended: 5 – 95 %,	In 0.1 % increments
Composition precision	< 0.2 % RSD or < 0.04 min SD, whichever is greater	At 1 mL/min; based on retention time at constant temperature

Performance Specifications Agilent 1220 Infinity II LC Autosampler

Table 5 Performance Specifications Agilent 1220 Infinity II LC Autosampler

Type	Specification	Comments
Injection range	0.1 – 100 µL in 0.1 µL increments	Up to 1500 µL with multiple draw (hardware modification required)
Precision	<1 % RSD of peak areas from 1 µL to 5 µL <0.25 % RSD of peak areas from 5 µL to 100 µL	Measured with caffeine
Replicate injections	1 – 99 from one vial	
Pressure range	0 – 60 MPa (0 – 600 bar, 0 – 8702 psi)	
Sample viscosity range	0.2 – 50 cp	
Sample capacity	100 × 2 mL vials in 1 tray 40 × 2 mL vials in ½ tray 15 × 6 mL vials in ½ tray (Agilent vials only)	
Carry over	Typically < 0.1 %, < 0.05 % with external needle cleaning	
Injection cycle time	Typically 50 s depending on draw speed and injection volume	
Minimum sample volume	1 µL from 5 µL sample in 100 µL microvial, or 1 µL from 10 µL sample in 300 µL microvial	

2 Site Requirements and Specifications

Performance Specifications

Performance Specifications Agilent 1220 Infinity II LC Column Oven

Table 6 Performance Specifications Agilent 1220 Infinity II LC Column Oven

Type	Specification	Comments
Temperature range	5 °C above ambient to 80 °C 10 °C above ambient to 80 °C (G4294B with DAD)	
Column capacity	1 column up to 25 cm	
Temperature stability	± 0.15 °C	Constant composition and flow rate
Temperature accuracy	± 0.8 °C	
Internal volume	6 µL	

Performance Specifications Agilent 1220 Infinity II LC VWD

Table 7 Performance Specifications Agilent 1220 Infinity II LC VWD

Type	Specification	Comments
Detection type	Double-beam photometer	
Light source	Deuterium lamp	
Maximum data rate	80 Hz	
Short term noise (ASTM)	<± 0.25·10 ⁻⁵ AU at 230 nm	See <i>Specification Conditions</i> below
Drift	< 1·10 ⁻⁴ AU/h at 230 nm	See <i>Specification Conditions</i> below
Linearity	> 2 AU (5 %) upper limit	See <i>Specification Conditions</i> below
Wavelength range	190 – 600 nm	
Wavelength accuracy	± 1 nm	Self-calibration with deuterium lines, verification with holmium oxide filter
Slit width	6.5 nm typical over whole wavelength range	
Flow cells	Standard: 14 µL volume, 10 mm cell path length and 40 bar (580 psi) pressure maximum High pressure: 14 µL volume, 10 mm cell path length and 400 bar (5800 psi) pressure maximum Micro: 1 µL volume, 5 mm cell path length and 40 bar (580 psi) pressure maximum Semi-micro: 5 µL volume, 6 mm cell path length and 40 bar (580 psi) pressure maximum	Can be repaired on component level

2 Site Requirements and Specifications

Performance Specifications

Performance Specifications Agilent 1220 Infinity II LC DAD

Table 8 Performance Specifications Agilent 1220 Infinity II LC DAD

Type	Specification	Comments
Detection type	1024-element diode array	
Light source	Deuterium and tungsten lamps	The UV-lamp is equipped with RFID tag that holds lamp typical information.
Number of signals	8	
Maximum sampling rate	80 Hz	
Short term noise (ASTM) Single and Multi-Wavelength	< ± 0.7·10 ⁻⁵ AU at 254 nm and 750 nm	see "Specification Conditions" below
Drift	< 0.9·10 ⁻³ AU/h at 254 nm	see "Specification Conditions" below
Linearity	> 2 AU (5 %)	see "Specification Conditions" below
Wavelength range	190 – 950 nm	
Wavelength accuracy	± 1 nm	Self-calibration with deuterium lines, verification with holmium oxide filter
Slit width	1, 2, 4 , 8, 16 nm	Programmable slit
Diode width	~ 1 nm	
Flow cell	Standard: 13 µL volume, 10 mm cell path length and 120 bar (1740 psi) pressure maximum	The flow cell is equipped with RFID tags that hold cell typical information. pH range 1.0 – 9.5
Time programmable	Wavelength, polarity, peak width, lamp bandwidth, autobalance, wavelength range, threshold, spectra storage mode	

Specification Conditions

ASTM: “Standard Practice for Variable Wavelength Photometric Detectors Used in Liquid Chromatography”.

VWD

Drift and noise: Standard flow cell, path length 10 mm, flow 1 mL/min LC-grade methanol, TC 2 s.

$$RT = 2.2 * TC$$

Linearity: Standard flow cell, path length 10 mm, measured with caffeine at 273 nm.

DAD

Drift and noise: Standard flow cell, path length 10 mm, flow 1 mL/min LC-grade methanol, TC 2 s. Measurements were performed at 254 nm or 750 nm with bandwidths of 4 nm (reference wavelengths 360 nm and 850 nm with bandwidths of 100 nm). Slit width 4 nm.

$$RT = 2.2 * TC$$

Linearity: Standard flow cell, path length 10 mm, measured with caffeine at 273 nm and a bandwidth of 4 nm.

NOTE

The specifications are based on the standard lamp (G1314-60100 for VWD, G2140-0820 for DAD) and may not be achieved when other lamp types or aged lamps are used.

ASTM drift tests require a temperature change below 2 °C/h (3.6 °F/h) over one hour period. Our published drift specification is based on these conditions. Larger ambient temperature changes will result in larger drift.

Better drift performance depends on better control of the temperature fluctuations. To realize the highest performance, minimize the frequency and the amplitude of the temperature changes to below 1 °C/h (1.8 °F/h). Turbulences around one minute or less can be ignored.

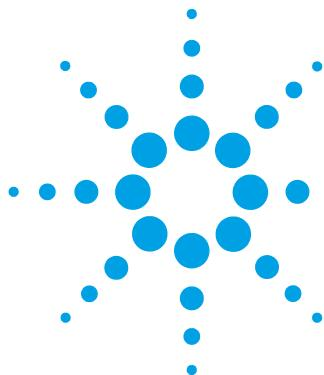
Performance tests should be done with a completely warmed up optical unit (> two hours). ASTM measurements require that the detector should be turned on at least 24 hours before start of testing.

2 Site Requirements and Specifications

Performance Specifications

Time Constant (TC) versus Response Time (RT)

According to ASTM E1657-98 "Standard Practice of Testing Variable-Wavelength Photometric Detectors Used in Liquid Chromatography" the time constant is converted to response time by multiplying by the factor 2.2.



3 Installation

Unpacking Your System	30
Delivery Checklist	30
Installing the Hardware	35
Installation Paths	35
Installing the Agilent 1220 Infinity II LC	37
Identifying the connections of the 1220 Infinity II LC	42
Connecting and Configuring the Instrument to the Chromatographic Data System	44
Connecting the Agilent 1220 Infinity II LC to the PC	45
The Instrument Utility / Lab Advisor Software	47
Configuration of the Instrument After an Upgrade Installation	48
Priming the System and Performing the 'Installation Check'	49
Performing a 'Checkout Run'	50
Remove the Restriction Capillary	51
Install a Column	53
Remove a Column	55
Install the Restriction Capillary	57

This chapter provides an overview on shipment content and installation.

NOTE

To install the Agilent 1220 Infinity II LC System, it is highly recommended to follow the installation instructions step by step.



Agilent Technologies

3 Installation

Unpacking Your System

Unpacking Your System

If the delivery packaging shows signs of external damage, please call your Agilent Technologies sales and service office immediately. Inform your service representative that the Agilent 1220 Infinity II LC may have been damaged during shipment.

CAUTION

Signs of damage

→ Do not attempt to install the Agilent 1220 Infinity II LC.

Delivery Checklist

Delivery Checklist

Ensure all parts and materials have been delivered with the Agilent 1220 Infinity II LC. The delivery checklist is shown below. Please report missing or damaged parts to your local Agilent Technologies sales and service office.

Table 9 Agilent 1220 Infinity II LC Checklist

Description	Quantity
Agilent 1220 Infinity II LC	1
Power cable	1
Flow cell	Installed
User Manual on Documentation Medium (part of the shipment - not module specific)	1 per order
Installation guide	1
Accessory kit (see below)	1

Accessory Kit Contents for G4286B

p/n	Description
G4286-68755	Accessory kit complete
0100-2562	Fitting, onepiece, fingertight
0890-1195	PTFE tubing, 0.052 in i.d.
0890-1711	Flexible tubing (to waste), 3 m
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5062-8535	Waste accessory kit
5188-2758	PTFE/silicone septa 16mm pre-silt 100/pk (delivered quantity is 0.010)
5190-1501	Syringe, 50.0 µL, FN, LC tip
9301-0411	Syringe, Plastic
9301-1337	Syringe adapter
9301-1377	Screw Cap Vial, clear, 6 mL 100/PK (delivered quantity is 0.010)
9301-1379	Screw caps for 6 mL vials 100/PK (delivered quantity is 0.010)
9301-1420	Solvent bottle, transparent
G1311-60003	Bottle-head assembly

3 Installation

Unpacking Your System

Accessory Kit Contents for G4288B/C

#	p/n	Description
1	G4288-68755	Accessory kit complete
1	0100-2562	Fitting, onepiece, fingertight
1	0890-1195	PTFE tubing, 0.052 in i.d.
1	0890-1711	Flexible tubing (to waste), 3 m
1	5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
1	5062-8535	Waste accessory kit
1	5188-2758	PTFE/silicone septa 16mm pre-silt 100/pk (delivered quantity is 0.010)
1	5190-1501	Syringe, 50.0 µL, FN, LC tip
1	9301-0411	Syringe, Plastic
1	9301-1337	Syringe adapter
1	9301-1377	Screw Cap Vial, clear, 6 mL 100/PK (delivered quantity is 0.010)
1	9301-1379	Screw caps for 6 mL vials 100/PK (delivered quantity is 0.010)
1	9301-1420	Solvent bottle, transparent
1	9301-1450	Solvent bottle, amber
2	G1311-60003	Bottle-head assembly

Accessory Kit Contents for G4290B/C, G4294B

p/n	Description
G4290-68755	Accessory kit complete
0100-2562	Fitting, onepiece, fingertight
0890-1195	PTFE tubing, 0.052 in i.d.
0890-1711	Flexible tubing (to waste), 3 m
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5062-8535	Waste accessory kit
9301-0411	Syringe, Plastic
9301-1337	Syringe adapter
9301-1420	Solvent bottle, transparent
9301-1450	Solvent bottle, amber
G1311-60003	Bottle-head assembly

3 Installation

Unpacking Your System

Optional Tool Kit for Agilent 1220 Infinity II LC

p/n	Description
G4296-68715	Tool kit complete
0100-1710	Mounting Tool for Tubing Connections
8710-0510	Wrench open 1/4 — 5/16 inch
8710-1924	Wrench open 14 mm
8720-0025	Wrench, 1/2 inch & 9/16 inch
01018-23702	Insert tool
8710-2392	Hex key 4 mm15 cm long T-handle
8710-2394	Hex key 9/64 inch 15 cm long T-handle
8710-2411	Hex key 3 mm12 cm long
8710-2412	Hex key 2.5 mm, 15 cm long, straight handle
8710-0899	Pozidriv screwdriver

Installing the Hardware

Installation Paths

Standard Installation Path

(No optional hardware upgrade is added to the module during installation)

Installation Path Including the Installation of Upgrade Kits

(Oven upgrade kit /Manual injector to ALS upgrade kit/ Isocratic to gradient upgrade kit)

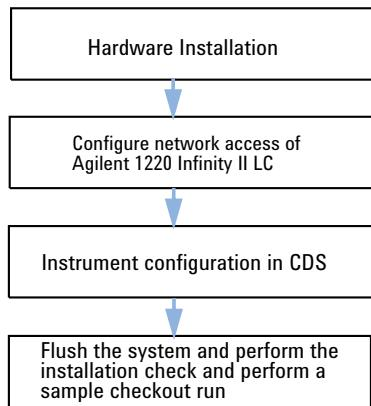
NOTE

The Installation of the gradient system upgrade kit and ALS upgrade kit must be done from Agilent-trained service personnel only.

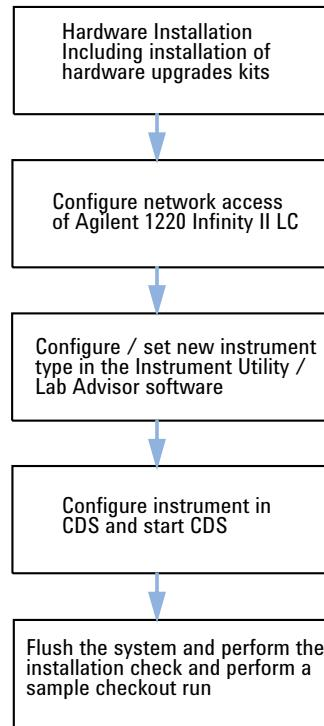
3 Installation

Installing the Hardware

Standard Installation Path



Installation Path Including the Installation of Upgrade Kits



NOTE

The installation of a Solvent Selection Valve (SSV) Option does not require the configuration of a new instrument type. The SSV needs to be configured in the CDS only.

Installing the Agilent 1220 Infinity II LC

- 1 Open the box and compare its content with the delivery checklist for completeness.
- 2 Place the instrument on top of the bench.
- 3 Remove both front covers (top and lower) by pressing the release buttons (on both sides).

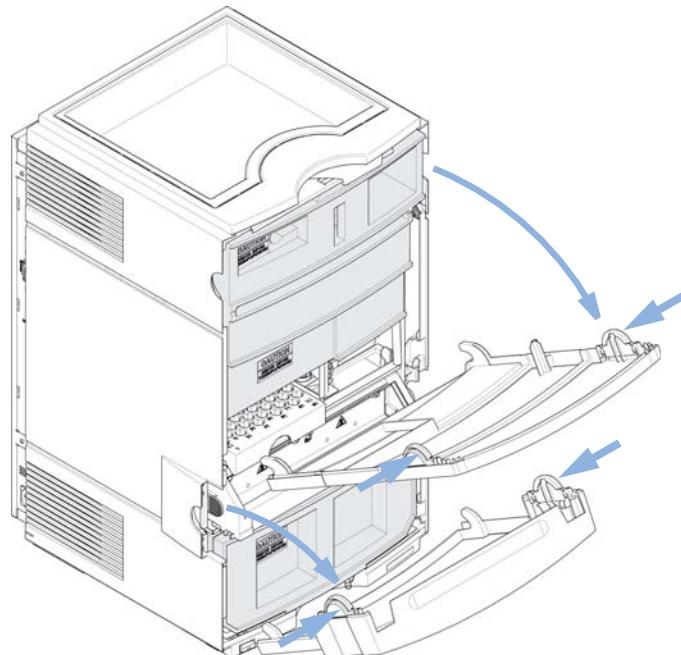


Figure 1 Front Cover Mechanism

3 Installation

Installing the Hardware

- 4 Remove the two transport foams.

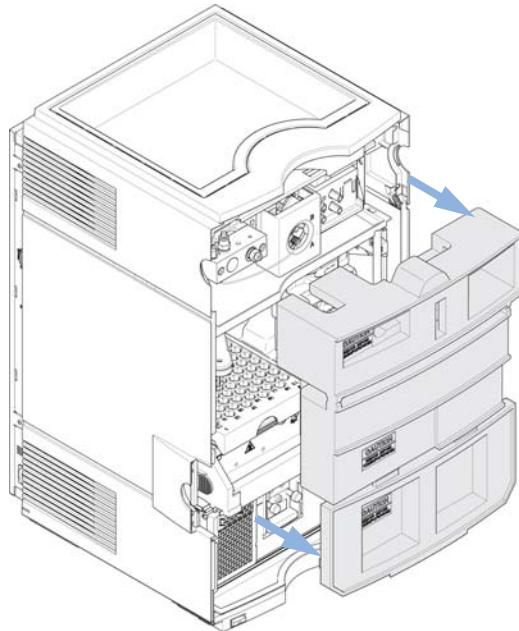


Figure 2 Remove the Transport Foam

[Figure 3](#) on page 39 shows the content of a fully equipped 1220 Infinity II LC system with removed front covers. (Module type shown G4290B)

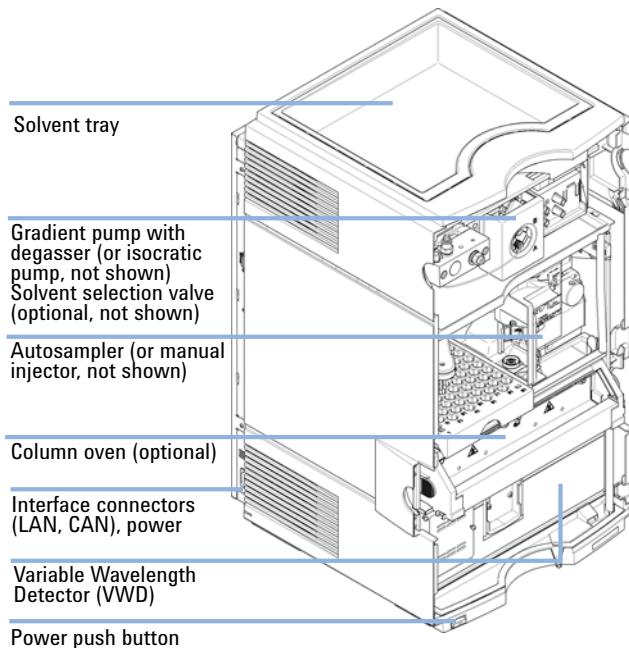


Figure 3 System Overview 1220 Infinity II LC

NOTE

Additional options or upgrade kits should be installed prior to all solvent path installations. Configuring your Agilent 1220 Infinity II LC Module Information can be found in the [“Configuration of the Instrument After an Upgrade Installation”](#) on page 48.

For further information about how to install the options and upgrade kits refer to the user manual.

- 5 Place the Solvent Bottle filled with 0.5 L HPLC grade water in the Solvent Tray.

3 Installation

Installing the Hardware

- 6 Place the Solvent Inlet Filter end of the Bottle Head Assembly in the Solvent Bottle (see picture below).

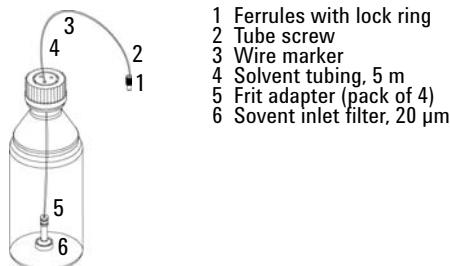


Figure 4 Bottle Head Assembly and Solvent Bottle

- 7 Prime the tubing using the Syringe (9301-044) and Syringe adapter (9301-1337) (part of the Accessory kit) until the tubing is completely filled with water.
- 8 Connect the Bottle Head Assembly connector (see [Figure 4](#) on page 40 Item 1+2) to:
 - the passive inlet valve (isocratic pump) or
 - the degasser inlet Channel A (gradient pump).
- 9 Connect the waste tube with the fitting attached (part of Accessory Kit) to the flow cell outlet and the other end to an appropriate solvent waste container (see [Figure 5](#) on page 41).
- 10 Attach the corrugated waste tube (part of Accessory Kit) to the VWD leak tray outlet adapter and guide it to a proper waste container (see [Figure 5](#) on page 41).
- 11 Connect the waste tube (part of Accessory Kit) to the purge valve outlet adapter and the other end to the waste container.
- 12 Connect the network connection between the instrument and your PC.

NOTE

More details about how to establish a network connection to the instrument can be found in "[Connecting the Agilent 1220 Infinity II LC to the PC](#)" on page 45 or "[LAN Configuration](#)" on page 59

- 13 Verify that the power push button at the front of the module (see [Figure 5](#) on page 41) stands off. Now connect the power cord to the instrument and the power line.

- 14** Before switching on the module check that all transport foams have been de-installed (see [Figure 2](#) on page 38). Then switch on the module via the power push button.

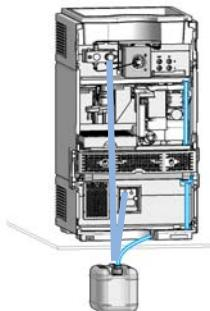


Figure 5 Leak and waste connections (example shows instrument with VWD)

3 Installation

Installing the Hardware

Identifying the connections of the 1220 Infinity II LC

Agilent 1220 Infinity II LC with VWD

Figure 6 on page 42 shows an overview of the possible connections on the 1220 Infinity II LC instrument with VWD.

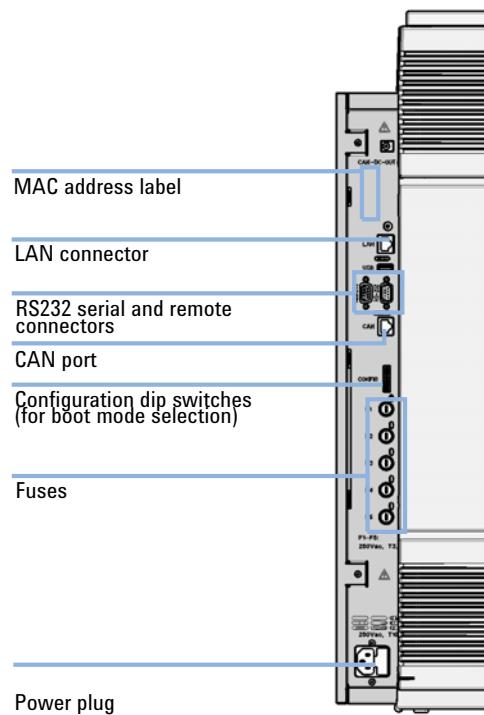


Figure 6 Connections of the Agilent 1220 Infinity II LC with VWD

Agilent 1220 Infinity II LC with DAD

Figure 7 on page 43 shows an overview of the possible connections on the 1220 Infinity II LC instrument with DAD. On the G4294B you must use the configuration dip switches on the DAD mainboard because this will be the communication host in this case. The short CAN cable is the communication connection between the DAD and the other 1220 modules.

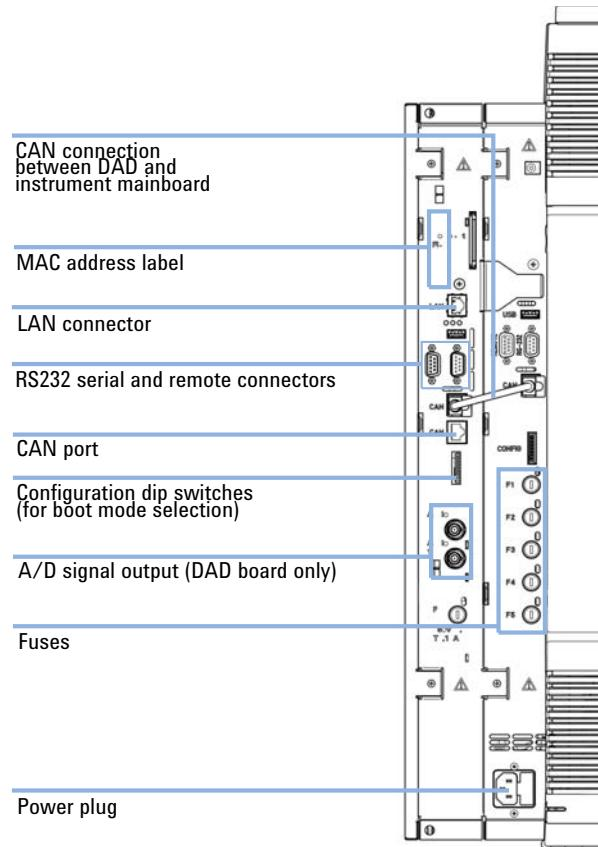


Figure 7 Connections of the Agilent 1220 Infinity II LC with DAD

3 Installation

Connecting and Configuring the Instrument to the Chromatographic Data System

Connecting and Configuring the Instrument to the Chromatographic Data System

The instrument is compatible to the following software for control and data evaluation:

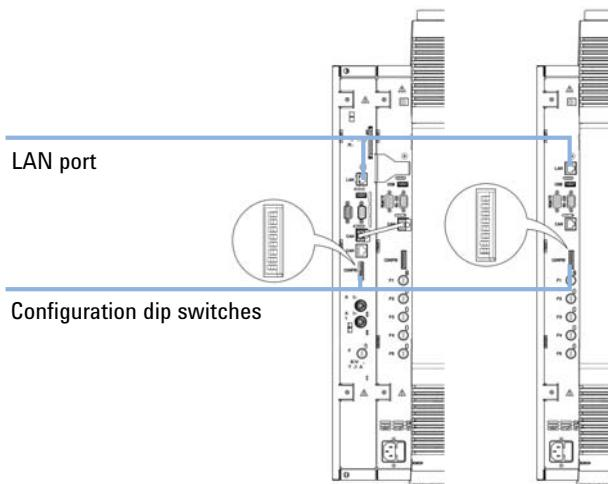
- Agilent ChemStation for LC
 - EZChrom Software
- 1 Install your Chromatographic Data System (CDS). Please refer to the installation documentation that has been delivered with the CDS.
 - 2 Start your CDS.
 - 3 At the instrument configuration screen enter the Instrument name (free of choice) and the Instrument type (*Agilent Compact LC*).
 - 4 For configuring the module choose **Auto Configuration**.

Connecting the Agilent 1220 Infinity II LC to the PC

The instrument is delivered from factory with the default network configuration settings. (Configuration dip switches 7 & 8 set to ON). This enables you to quickly connect it via the crossover patch cable (part of the accessory kit) with your PC.

Factory default IP address:

192.168.254.11



Instrument with DAD Instrument with VWD

Figure 8 Location of the Configuration Dip Switches and LAN Port

NOTE

On the G4294B use the configuration dip switches of the DAD extension board to configure your LAN connection.

3 Installation

Connecting the Agilent 1220 Infinity II LC to the PC

- 1 To connect the instrument with your PC using this default address configure the PC's network settings as followed:

IP: 192.168.254.10

Subnet Mask: 255.255.255.0

Default Gateway: N/A

- 2 Connect the crossover patch cable between the LAN-Port of the Agilent 1220 Infinity II LC ([Figure 8](#) on page 45) and the network connector of your PC.

NOTE

The crossover cable is only for the direct connection between the module and the PC. If you want to connect your Agilent 1220 Infinity II LC via a hub to the network you should contact your local network administrator.

If you want to connect the instrument to a network we strongly recommend consulting your local network administrator to provide you with a valid network address. For further details about LAN configuration of the instrument please refer to the *LAN Configuration* section of the *Installation chapter* in the user manual.

Here you find details about:

- TCP/IP parameter configuration
- Configuration Switches
- Using a fixed IP address
- How to configure an individual IP address

The Instrument Utility / Lab Advisor Software

During the installation process of the instrument this software is used for flushing the system and performing the System Installation check (see “[Priming the System and Performing the ‘Installation Check’](#)” on page 49).

- 1 Install the Instrument Utility or Lab Advisor software according to the Installation Procedure on the software CD.
- 2 Setup your Instrument to the software and **Connect it**.

NOTE

Whenever a hardware upgrade (Isocratic to Gradient, Column Oven, Autosampler upgrade) has been installed, it is necessary to re-configure your instrument in the Instrument Utility / Lab Advisor software. Therefore please follow the steps that are described in “[Configuration of the Instrument After an Upgrade Installation](#)” on page 48 before you proceed.

3 Installation

Configuration of the Instrument After an Upgrade Installation

Configuration of the Instrument After an Upgrade Installation

This step is only necessary if any of the following hardware upgrade kits has been installed to the instrument.

- G4297A - 1220 Infinity Oven kit
- G4298A - 1220 Infinity upgrade manual injector to ALS
- G4299A - 1220 Infinity upgrade isocratic to gradient pump

1 Connect the Instrument to the Instrument Utility / Lab Advisor software.

2 Configure your Instrument according to the applied hardware changes:

Software Revision B.01.04 and below:

- **Tools > Module Service Center** (any module)

Software Revision B.02.01 and higher:

- **Instrument Control > Controls (any module) > Convert Device Type**

(For example if you installed a G4297A - 1220 Infinity Oven kit then press

Add Oven

3 Disconnect from the software and re-boot your instrument.

4 Re-connect to the Instrument to the Instrument Utility / Lab Advisor software.

5 Now perform the **Installation Check** as described in “[Priming the System and Performing the ‘Installation Check’](#)” on page 49.

6 When starting the Chromatographic Data System use **Auto Configuration** to have your new hardware configuration configured to the CDS.

Priming the System and Performing the 'Installation Check'

The steps described below are performed using the Instrument Utility / Lab Advisor software.

- 1 Connect all channels with HPLC grade water and use the **Purge Pump** for flushing the solvent channels.

Software Revision B.01.04 and below:

- **Tools > Pump > Purge Pump**

Software Revision B.02.01 and higher:

- **Service & Diagnostics > Pump (Tools must be checked) > Purge Pump**

- 2 Prime all connected channels sufficiently until all channels are bubble free.

- 3 Use the **Instrument Control** functionality to prime your system with HPLC grade water to remove air out of the system.

Apply the following conditions:

- Purge Valve: closed
- Flow: 2 mL/min
- Time: 5 min/channel
- Set stroke: 100 µL

- 4 Perform the **Installation Check** from the **Service and Diagnostics** menu and print out the test result.

- 5 Create and print out a **Status Report**.

3 Installation

Performing a 'Checkout Run'

Performing a 'Checkout Run'

1 Start the Chromatographic Data System

2 Create a checkout method with the following parameters:

- Flow: 1 mL/min
- Inj. Volume: 20 µL (Autosampler)
- Oven temp: not controlled
- VWD Wavelength: 254 nm
- Runtime: 1 min

This checkout run is done with the factory installed 'restriction capillary' in place.

3 Prepare 1ml of a checkout sample (acetone for example) and put it on vial position 1 in the autosampler tray.

For Manual Injector configurations load 20 µL of checkout sample in the loop. Overfill the injection loop at least 3 times. (e.g. inject at least 60 µL at a 20 µL sample loop)

4 Start a single run.

As result a single peak should be visible.

5 Print out the report.

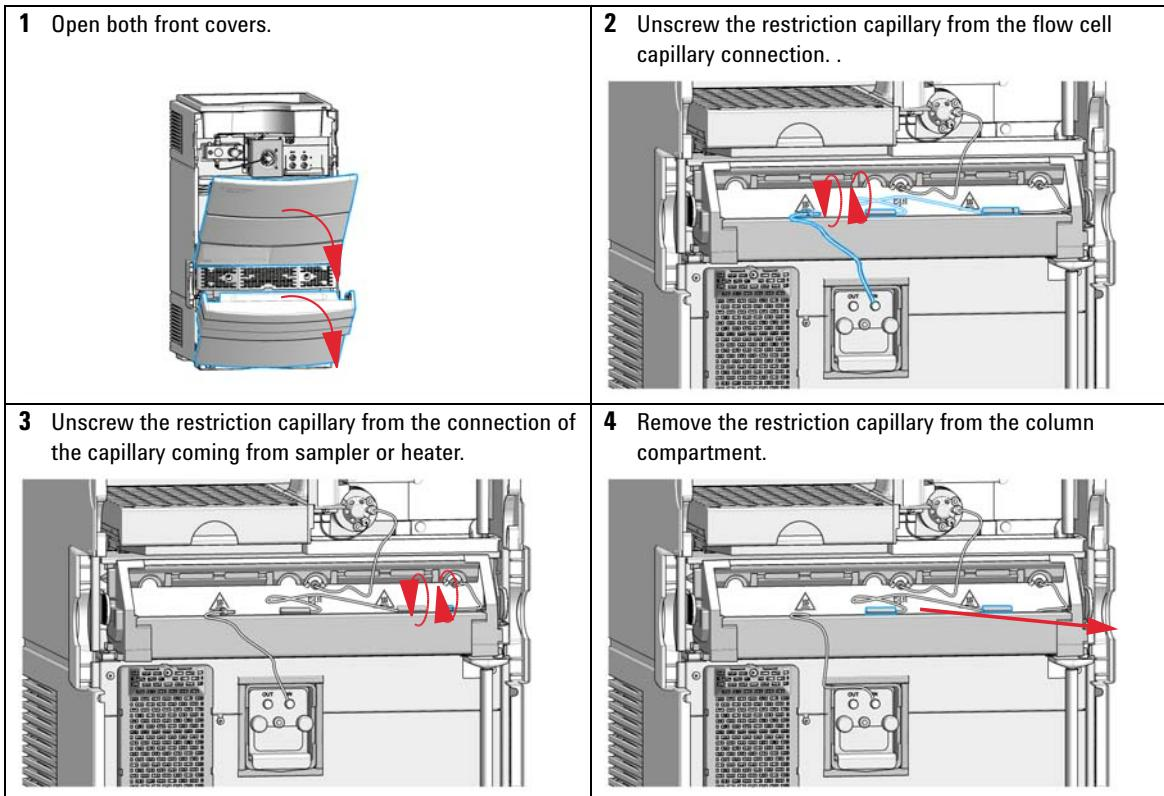
6 Store all created and printed out reports in a binder.

The instrument is installed.

Remove the Restriction Capillary

When For column installation

Tools required	p/n	Description
	8710-0510	Wrench open 1/4 — 5/16 inch

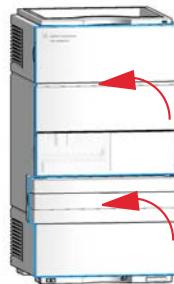


3 Installation

Remove the Restriction Capillary

5 Install a column, see “[Install a Column](#)” on page 53.

6 Close both front covers.



Install a Column

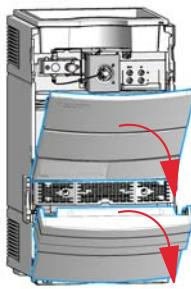
When Installing or changing columns

Tools required **p/n** **Description**
8710-0510 Wrench open 1/4 — 5/16 inch
Flexible wrench

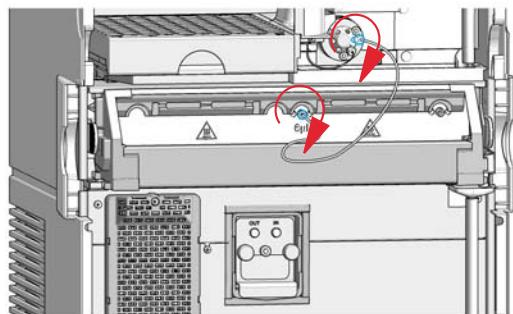
Parts required **Description**
Ordered Column

Preparations Remove the restriction capillary (if installed), see “[Remove the Restriction Capillary](#)” on page 51.

- 1** Open both front covers.



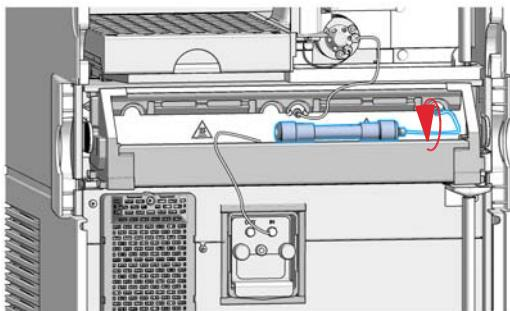
- 2** Connect the column inlet capillary to the injection valve (port 6).



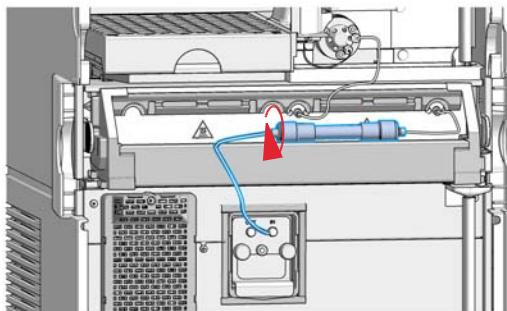
3 Installation

Install a Column

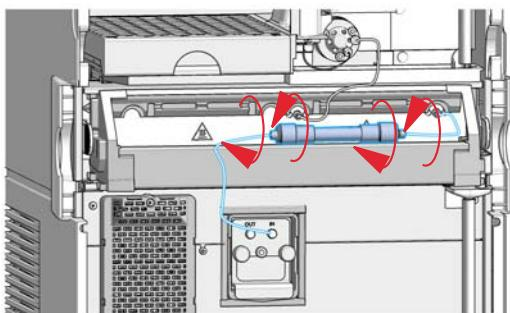
- 3** Connect the capillary coming from the injection valve to the column inlet.



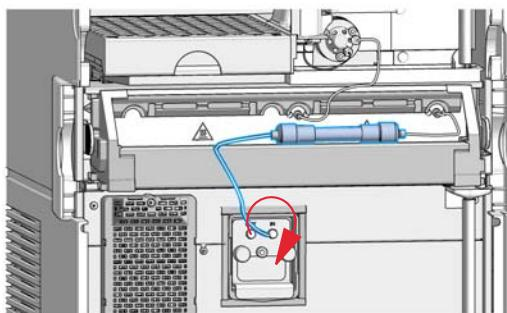
- 4** Connect the column outlet capillary to the column.



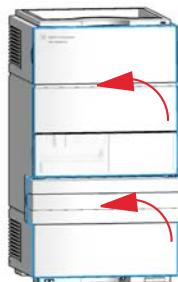
- 5** Use a wrench to counter the column while tightening the capillary fittings on both sides of the column.



- 6** Connect the column outlet capillary to the flow cell inlet at the detector.



- 7** Close both front covers.



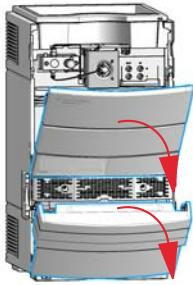
- 8** Configure the column.

Remove a Column

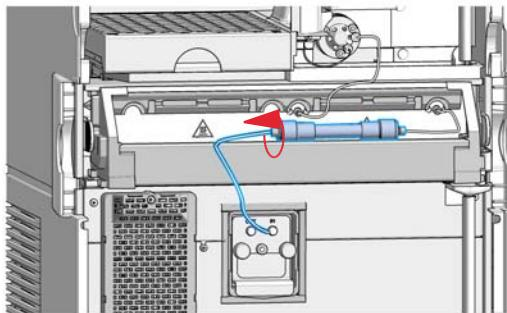
When Installing the restriction capillary or changing the column

Tools required	p/n	Description
	8710-0510	Wrench open 1/4 — 5/16 inch Flexible wrench

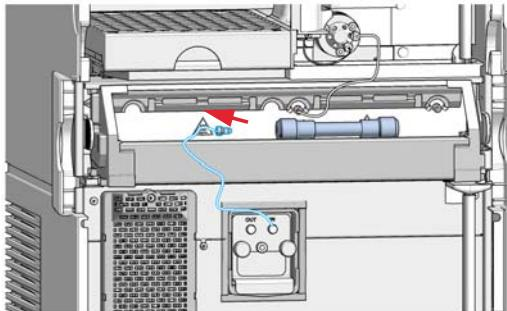
- 1** Open both front covers.



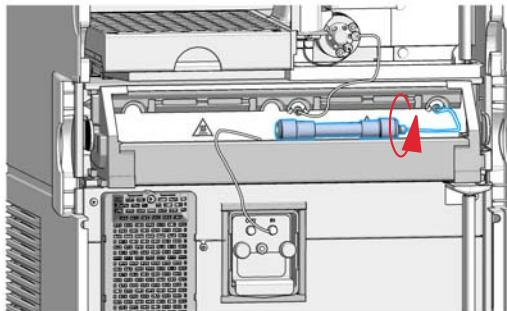
- 2** Unscrew the fitting at the column outlet.



- 3** Disconnect the column outlet capillary from the column.



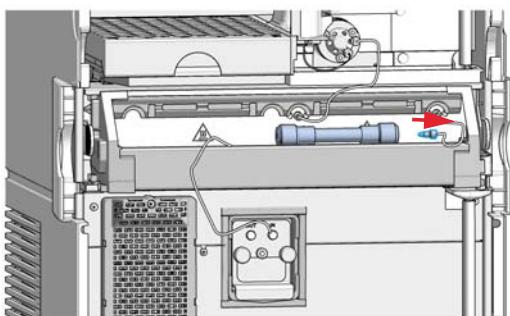
- 4** Unscrew the fitting at the column inlet.



3 Installation

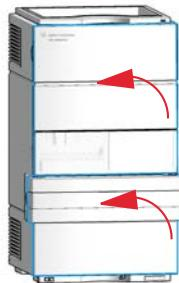
Remove a Column

- 5 Disconnect the column inlet capillary from the column.



- 6 Install the restriction capillary, see "Install the Restriction Capillary" on page 57.

- 7 Close both front covers.



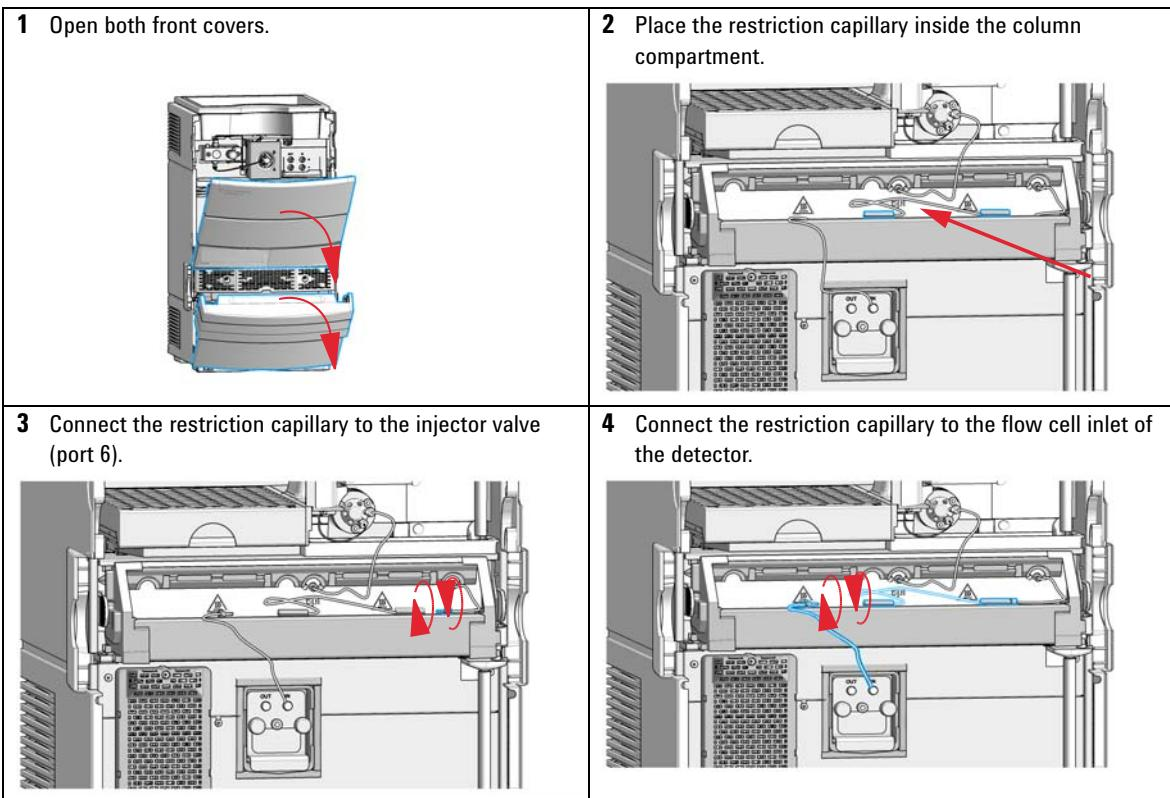
Install the Restriction Capillary

When

Tools required	p/n	Description
	8710-0510	Wrench open 1/4 — 5/16 inch

Parts required	Description
	Restriction capillary

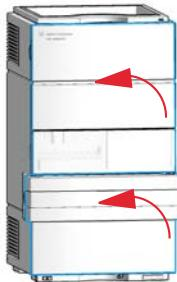
Preparations Remove the column, see “Remove a Column” on page 55.

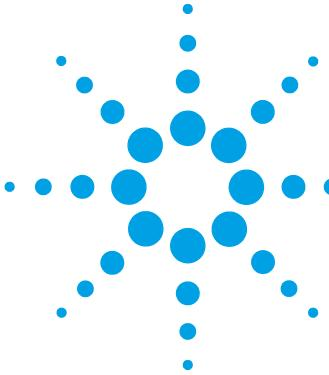


3 Installation

Install the Restriction Capillary

- 5 Close both front covers.





4

LAN Configuration

To do first	60
TCP/IP parameter configuration	62
Configuration Switches	63
Initialization mode selection	64
Dynamic Host Configuration Protocol (DHCP)	66
General Information (DHCP)	66
Setup (DHCP)	67
Link configuration selection	69
Automatic Configuration with BootP	70
About Agilent BootP Service	70
How BootP Service Works	71
Situation: Cannot Establish LAN Communication	71
Installation of BootP Service	71
Two Methods to Determine the MAC Address	73
Assigning IP Addresses Using the Agilent BootP Service	74
Changing the IP Address of an Instrument Using the Agilent BootP Service	77
Storing the settings permanently with Bootp	79
Manual Configuration	80
With Telnet	81

This chapter provides information on connecting the instrument to the Agilent ChemStation PC.



Agilent Technologies

4 LAN Configuration

To do first

To do first

The instrument has an on-board LAN communication interface.

- 1 Note the MAC (Media Access Control) address for further reference. The MAC or hardware address of the LAN interfaces is a world wide unique identifier. No other network device will have the same hardware address. The MAC address can be found on a label at the rear left side of the instrument next to the configuration switch.



Figure 9 MAC label

- 2** Connect the instrument's LAN interface to
- the PC network card using a crossover network cable (point-to-point) or
 - a hub or switch using a standard LAN cable.

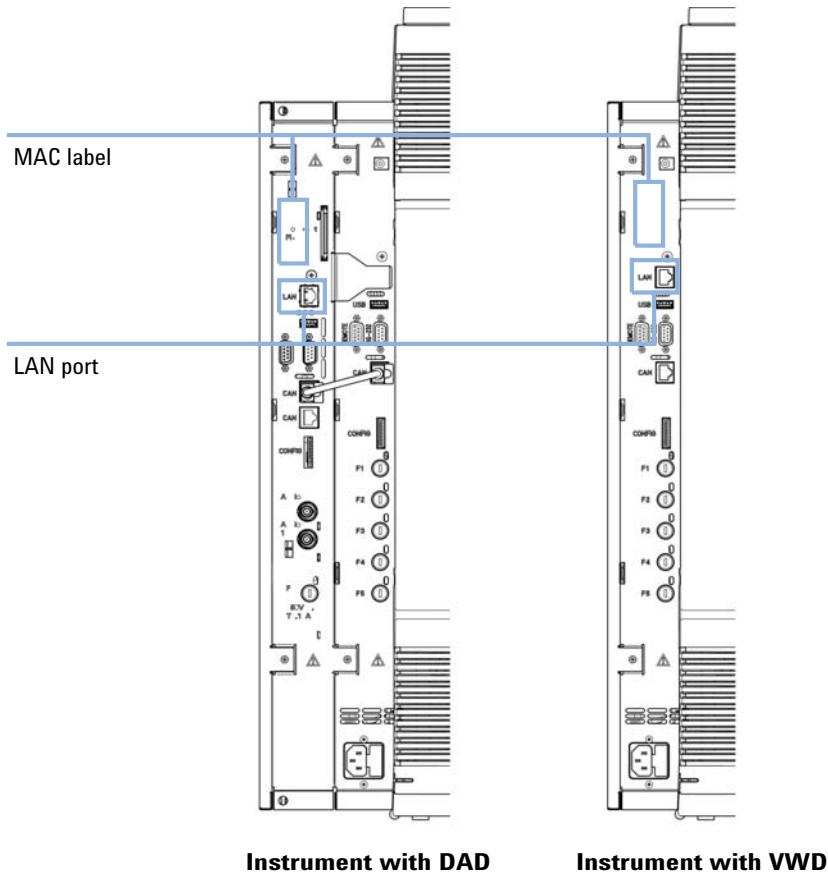


Figure 10 Location of LAN interface and MAC label

4 LAN Configuration

TCP/IP parameter configuration

TCP/IP parameter configuration

To operate properly in a network environment, the LAN interface must be configured with valid TCP/IP network parameters. These parameters are:

- IP address
- Subnet Mask
- Default Gateway

The TCP/IP parameters can be configured by the following methods:

- by automatically requesting the parameters from a network-based DHCP Server (using the so-called Dynamic Host Configuration Protocol). This mode requires a LAN-onboard Module or a G1369C LAN Interface card, see “[Setup \(DHCP\)](#)” on page 67
- by manually setting the parameters using Telnet

The LAN interface differentiates between several initialization modes. The initialization mode (short form ‘init mode’) defines how to determine the active TCP/IP parameters after power-on. The parameters may be derived from non-volatile memory or initialized with known default values. The initialization mode is selected by the configuration switch, see [Table 11](#) on page 64.

Configuration Switches

The configuration switch can be accessed at the rear left side of the instrument.

The instrument is shipped with switches 7 and 8 set to **ON**, which means that the instrument is set to a default fixed IP address: 192.168.254.11

NOTE

To configure the LAN, SW1 and SW2 must be set to **OFF**.

Table 10 Factory Default Settings

Initialization ('Init') Mode	Using Default, switches 7 and 8 set to ON .
Link Configuration	Speed and duplex mode determined by auto-negotiation

NOTE

For the G4294B, the Configuration switches on the DAD main board must be used for configuring the LAN access of the instrument. The switches on the main board must all be set to off.

4 LAN Configuration

Initialization mode selection

Initialization mode selection

The following initialization (init) modes are selectable:

Table 11 Initialization Mode Switches

	SW 6	SW 7	SW 8	Init Mode
	OFF	ON	OFF	Using Stored
	OFF	ON	ON	Using Default
	ON	OFF	OFF	DHCP ¹

¹ Requires firmware B.06.40 or above. Modules without LAN on board, see G1369C LAN Interface Card

Using Stored

When initialization mode **Using Stored** is selected, the parameters are taken from the non-volatile memory of the module. The TCP/IP connection will be established using these parameters. The parameters were configured previously by one of the described methods.

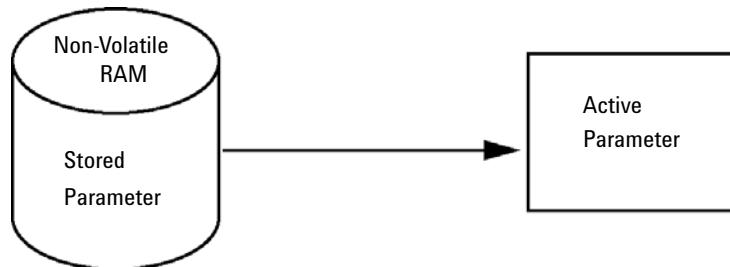


Figure 11 Using Stored (Principle)

Using Default

When **Using Default** is selected, the factory default parameters are taken instead. These parameters enable a TCP/IP connection to the LAN interface without further configuration, see [Table 12](#) on page 65.

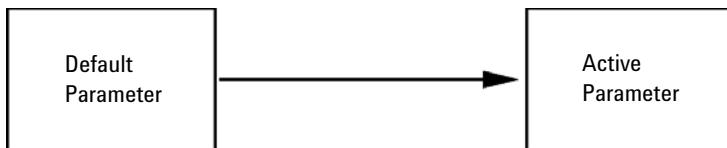


Figure 12 Using Default (Principle)

NOTE

Using the default address in your local area network may result in network problems. Take care and change it to a valid address immediately.

Table 12 Using Default Parameters

IP address:	192.168.254.11
Subnet Mask:	255.255.255.0
Default Gateway	not specified

Since the default IP address is a so-called local address, it will not be routed by any network device. Thus, the PC and the module must reside in the same subnet.

The user may open a Telnet session using the default IP address and change the parameters stored in the non-volatile memory of the module. He may then close the session, select the initialization mode Using Stored, power-on again and establish the TCP/IP connection using the new parameters.

When the module is wired to the PC directly (e.g. using a cross-over cable or a local hub), separated from the local area network, the user may simply keep the default parameters to establish the TCP/IP connection.

NOTE

In the **Using Default** mode, the parameters stored in the memory of the module are not cleared automatically. If not changed by the user, they are still available, when switching back to the mode Using Stored.

4 LAN Configuration

Dynamic Host Configuration Protocol (DHCP)

Dynamic Host Configuration Protocol (DHCP)

General Information (DHCP)

The Dynamic Host Configuration Protocol (DHCP) is an auto configuration protocol used on IP networks. The DHCP functionality is available on all Agilent HPLC modules with on-board LAN Interface or LAN Interface Card G1369C, and "B"-firmware (B.06.40 or above) or modules with "D"-firmware. All modules should use latest firmware from the same set.

When the initialization mode "DHCP" is selected, the card tries to download the parameters from a DHCP Server. The parameters obtained become the active parameters immediately. They are not stored to the non-volatile memory of the card.

Besides requesting the network parameters, the card also submits its hostname to the DHCP Server. The hostname equals the MAC address of the card, e.g. *0030d3177321*. It is the DHCP server's responsibility to forward the hostname/address information to the Domain Name Server. The card does not offer any services for hostname resolution (e.g. NetBIOS).



Figure 13 DHCP (Principle)

NOTE

- 1 It may take some time until the DHCP server has updated the DNS server with the hostname information.
- 2 It may be necessary to fully qualify the hostname with the DNS suffix, e.g. *0030d3177321.country.company.com*.
- 3 The DHCP server may reject the hostname proposed by the card and assign a name following local naming conventions.

Setup (DHCP)

The DHCP functionality is available on all Agilent HPLC modules with on-board LAN Interface or LAN Interface Card G1369C, and "B"-firmware (B.06.40 or above) or modules with "D"-firmware. All modules should use latest firmware from the same set.

- 1 Note the MAC address of the LAN interface (provided with G1369C LAN Interface Card or Main Board). This MAC address is on a label on the card or at the rear of the main board, e.g. *0030d3177321*.

On the Instant Pilot the MAC address can be found under **Details** in the LAN section.

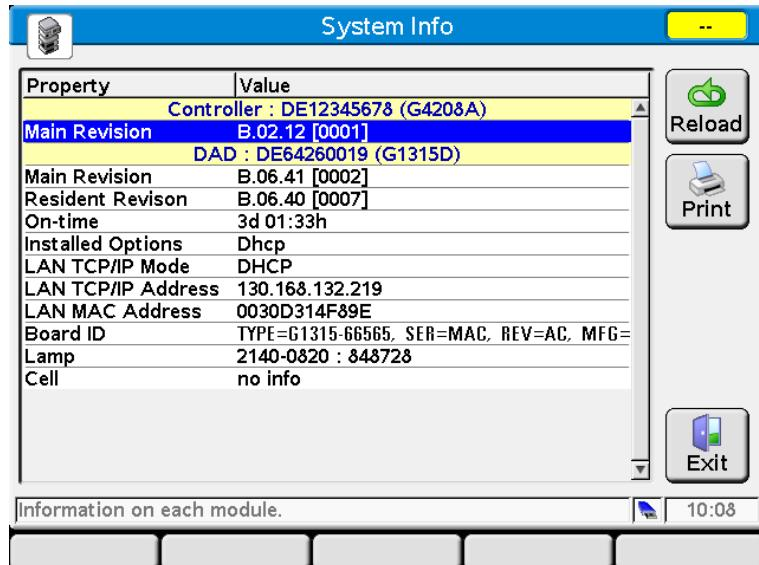


Figure 14 LAN Setting on Instant Pilot

4 LAN Configuration

Dynamic Host Configuration Protocol (DHCP)

- 2** Set the Configuration Switch to DHCP either on the G1369C LAN Interface Card or the main board of above mentioned modules.

Table 13 G1369C LAN Interface Card (configuration switch on the card)

SW 4	SW 5	SW 6	SW 7	SW 8	Initialization Mode
ON	OFF	OFF	OFF	OFF	DHCP

Table 14 LC Modules with 8-bit configuration switch (B-firmware) (configuration switch at rear of the instrument)

SW 6	SW 7	SW 8	Initialization Mode
ON	OFF	OFF	DHCP

- 3** Turn on the module that hosts the LAN interface.
- 4** Configure your Control Software (e.g. OpenLAB CDS ChemStation Edition, Lab Advisor, Firmware Update Tool) and use MAC address as host name, e.g. *0030d3177321*.

The LC system should become visible in the control software (see Note in section “[General Information \(DHCP\)](#)” on page 66).

Link configuration selection

The LAN interface supports 10 or 100 Mbps operation in full- or half-duplex modes. In most cases, full-duplex is supported when the connecting network device - such as a network switch or hub - supports IEEE 802.3u auto-negotiation specifications.

When connecting to network devices that do not support auto-negotiation, the LAN interface will configure itself for 10- or 100-Mbps half-duplex operation.

For example, when connected to a non-negotiating 10-Mbps hub, the LAN interface will be automatically set to operate at 10-Mbps half-duplex.

If the module is not able to connect to the network through auto-negotiation, you can manually set the link operating mode using link configuration switches on the module.

Table 15 Link Configuration Switches

	SW 3	SW 4	SW 5	Link Configuration
	OFF	-	-	speed and duplex mode determined by auto-negotiation
	ON	OFF	OFF	manually set to 10 Mbps, half-duplex
	ON	OFF	ON	manually set to 10 Mbps, full-duplex
	ON	ON	OFF	manually set to 100 Mbps, half-duplex
	ON	ON	ON	manually set to 100 Mbps, full-duplex

4 LAN Configuration

Automatic Configuration with BootP

Automatic Configuration with BootP

NOTE

All examples shown in this chapter will not work in your environment. You need your own IP-, Subnet-Mask- and Gateway addresses.

NOTE

Assure that the detector configuration switch is set properly. The setting should be either **BootP** or **BootP & Store**, see [Table 11](#) on page 64.

NOTE

Assure that the detector connected to the network is powered off.

NOTE

If the Agilent BootP Service program is not already installed on your PC, then install it from your Agilent ChemStation DVD, located in folder **BootP**.

About Agilent BootP Service

The Agilent BootP Service is used to assign the LAN Interface with an IP address.

The Agilent BootP Service is provided on the ChemStation DVD. The Agilent BootP Service is installed on a server or PC on the LAN to provide central administration of IP addresses for Agilent instruments on a LAN. The BootP service must be running TCP/IP network protocol and cannot run a DHCP server.

How BootP Service Works

When an instrument is powered on, an LAN Interface in the instrument broadcasts a request for an IP address or host name and provides its hardware MAC address as an identifier. The Agilent BootP Service answers this request and passes a previously defined IP address and host name associated with the hardware MAC address to the requesting instrument.

The instrument receives its IP address and host name and maintains the IP address as long as it is powered on. Powering down the instrument causes it to lose its IP address, so the Agilent BootP Service must be running every time the instrument powers up. If the Agilent BootP Service runs in the background, the instrument will receive its IP address on power-up.

The Agilent LAN Interface can be set to store the IP address and will not lose the IP address if power cycled.

Situation: Cannot Establish LAN Communication

If a LAN communication with BootP service cannot be established, check the following on the PC:

- Is the BootP service started? During installation of BootP, the service is not started automatically.
- Does the Firewall block the BootP service? Add the BootP service as an exception.
- Is the LAN Interface using the BootP-mode instead of "Using Stored" or "Using Default" modes?

Installation of BootP Service

Before installing and configuring the Agilent BootP Service, be sure to have the IP addresses of the computer and instruments on hand.

- 1 Log on as Administrator or other user with Administrator privileges.
- 2 Close all Windows programs.
- 3 Insert the Agilent ChemStation software DVD into the drive. If the setup program starts automatically, click **Cancel** to stop it.

4 LAN Configuration

Automatic Configuration with BootP

- 4 Open Windows Explorer.
- 5 Go to the BootP directory on the Agilent ChemStation DVD and double-click **BootPPackage.msi**.
- 6 If necessary, click the **Agilent BootP Service...** icon in the task bar.
- 7 The **Welcome** screen of the **Agilent BootP Service Setup Wizard** appears. Click **Next**.
- 8 The **End-User License Agreement** screen appears. Read the terms, indicate acceptance, then click **Next**.
- 9 The **Destination Folder** selection screen appears. Install BootP to the default folder or click **Browse** to choose another location. Click **Next**.

The default location for installation is:

C:\Program Files\Agilent\BootPService\

- 10 Click **Install** to begin installation.

- 11 Files load; when finished, the **BootP Settings** screen appears.

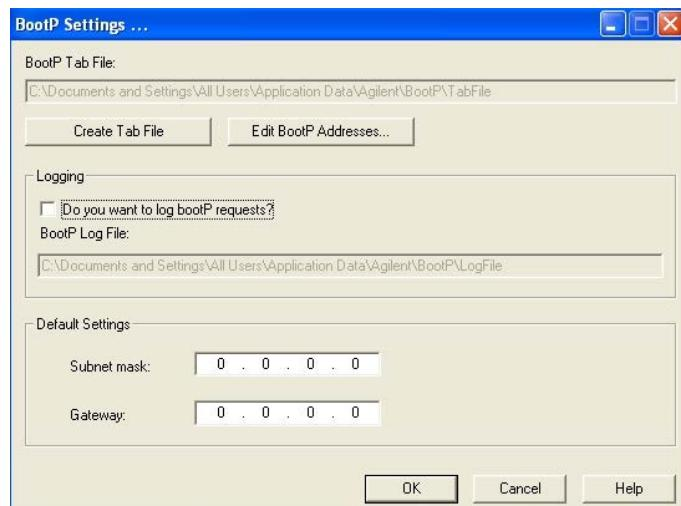


Figure 15 BootP Settings screen

- 12 In the **Default Settings** part of the screen, if known, you can enter the subnet mask and gateway.

Defaults can be used:

- The default subnet mask is 255.255.255.0

- The default gateway is 192.168.254.11

- 13 On the **BootP Settings** screen, click **OK**. The **Agilent BootP Service Setup** screen indicates completion.
- 14 Click **Finish** to exit the **Agilent BootP Service Setup** screen.
- 15 Remove the DVD from the drive.

This completes installation.

- 16 Start BootP Service in the Windows® services: On the Windows® desktop click right on **Computer** icon, select **Manage > Services and Applications > Services**. Select the **Agilent BootP Service** and click **Start**.

Two Methods to Determine the MAC Address

Enabling logging to discover the MAC address using BootP

If you want to see the MAC address, select the **Do you want to log BootP requests?** check box.

- 1 Open BootP Settings from **Start > All Programs > Agilent BootP Service > EditBootPSettings**.
- 2 In **BootP Settings...** check **Do you want to log BootP requests?** to enable logging.

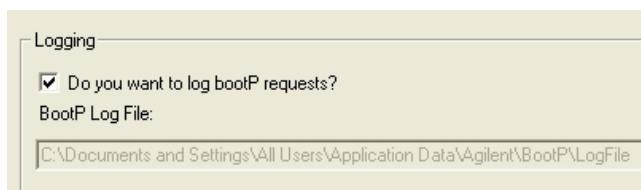


Figure 16 Enable BootP logging

The log file is located in

C:\Documents and Settings\All Users\Application Data\Agilent\BootP\LogFile

It contains a MAC address entry for each device that requests configuration information from BootP.

- 3 Click **OK** to save the values or **Cancel** to discard them. The editing ends.
- 4 After each modification of the BootP settings (i.e. **EditBootPSettings**) a stop or start of the BootP service is required for the BootP service to accept

4 LAN Configuration

Automatic Configuration with BootP

changes. See “[Stopping the Agilent BootP Service](#)” on page 77 or “[Restarting the Agilent BootP Service](#)” on page 78.

- 5 Uncheck the **Do you want to log BootP requests?** box after configuring instruments; otherwise, the log file will quickly fill up disk space.

Determining the MAC address directly from the LAN Interface card label

- 1 Turn off the instrument.
- 2 Read the MAC address from the label and record it.
The MAC address is printed on a label on the rear of the module. It is the number below the barcode and after the colon (:) and usually begins with the letters AD, see [Figure 9](#) on page 60 and [Figure 10](#) on page 61.
- 3 Turn on the instrument.

Assigning IP Addresses Using the Agilent BootP Service

The Agilent BootP Service assigns the Hardware MAC address of the instrument to an IP address.

Determining the MAC address of the instrument using BootP Service

- 1 Power cycle the Instrument.
- 2 After the instrument completes self-test, open the log file of the BootP Service using Notepad.
 - The default location for the logfile is C:\Documents and Settings\All Users\Application Data\Agilent\BootP\LogFile.
 - The logfile will not be updated if it is open.

The contents will be similar to the following:

02/25/10 15:30:49 PM

Status: BootP Request received at outermost layer

Status: BootP Request received from hardware address: 0010835675AC

Error: Hardware address not found in BootPTAB: 0010835675AC

Status: BootP Request finished processing at outermost layer

- 3 Record the hardware (MAC) address (for example, 0010835675AC).
- 4 The Error means the MAC address has not been assigned an IP address and the Tab File does not have this entry. The MAC address is saved to the Tab File when an IP address is assigned.
- 5 Close the log file before turning on another instrument.
- 6 Uncheck the **Do you want to log BootP requests?** box after configuring instruments to avoid having the logfile use up excessive disk space.

Adding each instrument to the network using BootP

- 1 Follow **Start > All Programs > Agilent BootP Service** and select **Edit BootP Settings**. The BootP Settings screen appears.
- 2 Uncheck the **Do you want to log BootP requests?** once all instruments have been added.
The **Do you want to log BootP requests?** box must be unchecked when you have finished configuring instruments; otherwise, the log file will quickly fill up disk space.
- 3 Click **Edit BootP Addresses...** The **Edit BootP Addresses** screen appears.
- 4 Click **Add...** The **Add BootP Entry** screen appears.

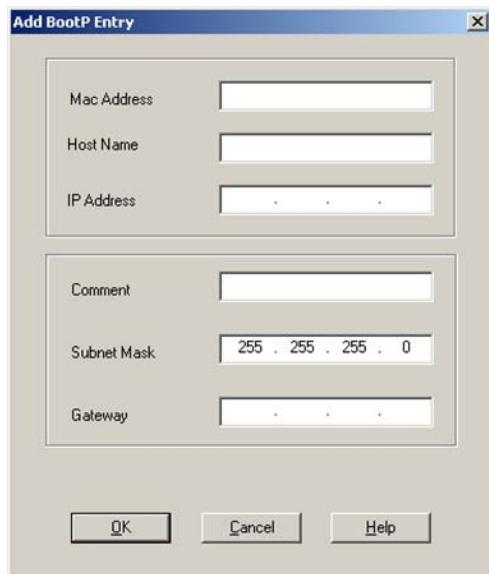


Figure 17 Enable BootP logging

4 LAN Configuration

Automatic Configuration with BootP

5 Make these entries for the instrument:

- MAC address
- Host name, Enter a Hostname of your choice.
The Host Name must begin with "alpha" characters (i.e. LC1260)
- IP address
- Comment (optional)
- Subnet mask
- Gateway address (optional)

The configuration information entered is saved in the Tab File.

6 Click **OK**.

7 Leave **Edit BootP Addresses** by pressing **Close**.

8 Exit **BootP Settings** by pressing **OK**.

9 After each modification of the BootP settings (i.e. EditBootPSettings) a stop or start of the BootP service is required for the BootP service to accept changes. See “[Stopping the Agilent BootP Service](#)” on page 77 or “[Restarting the Agilent BootP Service](#)” on page 78.

10 Power cycle the Instrument.

OR

If you changed the IP address, power cycle the instrument for the changes to take effect.

11 Use the PING utility to verify connectivity by opening a command window and typing:

Ping 192.168.254.11 for example.

The Tab File is located at

C:\Documents and Settings\All Users\Application Data\Agilent\BootP\TabFile

Changing the IP Address of an Instrument Using the Agilent BootP Service

Agilent BootP Service starts automatically when your PC reboots. To change Agilent BootP Service settings, you must stop the service, make the changes, and then restart the service.

Stopping the Agilent BootP Service

- 1 From the Windows control panel, select **Administrative Tools > Services**. The **Services** screen appears.

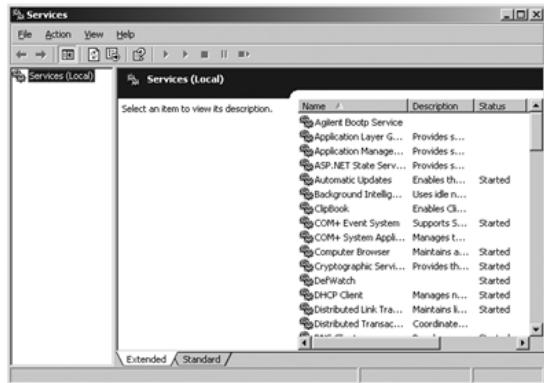


Figure 18 Windows Services screen

- 2 Right-click **Agilent BootP Service**.
- 3 Select **Stop**.
- 4 Close the **Services and Administrative Tools** screen.

Editing the IP address and other parameters in EditBootPSettings

- 1 Select **Start > All Programs > Agilent BootP Service** and select **Edit BootP Settings**. The **BootP Settings** screen appears.
- 2 When the **BootP Settings** screen is first opened, it shows the default settings from installation.

4 LAN Configuration

Automatic Configuration with BootP

- 3 Press **Edit BootP Addresses...** to edit the Tab File.

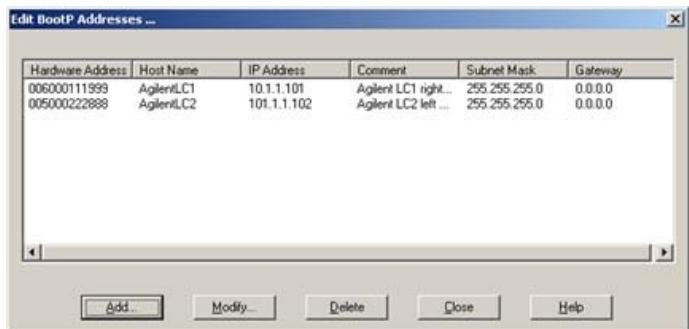


Figure 19 Edit BootP Addresses screen

- 4 In the **Edit BootP Addresses...** screen press **Add...** to create a new entry or select an existing line from the table and press **Modify...** or **Delete** to change the IP address, comment, subnet mask, for example, in the Tab File.
If you change the IP address, it will be necessary to power cycle the instrument for the changes to take effect.
- 5 Leave **Edit BootP Addresses...** by pressing **Close**.
- 6 Exit BootP Settings by pressing OK.

Restarting the Agilent BootP Service

- 1 In the Windows control panel, select **Administrative Tools > Services**. The **Services** screen appears, see [Figure 18](#) on page 77.
- 2 Right-click **Agilent BootP Service** and select **Start**.
- 3 Close the **Services and Administrative Tools** screens.

Storing the settings permanently with Bootp

If you want to change parameters of the module using the Bootp follow the instructions below.

- 1 Turn off the module.
- 2 Change the module's settings of the Configuration Switch to “*Bootp & Store*” mode, see [Table 11](#) on page 64.
- 3 Start the Agilent Bootp Service and open its window.
- 4 If required, modify the parameters for the module according to your needs using the existing configuration.
- 5 Press **OK** to exit the Bootp Manager.
- 6 Now turn on the module and view the Bootp Server window. After some time the Agilent Bootp Service will display the request from the LAN interface. The parameters are now stored permanently in the non-volatile memory of the module.
- 7 Close the Agilent Bootp Service and turn off the module.
- 8 Change the settings of the module's Configuration Switch to “*Using Stored*” mode, see [Table 11](#) on page 64.
- 9 Power cycle the module. The module can be accessed now via LAN without the Agilent Bootp Service.

4 LAN Configuration

Manual Configuration

Manual Configuration

Manual configuration only alters the set of parameters stored in the non-volatile memory of the module. It never affects the currently active parameters. Therefore, manual configuration can be done at any time. A power cycle is mandatory to make the stored parameters become the active parameters, given that the initialization mode selection switches are allowing it.

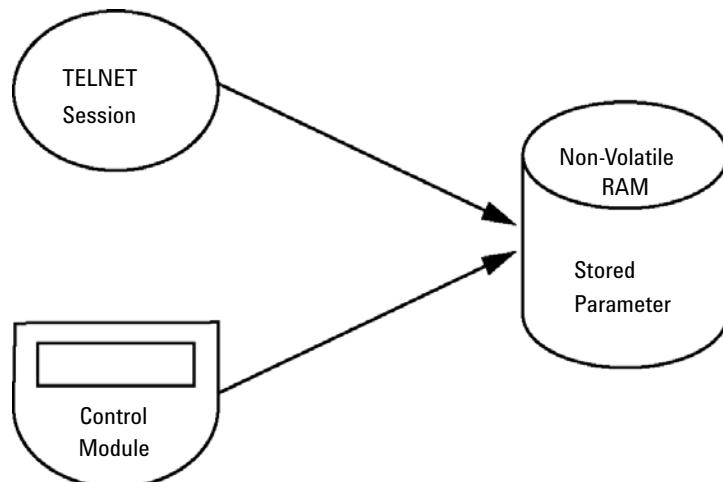


Figure 20 Manual Configuration (Principle)

With Telnet

Whenever a TCP/IP connection to the module is possible (TCP/IP parameters set by any method), the parameters may be altered by opening a Telnet session.

- 1 Open the system (DOS) prompt window by clicking on Windows **START** button and select “**Run...**”. Type “cmd” and press OK.
- 2 Type the following at the system (DOS) prompt:
 - **c:\>telnet <IP address>** or
 - **c:\>telnet <host name>**

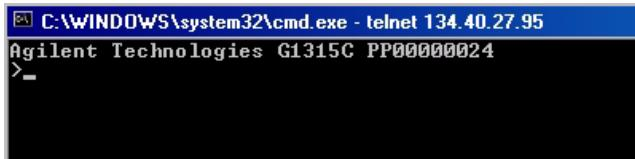


```
C:\WINDOWS\system32\cmd.exe
C:\>telnet 134.40.27.95
```

Figure 21 Telnet - Starting a session

where <IP address> may be the assigned address from a Bootp cycle, a configuration session with the Handheld Controller, or the default IP address (see “[Configuration Switches](#)” on page 63).

When the connection was established successfully, the module responds with the following:



```
C:\WINDOWS\system32\cmd.exe - telnet 134.40.27.95
Agilent Technologies G1315C PP00000024
>_
```

Figure 22 A connection to the module is made

- 3 Type
? and press enter to see the available commands.

4 LAN Configuration

Manual Configuration

command syntax	description
?	display help info
/	display current LAN settings
ip <x.x.x.x>	set IP Address
sm <x.x.x.x>	set Subnet Mask
gw <x.x.x.x>	set Default Gateway
exit	exit shell
>	

Figure 23 Telnet Commands

Table 16 Telnet Commands

Value	Description
?	displays syntax and descriptions of commands
/	displays current LAN settings
ip <x.x.x.x>	sets new ip address
sm <x.x.x.x>	sets new subnet mask
gw <x.x.x.x>	sets new default gateway
exit	exits shell and saves all changes

4 To change a parameter follows the style:

- parameter value, for example:
`ip 134.40.27.230`

Then press [Enter], where parameter refers to the configuration parameter you are defining, and value refers to the definitions you are assigning to that parameter. Each parameter entry is followed by a carriage return.

5 Use the "/" and press Enter to list the current settings.

```
C:\WINDOWS\system32\cmd.exe - telnet 134.40.27.95
>/
LAN Status Page
-----
MAC Address : 0030D30A0838
-----
Init Mode   : Using Stored
-----
TCP/IP Properties
- active -
IP Address  : 134.40.27.95
Subnet Mask : 255.255.248.0
Def. Gateway : 134.40.24.1
-----
TCP/IP Status : Ready
-----
Controllers  : no connections
>_
```

information about the LAN interface
MAC address, initialization mode
Initialization mode is Using Stored
active TCP/IP settings
TCP/IP status - here ready
connected to PC with controller software (e.g. Agilent ChemStation), here not connected

Figure 24 Telnet - Current settings in "Using Stored" mode

6 Change the IP address (in this example 134.40.27.99) and type "/" to list current settings.

```
C:\WINDOWS\system32\cmd.exe - telnet 134.40.27.95
>ip 134.40.27.99
>/
LAN Status Page
-----
MAC Address : 0030D30A0838
-----
Init Mode   : Using Stored
-----
TCP/IP Properties
- active -
IP Address  : 134.40.27.95
Subnet Mask : 255.255.248.0
Def. Gateway : 134.40.24.1
- stored -
IP Address  : 134.40.27.99
Subnet Mask : 255.255.248.0
Def. Gateway : 134.40.24.1
-----
TCP/IP Status : Ready
-----
Controllers  : no connections
>_
```

change of IP setting to
Initialization mode is Using Stored
active TCP/IP settings
stored TCP/IP settings in non-volatile memory

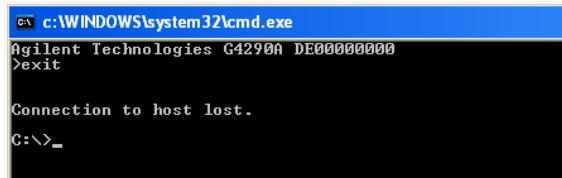
connected to PC with controller software (e.g. Agilent ChemStation), here not connected

Figure 25 Telnet - Change IP settings

4 LAN Configuration

Manual Configuration

- When you have finished typing the configuration parameters, type **exit** and press **Enter** to exit with storing parameters.



```
c:\WINDOWS\system32\cmd.exe
Agilent Technologies G4290A DE00000000
>exit

Connection to host lost.

C:>_
```

Figure 26 Closing the Telnet Session

NOTE

If the Initialization Mode Switch is changed now to “Using Stored” mode, the instrument will take the stored settings when the module is re-booted. In the example above it would be 134.40.27.99.

5

Solvent Delivery System Description

Overview 86

Degasser 87

Principles of Operation 88

Compressibility Compensation 92

Variable Stroke Volume 95

Using the Pump 96

This chapter provides an overview on the operational principles of the solvent delivery system (pump and optional degasser).



Agilent Technologies

5 Solvent Delivery System Description

Overview

Overview

The pump is based on a two-channel, dual-plunger in-series design that provides all essential functions that a solvent delivery system has to fulfill. Metering of solvent and delivery to the high-pressure side are performed by one pump assembly that can generate a pressure up to 600 bar.

The solvents are degassed by a vacuum degasser, and solvent compositions are generated on the low-pressure side by a high-speed proportioning valve. The dual-channel gradient pump includes a built-in dual-channel online vacuum degasser. The isocratic pump of the Agilent 1220 Infinity II LC has no degasser.

The pump assembly includes a pump head with an inlet valve and an outlet valve. A damping unit is connected between the two plunger chambers. A purge valve, including a PTFE frit, is fitted at the pump outlet for convenient priming of the pump head.

Degasser

The dual-channel gradient pump comes with a built-in online degasser. The degasser is switched on automatically when the pump is switched on, even if the flow is set to 0 mL/min. A constant vacuum of 75 Torr (100 mbar) is created in the vacuum chamber of the two channels. The solvent flows through a PTFE AF tube, with an internal volume of 1.5 mL/channel inside the vacuum chamber.

5 Solvent Delivery System Description

Principles of Operation

Principles of Operation

The liquid runs from the solvent reservoir through the degasser to the DCGV, and from there to the inlet valve. The pump assembly comprises two substantially identical plunger/chamber units. Both plunger/chamber units comprise a ball-screw drive and a pump head containing one reciprocating sapphire plunger.

A servo-controlled variable-reluctance motor drives the two ball-screw drives in opposite directions. The gears for the ball-screw drives have different circumferences (ratio 2:1), allowing the first plunger to move at twice the speed of the second plunger. The solvent enters the pump head close to the bottom limit and leaves it at its top. The outer diameter of the plunger is smaller than the inner diameter of the pump head chamber, allowing the solvent to fill the gap in between. The first plunger has a stroke volume in the range of 20 – 100 µL depending on the flow rate. The microprocessor controls all flow rates in a range of 1 µL/min–10 mL/min. The inlet of the first plunger/chamber unit is connected to the inlet valve, which is opened or closed allowing solvent to be drawn into the first plunger pump unit.

The outlet of the first plunger/chamber unit is connected through the outlet ball valve and the damping unit to the inlet of the second plunger/chamber unit. The outlet of the purge valve assembly is then connected to the chromatographic system.

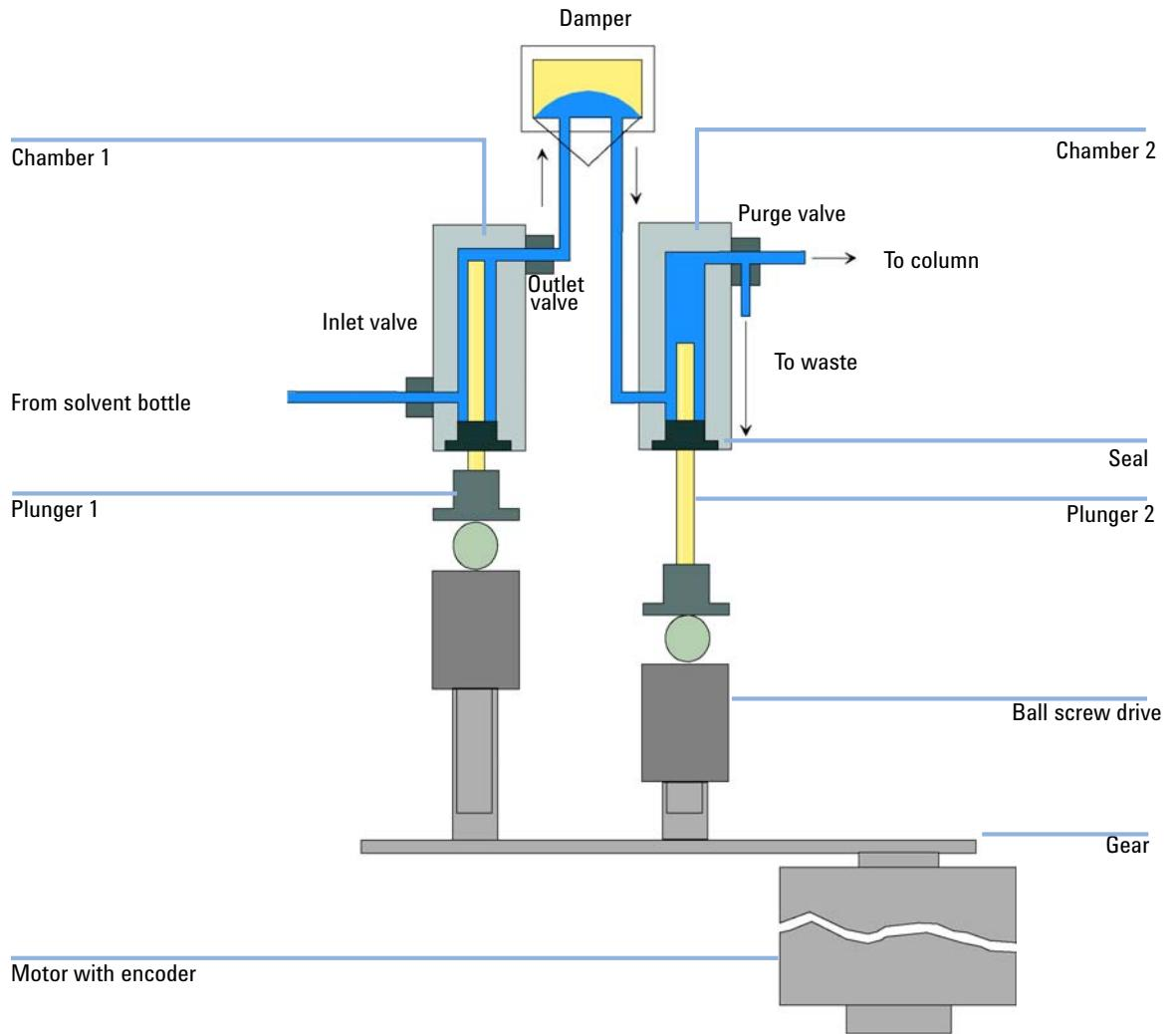


Figure 27 Principle of the pump

When turned on, the pump runs through an initialization procedure to determine the upper dead center of the first plunger. The first plunger moves slowly upwards into the mechanical stop of chamber, and from there it moves back a predetermined distance. The controller stores this plunger position in memory. After this initialization, the pump starts operation with the set

5 Solvent Delivery System Description

Principles of Operation

parameters. The inlet valve is opened and the down-moving plunger draws solvent into the first chamber. At the same time, the second plunger moves upwards, delivering into the system. After a controller-defined stroke length (depending on the flow rate), the drive motor is stopped and the inlet valve is closed. The motor direction is reversed and moves the first plunger up until it reaches the stored upper limit and at the same time moves the second plunger downwards. The sequence then starts again, moving the plungers up and down between the two limits. During the up movement of the first plunger, the solvent in the chamber is pushed through the outlet ball valve into the second chamber. The second plunger draws in half of the volume displaced by the first plunger and the remaining half volume is directly delivered into the system. During the drawing stroke of the first plunger, the second plunger delivers the drawn volume into the system.

For solvent compositions from the solvent bottles A and B, the controller divides the length of the intake stroke into certain fractions in which the gradient valve connects the specified solvent channel to the pump input.

Table 17 Isocratic pump details

Dead volume	800 – 1100 µL, depending on back pressure
Materials in contact with mobile phase	
Pump head	SST, gold, sapphire, ceramic
Active inlet valve	SST, gold, sapphire, ruby, ceramic, PTFE
Outlet valve	SST, gold, sapphire, ruby
Adapter	SST, gold
Purge valve	SST, gold, PTFE, ceramic, PEEK
Degasser chamber	TFE/PDD copolymer, FEP, PEEK, PPS

Table 18 Gradient pump details

Delay volume	800 – 1100 µL, dependent on back pressure
Materials in contact with mobile phase	
MCGV	PTFE
Pump head	SST, gold, sapphire, ceramic
Active inlet valve	SST, gold, sapphire, ruby, ceramic, PTFE
Outlet valve	SST, gold, sapphire, ruby
Adapter	SST, gold
Purge valve	SST, gold, PTFE, ceramic, PEEK
Damping unit	Gold, SST
Degasser chamber	TFE/PDD copolymer, FEP, PEEK, PPS

5 Solvent Delivery System Description

Compressibility Compensation

Compressibility Compensation

Principles of compressibility compensation

The compressibility of the solvents in use affects retention-time stability when the back pressure in the system changes (for example, ageing of the column). To minimize this effect, the pump provides a compressibility compensation feature that optimizes the flow stability according to the solvent type. The compressibility compensation is set to a default value and can be changed through the user interface.

Without compressibility compensation, the following happens during a stroke of the first plunger: the pressure in the plunger chamber increases and the volume in the chamber is compressed, depending on backpressure and solvent type. The volume displaced into the system is reduced by the compressed volume.

When a compressibility value is set, the processor calculates a compensation volume that is depending on the backpressure in the system and the selected compressibility. This compensation volume is added to the normal stroke volume and compensates for the previously described *loss* of volume during the delivery stroke of the first plunger.

Optimizing the compressibility compensation setting

The default compressibility compensation setting is $46 \cdot 10^{-6}$ /bar. This setting represents an average value. Under normal conditions, the default setting reduces the pressure pulsation to values (below 1 % of system pressure) that are sufficient for most applications and for all gradient analyses. For applications using sensitive detectors, the compressibility settings can be optimized by using the values for the various solvents. If the solvent in use is not listed in the compressibility tables, when using isocratic mixtures of solvents and if the default settings are not sufficient for your application, the following procedure can be used to optimize the compressibility settings.

NOTE

When using mixtures of solvents, it is not possible to calculate the compressibility of the mixture by interpolating the compressibility values of the pure solvents used in that mixture or by applying any other calculation. In these cases, the following empirical procedure has to be applied to optimize your compressibility setting.

- 1 Start the pump with the required flow rate.
- 2 Before starting the optimization procedure, the flow must be stable. Use degassed solvent only. Check the tightness of the system with the pressure test.
- 3 Your pump must be connected to control software with which the pressure and %-ripple can be monitored.

5 Solvent Delivery System Description

Compressibility Compensation

- 4 The compressibility compensation setting that generates the smallest pressure ripple is the optimum value for your solvent composition.

Table 19 Solvent Compressibility

Solvent (pure)	Compressibility ($1 \cdot 10^{-6}$ /bar)
Acetone	126
Acetonitrile	115
Benzene	95
Carbon tetrachloride	110
Chloroform	100
Cyclohexane	118
Ethanol	114
Ethyl acetate	104
Heptane	120
Hexane	150
Isobutanol	100
Isopropanol	100
Methanol	120
1-Propanol	100
Toluene	87
Water	46

Variable Stroke Volume

Due to the compression of the pump-chamber volume, each plunger stroke of the pump generates a small pressure pulsation, influencing the flow ripple of the pump. The amplitude of the pressure pulsation is dependent mainly on the stroke volume and the compressibility compensation for the solvent in use. Small stroke volumes generate pressure pulsations of smaller amplitude than larger stroke volumes at the same flow rate. In addition, the frequency of the pressure pulsations are higher. This decreases the influence of flow pulsations on quantitative results.

In gradient mode, smaller stroke volumes result in less flow ripple and improve composition ripple.

The pump uses a processor-controlled spindle system to drive its plungers. The normal stroke volume is optimized for the selected flow rate. Low flow rates use a small stroke volume, while higher flow rates use a larger stroke volume.

When the stroke volume for the pump is set to AUTO mode, the stroke is optimized for the flow rate in use. A change to larger stroke volumes is possible but not recommended.

5 Solvent Delivery System Description

Using the Pump

Using the Pump

Hints for successful use of the Agilent 1220 Infinity II LC pump

- When using salt solutions and organic solvents in the Agilent 1220 Infinity II LC pump, it is recommended to connect the salt solution to one of the lower gradient valve ports and the organic solvent to one of the upper gradient valve port. It is best to have the organic channel directly above the salt solution channel. Regular flushing with water of all DCGV channels is recommended to remove all possible salt deposits in the valve ports.
- Before operating the pump, flush the vacuum degasser (optional) with at least two volumes (3 mL), especially when the pump has been turned off for some time (for example, overnight) and volatile solvent mixtures are used in the channels.
- Prevent blocking of solvent inlet filters (never use the pump without solvent inlet filter). Growth of algae should be avoided.
- Check the purge valve frit and column frit regularly. A blocked purge valve frit can be identified by black or yellow layers on its surface, or by a pressure greater than 10 bar when pumping distilled water at a rate of 5 mL/min with an open purge valve.
- When using the pump at low flow rates (for example, 0.2 mL/min), check all 1/16-inch fittings for any signs of leaks.
- When exchanging the pump seals, also exchange the purge valve frit.
- When using buffer solutions, flush the system with water before switching it off.
- Check the pump plungers for scratches when changing the plunger seals. Scratched plungers will lead to micro leaks and will decrease the lifetime of the seal.
- After changing the plunger seals, pressurize the system according to the wear-in procedure.

Preventing blockage of solvent filters

Contaminated solvents or algae growth in the solvent bottle will reduce the lifetime of the solvent filter and will influence the performance of the pump. This is especially true for aqueous solvents or phosphate buffers (pH 4 – 7). The following suggestions will prolong the lifetime of the solvent filter and will maintain the performance of the pump.

- Use a sterile, if possible amber, solvent bottle to slow down algae growth.
- Filter solvents through filters or membranes that remove algae.
- Exchange solvents every two days, or refilter.
- If the application permits, add 0.0001 – 0.001 M sodium azide to the solvent.
- Place a layer of argon on top of your solvent.
- Avoid exposing the solvent bottle to direct sunlight.

NOTE

Never use the system without a solvent filter installed.

5 Solvent Delivery System Description

Using the Pump

6

Injection System Description

Manual Injector 100

The Injection Seal 100

Injecting Sample 101

Needles 102

Autosampler 103

Sampling Sequence 104

Injection Sequence 105

Sampling Unit 107

Needle-Drive 107

Analytical head 108

Injection Valve 108

Transport Assembly 109

Supported Trays for the Autosampler 110

Choice of Vials and Caps 111

This chapter provides an overview of the operational principles of the injection systems: manual injector and autosampler.



Agilent Technologies

6 Injection System Description

Manual Injector

Manual Injector

The Agilent 1220 Infinity II LC manual injector uses a Rheodyne, 6-port sample injection valve (5067-4202). Sample is loaded into the external 20 μL sample loop through the injection port at the front of the valve. The valve has a PEEK™ injection seal. A make-before-break passage in the stator ensures flow is not interrupted when the valve is switched between the INJECT and LOAD positions, and back again.

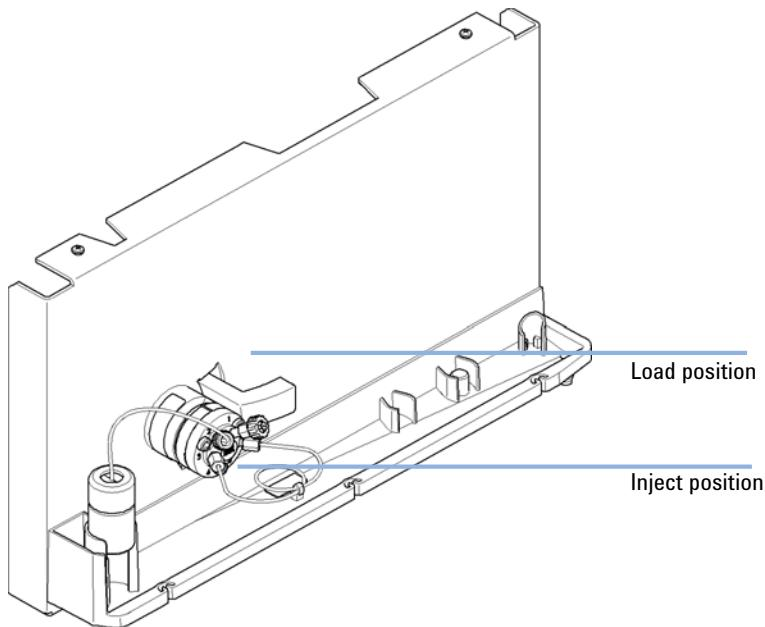


Figure 28 Rheodyne 6-port sample injection valve

The Injection Seal

The manual injector is supplied with a PEEK™ injection seal as standard.

Injecting Sample

WARNING

Ejection of mobile phase

When using sample loops larger than 100 µL, mobile phase may be ejected from the needle port as the mobile phase in the sample loop decompresses.

- Please observe appropriate safety procedures (for example, goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the solvent vendor, especially when toxic or hazardous solvents are used.

LOAD Position

In the LOAD position (see [Figure 29](#) on page 101), the pump is connected directly to the column (ports 4 and 5 connected), and the needle port is connected to the sample loop. At least 2 to 3 sample-loop volumes (more if better precision is required) of sample should be injected through the needle port to provide good precision. The sample fills the loop, and excess sample is expelled through the vent tube connected to port 2.

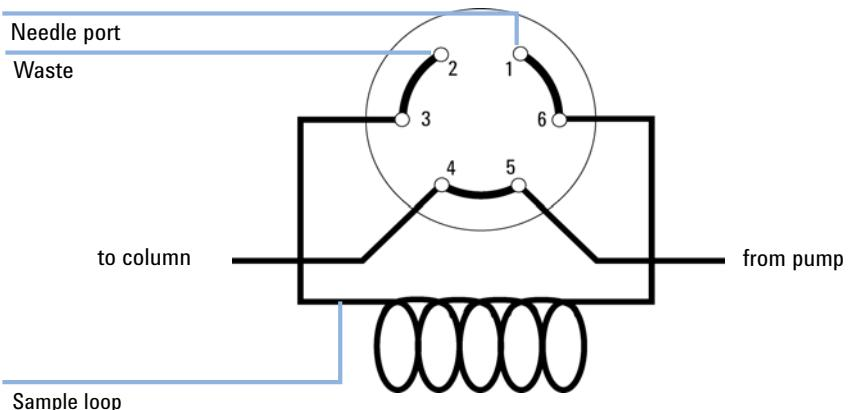


Figure 29 LOAD Position

6 Injection System Description

Manual Injector

INJECT Position

In the INJECT position (see [Figure 30](#) on page 102), the pump is connected to the sample loop (ports 5 and 6 connected). All of the sample is washed out of the loop onto the column. The needle port is connected to the vent tube (port 2).

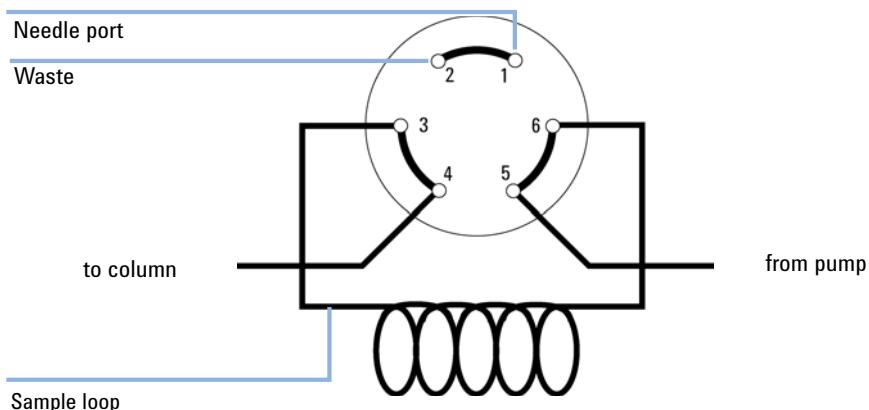


Figure 30 INJECT Position

Needles

CAUTION

Needle can damage valve

→ Always use the correct needle size.

Use needles with 0.028-inch outer diameter (22 gauge) × 2-inch long needle, without electro-taper, and with 90° point style (square tip).

Autosampler

Three sample-rack sizes are available for the autosampler. The standard full-size rack holds $100 \times 1.8\text{ mL}$ vials, while the two half-size racks provide space for $40 \times 1.8\text{ mL}$ and $15 \times 6\text{ mL}$ vials respectively. Any two half-size rack trays can be installed in the autosampler simultaneously. The analytical head device provides injection volumes from $0.1 - 100\text{ }\mu\text{L}$.

The autosamplers transport mechanism uses an X-Z-Theta movement to optimize vial pick-up and return. Vials are picked up by the gripper arm, and positioned below the sampling unit. The gripper transport mechanism and sampling unit are driven by motors. Movement is monitored by optical sensors and optical encoders to ensure correct operation. The metering device is always flushed after injection to ensure minimum carry-over.

The six-port injection valve unit (only 5 ports are used) is driven by a high-speed hybrid stepper motor. During the sampling sequence, the valve unit bypasses the autosampler, and directly connects the flow from the pump to the column. During injection and analysis, the valve unit directs the flow through the autosamplers which ensures that the sample is injected completely into the column, and that any sample residue is removed from the metering unit and needle from before the next sampling sequence begins.

6 Injection System Description

Autosampler

Sampling Sequence

The movements of the autosampler components during the sampling sequence are monitored continuously by the processor. The processor defines specific time windows and mechanical ranges for each movement. If a specific step of the sampling sequence can't be completed successfully, an error message is generated.

Solvent is bypassed from the autosamplers by the injection valve during the sampling sequence. The sample vial is selected by a gripper arm from a static sample rack. The gripper arm places the sample vial below the injection needle. The required volume of sample is drawn into the sample loop by the metering device. Sample is applied to the column when the injection valve returns to the mainpass position at the end of the sampling sequence.

The sampling sequence occurs in the following order:

- 1** The injection valve switches to the bypass position.
- 2** The plunger of the metering device moves to the initialization position.
- 3** The gripper arm selects the vial. At the same time, the needle lifts out of the seat.
- 4** The gripper arm places the vial below the needle.
- 5** The needle lowers into the vial.
- 6** The metering device draws the defined sample volume.
- 7** The needle lifts out of the vial.
- 8** If the automated needle wash is selected, the gripper arm replaces the sample vial, positions the wash vial below the needle, lowers the needle into the vial, then lifts the needle out of the wash vial.
- 9** The gripper arm checks if the safety flap is in position.
- 10** The gripper arm replaces the vial. Simultaneously, the needle lowers into the seat.
- 11** The injection valve switches to the mainpass position.

Injection Sequence

Before the start of the injection sequence, and during an analysis, the injection valve is in the mainpass position. In this position, the mobile phase flows through the autosamplers metering device, sample loop, and needle, ensuring all parts in contact with sample are flushed during the run, thus minimizing carry-over.

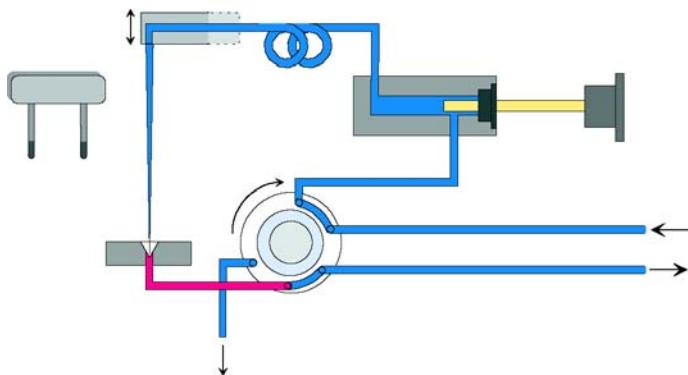


Figure 31 Mainpass Position

When the sample sequence begins, the valve unit switches to the bypass position. Solvent from the pump enters the valve unit at port 1, and flows directly to the column through port 6.

6 Injection System Description

Autosampler

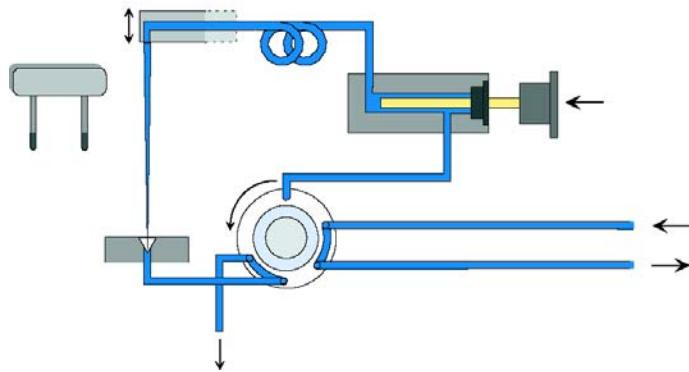


Figure 32 Bypass Position

Next, the needle is raised, and the vial is positioned below the needle. The needle moves down into the vial, and the metering unit draws the sample into the sample loop.

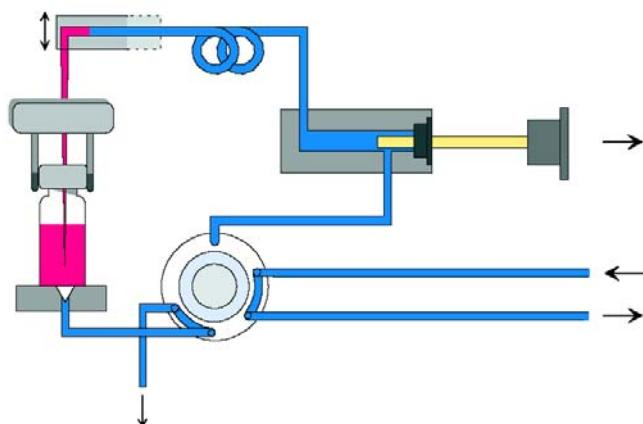


Figure 33 Drawing the Sample

When the metering unit has drawn the required volume of sample into the sample loop, the needle is raised, and the vial is replaced in the sample tray. The needle is lowered into the needle seat, and the injection valve switches back to the mainpass position, flushing the sample onto the column.

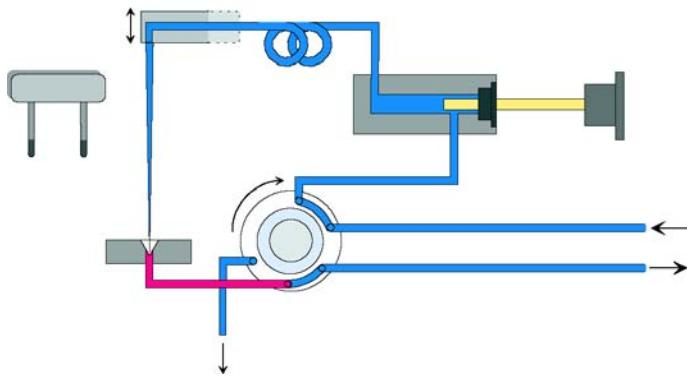


Figure 34 Mainpass Position (Sample Injection)

Sampling Unit

The sampling unit comprises three main assemblies: needle drive, metering device, and injection valve.

NOTE

The replacement sampling unit excludes the injection valve and metering head assemblies.

Needle-Drive

The needle movement is driven by a stepper motor connected to the spindle assembly by a toothed belt. The circular motion of the motor is converted to linear motion by the drive nut on the spindle assembly. The upper and lower needle positions are detected by reflection sensors on the sampling unit flex

6 Injection System Description

Autosampler

board, while the needle-in-vial position is determined by counting the motor steps from the upper needle-sensor position.

Analytical head

The analytical head is driven by the stepper motor connected to the drive shaft by a toothed belt. The drive nut on the spindle converts the circular movement of the spindle to linear motion. The drive nut pushes the sapphire plunger against the tension of the spring into the analytical head. The base of the plunger sits on the large bearing of the drive nut, which ensures the plunger is always centered. A ceramic ring guides the movement of the plunger in the analytical head. The home position of the plunger is sensed by an infra-red sensor on the sampling unit flex board, while the sample volume is determined by counting the number of steps from the home position. The backward movement of the plunger (driven by the spring) draws sample from the vial.

Table 20 Analytical head technical data

Standard (100 µL)	
Number of steps	15000
Volume resolution	7 nL/motor step
Maximum stroke	100 µL
Pressure limit	600 bar
Plunger material	Sapphire

Injection Valve

The two-position 6-port injection valve is driven by a stepper motor. Only five of the six ports are used (port 3 is not used). A lever/slider mechanism transfers the movement of the stepper motor to the injection valve. Two microswitches monitor switching of the valve (bypass and mainpass end positions).

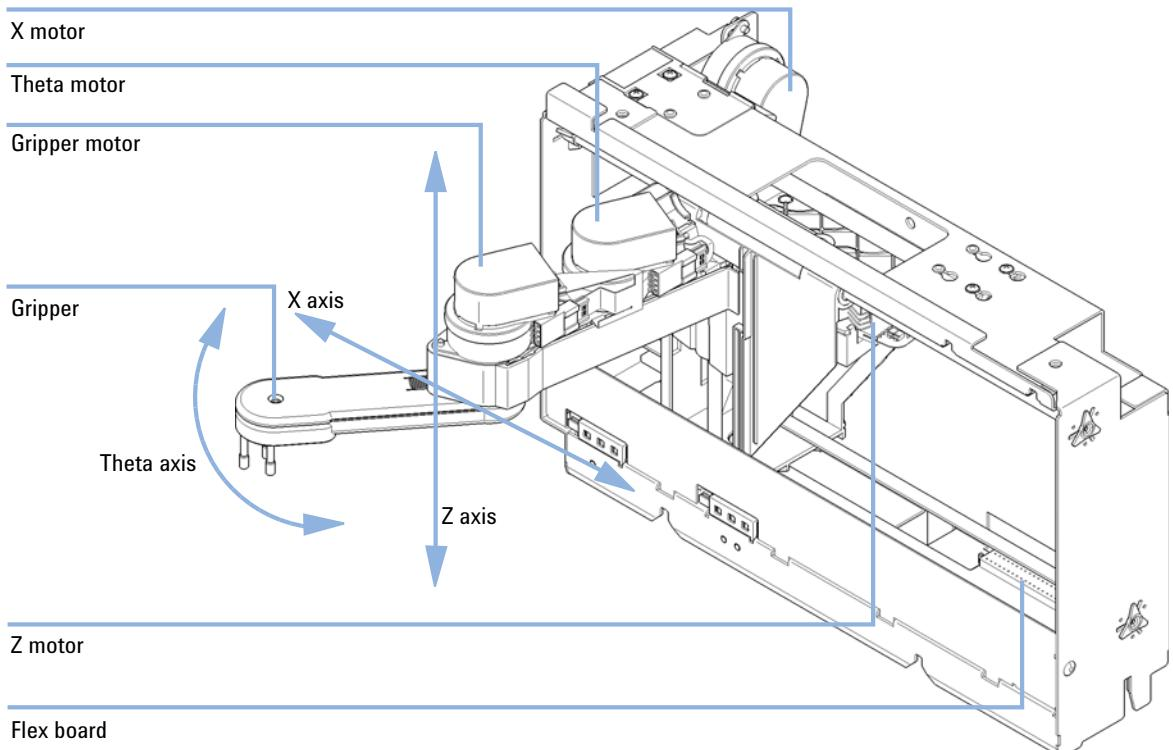
No valve adjustments are required after replacing internal components.

Table 21 Injection valve technical data

	Standard
Motor type	4 V, 1.2 A stepper motor
Seal material	Vespel™ (Tefzel™ available)
Number of ports	6
Switching time	< 150 ms

Transport Assembly

The transport unit comprises an X-axis slide (left-right motion), a Z-axis arm (up-down motion), and a gripper assembly (rotation and vial-gripping).

**Figure 35** Transport Assembly

6 Injection System Description

Autosampler

The transport assembly uses four stepper motors driven in closed-loop mode for accurate positioning of the gripper assembly for sample-vial transport. The rotational movement of the motors is converted to linear motion (X- and Z-axes) by toothed belts connected to the drive spindles. The rotation (theta axes) of the gripper assembly is transferred from the motor by a toothed belt and series of gears. The opening and closing of the gripper fingers are driven by a stepper motor linked by a toothed belt to the planetary gearing inside the gripper assembly.

The stepper motor positions are determined by the optical encoders mounted onto the stepper-motor housing. The encoders monitor the position of the motors continually, and correct for position errors automatically (e.g. if the gripper is accidentally moved out of position when loading vials into the vial tray). The initialization positions of the moving components are sensed by reflection sensors mounted on the flex board. These positions are used by the processor to calculate the actual motor position. An additional six reflection sensors for tray recognition are mounted on the flex board at the front of the assembly.

Supported Trays for the Autosampler

p/n	Description
G1313-44510	Tray for 100 x 2 mL vials
G1313-44513	Halftray for 15 x 6 mL vials
G1313-44512	Halftray for 40 x 2 mL vials

Half-tray combinations

Half-trays can be installed in any combination enabling both 2 mL- and 6 mL-vials to be used simultaneously.

Numbering of vial positions

The standard 100-vial tray has vial positions 1 to 100. However, when using two half-trays, the numbering convention is slightly different. The vial positions of the right-hand half tray begin at position 101 as follows:

Left-hand 40-position tray: 1 – 40

Left-hand 15-position tray: 1 – 15

Right-hand 40-position tray: 101 – 140

Right-hand 15-position tray: 101 – 115

Choice of Vials and Caps

For reliable operation, vials used with the Agilent 1220 Infinity II LC autosampler must not have tapered shoulders or caps that are wider than the body of the vial. The vials and caps shown with their part numbers in the tables below have been successfully tested using a minimum of 15000 injections with the Agilent 1220 Infinity II LC autosampler.

Crimp top vials

p/n	Description
5181-3375	Crimp Top Vial, 2 mL, clear glass, 100/Pack
5183-4491	Crimp Top Vial, 2 mL, clear glass, 1000/Pack
5182-0543	Crimp Top Vial, 2 mL, clear glass, write-on spot, 100/Pack
5183-4492	Crimp Top Vial, 2 mL, clear glass, write-on spot, 1000/Pack
5183-4494	Crimp Top Vial, 2 mL, clear glass, write-on spot, 100/Pack (silanized)
5181-3376	Crimp Top Vial, 2 mL, amber glass, write-on spot, 100/Pack
5183-4493	Crimp Top Vial, 2 mL, amber glass, write-on spot, 1000/Pack
5183-4495	Crimp Top Vial, 2 mL, amber glass, write-on spot, 100/Pack (silanized)
5182-0567	Crimp Top Vial, 1 mL, polypropylene, wide opening, 100/Pack
5183-4496	Crimp Top Vial, 1 mL, polypropylene, wide opening, 100/Pack (silanized)
9301-0978	crimp top vial, 0.3 mL, polypropylene, wide opening, 1000/pack

6 Injection System Description

Autosampler

Snap top vials

p/n	Description
5182-0544	Snap Top Vial, 2 mL, clear glass, 100/Pack
5183-4504	Snap Top Vial, 2 mL, clear glass, 1000/Pack
5183-4507	Snap Top Vial, 2 mL, clear glass, 100/Pack (silanized)
5182-0546	Snap Top Vial, 2 mL, clear glass, write-on spot, 100/Pack
5183-4505	Snap Top Vial, 2 mL, clear glass, write-on spot, 1000/Pack
5183-4508	Snap Top Vial, 2 mL, clear glass, write-on spot, 100/Pack (silanized)
5182-0545	Snap Top Vial, 2 mL, amber glass, write-on spot, 100/Pack
5183-4506	Snap Top Vial, 2 mL, amber glass, write-on spot, 1000/Pack
5183-4509	Snap Top Vial, 2 mL, amber glass, write-on spot, 100/Pack (silanized)

Screw top vials

p/n	Description
5182-0714	Screw Cap Vials, 2 mL, clear glass, 100/Pack
5183-2067	Screw Top Vial, 2 mL, clear glass, 1000/Pack
5183-2070	Screw Top Vial, 2 mL, clear glass, 100/Pack (silanized)
5182-0715	Screw Top Vial, 2 mL, clear glass, write-on spot, 100/Pack
5183-2068	Screw Top Vial, 2 mL, clear glass, write-on spot, 1000/Pack
5183-2071	Screw Top Vial, 2 mL, clear glass, write-on spot, 100/Pack (silanized)
5182-0716	Screw Cap Vial, 2 mL, amber glass, write-on spot, 100/pk
5183-2069	Screw Top Vial, 2 mL, amber glass, write-on spot, 1000/Pack
5183-2072	Screw Top Vial, 2 mL, amber glass, write-on spot, 100/Pack (silanized)

Crimp caps

p/n	Description
5181-1210	Crimp Cap, silver aluminum, septum (clear PTFE/red rubber), 100/Pack
5183-4498	Crimp Cap, silver aluminum, septum (clear PTFE/red rubber), 1000/Pack
5181-1215	Crimp Cap, blue aluminum, septum (clear PTFE/red rubber), 100/Pack
5181-1216	Crimp Cap, green aluminum, septum (clear PTFE/red rubber), 100/Pack
5181-1217	Crimp Cap, red aluminum, septum (clear PTFE/red rubber), 100/Pack

Snap caps

p/n	Description
5182-0550	Snap Cap, clear polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-3458	Snap Cap, blue polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-3457	Snap Cap, green polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-3459	Snap Cap, red polypropylene, septum (clear PTFE/red rubber), 100/Pack

6 Injection System Description

Autosampler

Screw caps

p/n	Description
5182-0717	Screw Cap, blue polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-0718	Screw Cap, green polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-0719	Screw Cap, red polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-0720	Screw Cap, blue polypropylene, septum (clear PTFE/silicone), 100/Pack
5182-0721	Screw Cap, green polypropylene, septum (clear PTFE/silicone), 100/Pack
5182-0722	Screw Cap, red polypropylene, septum (clear PTFE/silicone), 100/Pack

7

Column Oven Description

Column Oven 116

This chapter provides an overview of the operational principles of the column oven.



Agilent Technologies

115

7 Column Oven Description

Column Oven

Column Oven

The column oven is based on a resistor heater matt with two thermal sensors to provide constant temperature in the whole column area. A built-in over temperature cut off fuse inhibits overheating.

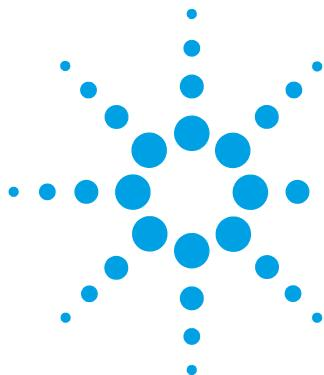
The inner volume of the oven capillary is 6 µL.

Maximum column length is 25 cm (10 inch).

Operational range is at least 10 °C up to 80 °C, max. specified flow rate is 5 mL/min at 60 °C.

NOTE

Never operate the column oven with open front cover, to ensure a correct column temperature always operate with closed front cover. The counterpart of the oven isolation is fixed at the inner side of the front cover.



8

Detector Description

Detector Types	118
Agilent 1220 Infinity II LC Variable Wavelength Detector (VWD) Detector	119
Agilent 1220 Infinity II LC Diode Array Detector (DAD)	120
Introduction to the Detector	120
Optical System	120
Peak width (response time)	123
Sample and Reference Wavelength and Bandwidth	124
Slit Width	127
Optimizing Spectral Acquisition (DAD only)	129
Margin for Negative Absorbance	130
Optimizing Selectivity	130
Spectrum Settings (DAD only)	134
Match the Flow Cell to the Column	136

This chapter provides an overview of the operational principles of the detector.



Agilent Technologies

8 Detector Description

Detector Types

Detector Types

There are two different detector types available for the Agilent 1220 Infinity II LC System:

- Variable Wavelength Detector (VWD, used in G4286B, G4288B/C, G4290B/C), optical unit of the G1314F VWD
- Diode Array Detector (DAD, used in G4294B), optical unit of the G1315C DAD

Agilent 1220 Infinity II LC Variable Wavelength Detector (VWD)

Detector

The Agilent 1220 Infinity II LC variable wavelength detector is designed for highest optical performance, GLP compliance and easy maintenance, with:

- Deuterium lamp for highest intensity and lowest detection limit over a wavelength range of 190 – 600 nm,
- Optional flow-cell cartridges (standard: 10 mm 14 µL, high pressure: 10 mm 14 µL, micro: 3 mm 2 µL, semi-micro: 6 mm 5 µL) are available and can be used depending on the application needs,
- Easy front access to lamp and flow cell for fast replacement, and
- Built-in holmium oxide filter for fast wavelength accuracy verification.

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

Introduction to the Detector

The detector is designed for highest optical performance, GLP compliance and easy maintenance. It includes the following features:

- 80 Hz data acquisition rate for (ultra-) fast LC applications
- RFID tags for all flow cells and UV-lamps provides traceable information about these assemblies
- Long-life deuterium with RFID tag and tungsten lamps for highest intensity and lowest detection limit over a wavelength range of 190 – 950 nm
- No loss in sensitivity for up to eight wavelengths simultaneous
- Programmable slit from 1 – 16 nm for complete optimization of sensitivity, linearity and spectral resolution
- Optional flow-cell cartridges with RFID tag (standard 10 mm¹ 13 µL, semi-micro 6 mm⁵ µL, micro 3 mm² µL, 80 nL, 500 nL, 10 mm, high pressure 10 mm¹.7 µL and prep-cells) are available and can be used depending on the application needs.
- Easy front access to lamps and flow cell for fast replacement
- Built-in holmium oxide filter for fast wavelength accuracy verification
- Built-in temperature control for improved baseline stability
- Additional diagnostic signals for temperature and lamp voltage monitoring

For specifications, see “[Performance Specifications Agilent 1220 Infinity II LC DAD](#)” on page 26.

Optical System

The optical system of the detector is shown in Figure below. Its illumination source is a combination of a deuterium-arc-discharge lamp for the ultraviolet (UV) wavelength range and a tungsten lamp for the visible (VIS) and short-wave near-infrared (SWNIR) wavelength range. The image of the filament of the tungsten lamp is focused on the discharge aperture of the

deuterium lamp by means of a special rear-access lamp design which allows both light sources to be optically combined and share a common axis to the source lens. The achromat (source lens) forms a single, focused beam of light through the flow cell. Each cell room and lamp are separated by a quartz window which can be cleaned or replaced. In the spectrograph, light is being dispersed onto the diode array by a holographic grating. This allows simultaneous access to all wavelength information.

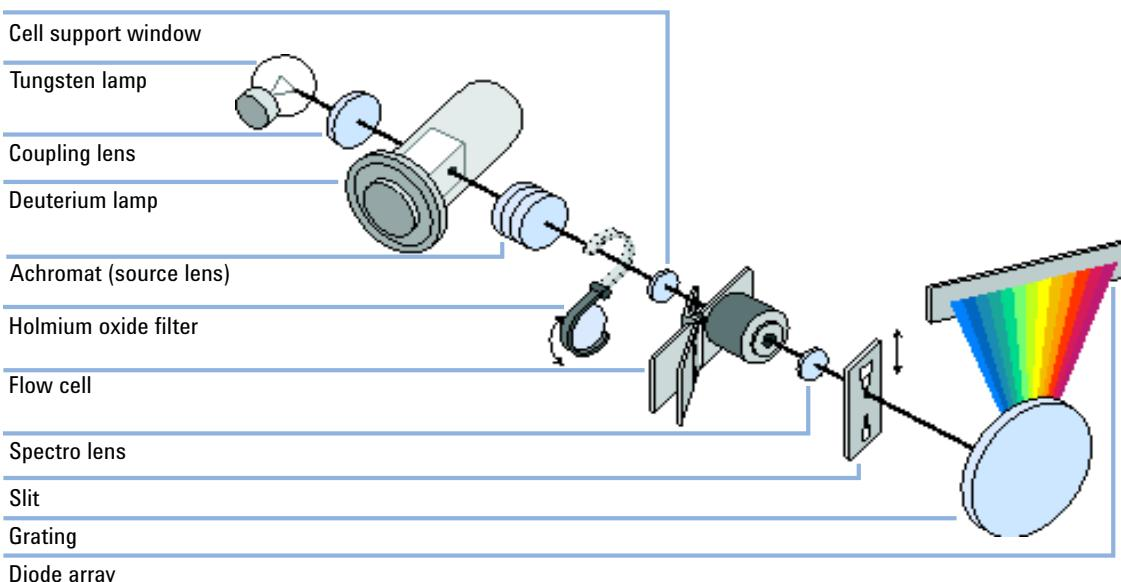


Figure 36 Optical System of the Detector

Lamps The light source for the UV-wavelength range is a deuterium lamp with a shine-through aperture. As a result of plasma discharge in low-pressure deuterium gas, the lamp emits light over the 190 nm to approximately 800 nm wavelength range. The light source for the visible and SWNIR wavelength range is a low noise tungsten lamp. This lamp emits light over the wavelength range 470 – 950 nm.

Achromat (Source Lens) The achromat receives the light from both lamps and focuses it so that the beam passes through the flow cell.

Holmium Oxide Filter The holmium oxide filter is electromechanically actuated. During the holmium filter test it moves into the light path.

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

Cell Support Window	The cell support window assembly separates the holmium filter area from the flow cell area.
Flow Cell Compartment	The optical unit has a flow cell compartment for easy access to flow cells. A variety of optional flow cells can be inserted using the same quick, simple mounting system. The flow cell can be removed to check the optical and electronic performance of the detector without having influences from the flow cell.
Spectrograph	<p>The spectrograph material is ceramic to reduce thermal effects to a minimum. The spectrograph consists of the spectrograph lens, the variable entrance slit, the grating and the photodiode array with front-end electronics. The spectrograph lens refocuses the light beam after it has passed through the flow cell. The sampling interval of the diode array is < 1 nm over the wavelength range 190 – 950 nm. Depending on the wavelength this varies from 1.0 to 1.25 diodes per nanometer (for example a diode every 0.8 to 1 nm).</p> <p>For a small wavelength range, the small non-linearity could be neglected. With the wavelength range from 190 – 950 nm a new approach is required to achieve wavelength accuracy over the full range. Each spectrograph is calibrated individually. The calibration data is stored in the spectrograph on an EEPROM. Based on these data, the built-in processors calculate absorbance data with linear intervals (1.0, 2.0, ...) between data points. This results in an excellent wavelength accuracy and instrument-to-instrument reproducibility.</p>
Variable Entrance Slit System	The micro-slit system makes use of the mechanical properties of silicon combined with the precise structuring capabilities of bulk micro-machining. It combines the required optical functions – slit and shutter – in a simple and compact component. The slit width is directly controlled by the micro-processor of the instrument and can be set as method parameter.
Grating	The combination of dispersion and spectral imaging is accomplished by using a concave holographic grating. The grating separates the light beam into all its component wavelengths and reflects the light onto the photodiode array.
Diode Array	The diode array is a series of 1024 individual photodiodes and control circuits located on a ceramic carrier. With a wavelength range from 190 – 950 nm the sampling interval is < 1 nm.

Peak width (response time)

Response time describes how fast the detector signal follows a sudden change of absorbance in the flow cell. The detector uses digital filters to adapt response time to the width of the peaks in your chromatogram. These filters do not affect peak area nor peak symmetry. When set correctly, such filters reduce baseline noise significantly ([Figure 37](#) on page 123), but reduce peak height only slightly. In addition, these filters reduce the data rate to allow optimum integration and display of your peaks and to minimize disk space required to store chromatograms and spectra.

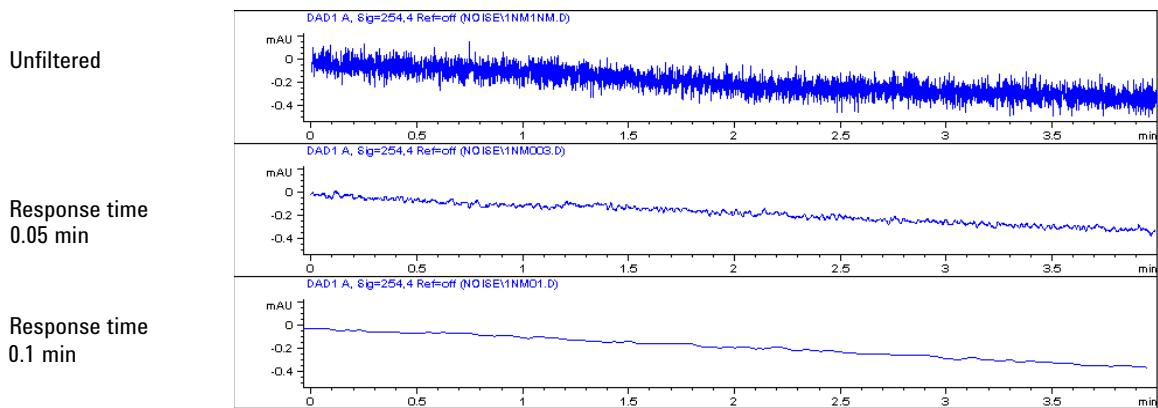


Figure 37 Influence of Response Time on Signal and Noise

[Table 22](#) on page 124 lists the filter choices of the detector. To get optimum results, set peak width as close as possible to a narrow peak of interest in your chromatogram. Response time will be approximately 1/3 of the peak width, resulting in less than 5 % peak-height reduction and less than 5 % additional peak dispersion. Decreasing the peak width setting in the detector will result in less than 5 % gain in peak height but baseline noise will increase by a factor of 1.4 for a factor of 2 response-time reduction. Increasing peak width (response time) by factor of two from the recommended setting (over-filtering) will reduce peak height by about 20 % and reduce baseline noise by a factor of 1.4. This gives you the best possible signal-to-noise ratio, but may affect peak resolution.

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

Table 22 Peak Width — Response Time — Data Rate

Peak Width [minutes]	Response Time [seconds]	Data Rate [Hz]
<0.0025	0.025	80
>0.0025	0.05	80
>0.005	0.1	40
>0.01	0.2	20
>0.03	0.5	10
>0.05	1.0	5
>0.10	2.0	2.5
>0.20	4.0	1.25
>0.40	8.0	0.62
>0.85	16.0	0.31

Sample and Reference Wavelength and Bandwidth

The detector measures absorbance simultaneously at wavelengths from 190 to 950 nm. Two lamps provide good sensitivity over the whole wavelength range. The deuterium discharge lamp provides the energy for the UV range (190 to 400 nm) and the tungsten lamp emits light from 400 to 950 nm for the visible and short wave near infrared.

If you know little about the analytes in your sample, use both lamps and store all spectra over the full wavelength range. This provides full information but fills up your disk space rather quickly. Spectra can be used to check a peak's purity and identity. Spectral information is also useful to optimize wavelength settings for your chromatographic signal.

The detector can compute and store at run time up to 8 signals with these properties:

- sample wavelength, the center of a wavelength band with the width of sample bandwidth (BW), and optionally
- reference wavelength, the center of a wavelength band with the width of reference bandwidth.

The signals comprises a series of data points over time, with the average absorbance in the sample wavelength band minus the average absorbance of the reference wavelength band.

Signal A in the detector default method is set to sample 250,100, reference 360,100, that is, the average absorbance from 200 – 300 nm minus the average absorbance from 300 – 400 nm. As all analytes show higher absorbance at 200 – 300 nm than at 300 – 400 nm, this signal will show you virtually every compound which can be detected by UV absorbance.

Many compounds show absorbance bands in the spectrum. [Figure 38](#) on page 125 shows the spectrum of anisic acid as an example.

To optimize for lowest possible detectable concentrations of anisic acid, set the sample wavelength to the peak of the absorbance band (that is, 252 nm) and the sample bandwidth to the width of the absorbance band (that is, 30 nm). A reference of 360,100 is adequate. Anisic acid does not absorb in this range.

If you work with high concentrations, you may get better linearity above 1.5 AU by setting the sample wavelength to a valley in the spectrum, like 225 nm for anisic acid.

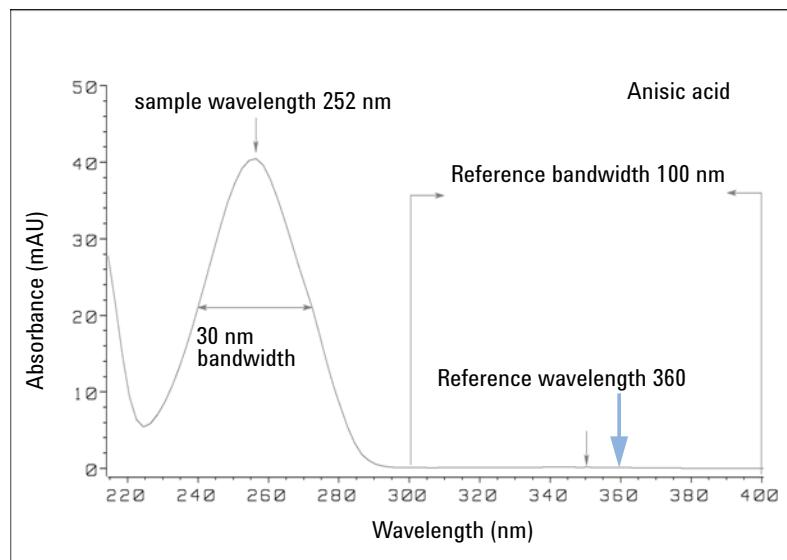


Figure 38 Optimization of Wavelength Setting

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

A wide bandwidth has the advantage of reducing noise by averaging over a wavelength range – compared to a 4 nm bandwidth, the baseline noise is reduced by a factor of approximately 2.5, whereas the signal is about 75 % of a 4 nm wide band. The signal-to-noise ratio for a 30 nm bandwidth is twice that for a 4 nm bandwidth in our example.

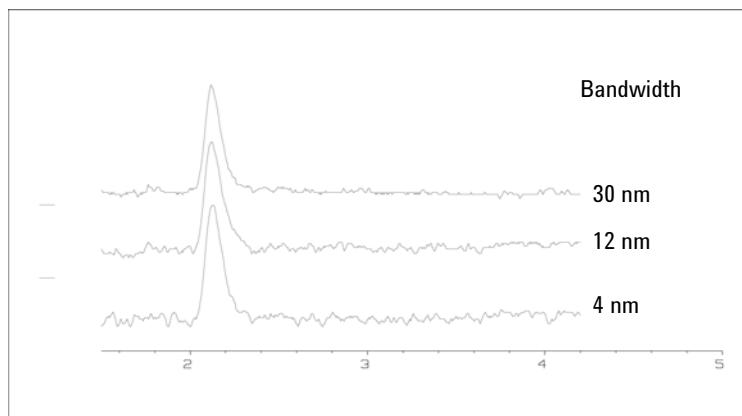


Figure 39 Influence of Bandwidth on Signal and Noise

Because the detector averages absorbance values that are calculated for each wavelength, using a wide bandwidth does not negatively impact linearity.

The use of a reference wavelength is highly recommended to further reduce baseline drift and wander induced by room temperature fluctuations or refractive index changes during a gradient.

An example of the reduction of baseline drifts is shown in [Figure 40](#) on page 127 for PTH-amino acids. Without a reference wavelength, the chromatogram drifts downwards due to refractive index changes induced by the gradient. This is almost completely eliminated by using a reference wavelength. With this technique, PTH-amino acids can be quantified in the low picomole range even in a gradient analysis.

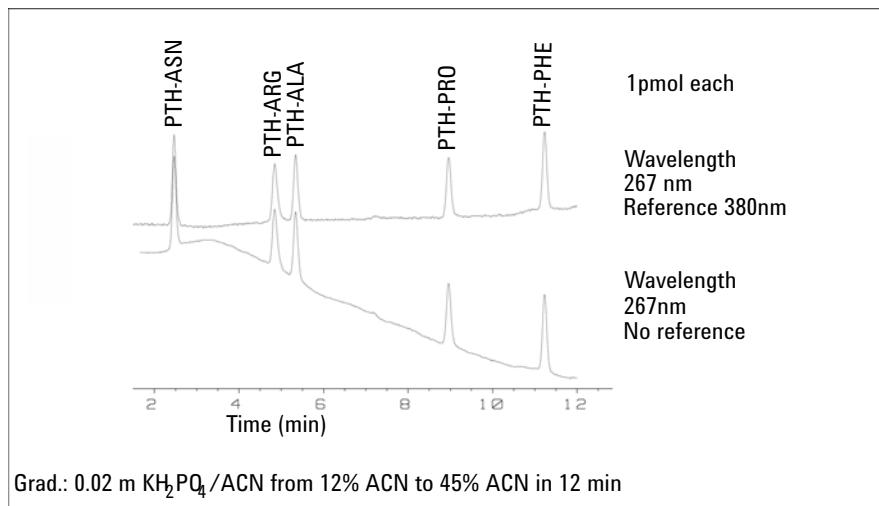


Figure 40 Gradient Analysis of PTH-Amino Acids (1 pmol each), with and without Reference

Slit Width

The detector has a variable slit at the entrance of the spectrograph. This is an effective tool to adapt the detector to changing demand of different analytical problems.

A narrow slit provides spectral resolution for analytes with very fine structures in the absorbance spectrum. An example of such a spectrum is benzene. The five main absorbance bands (fingers) are only 2.5 nm wide and just 6 nm apart from each other.

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

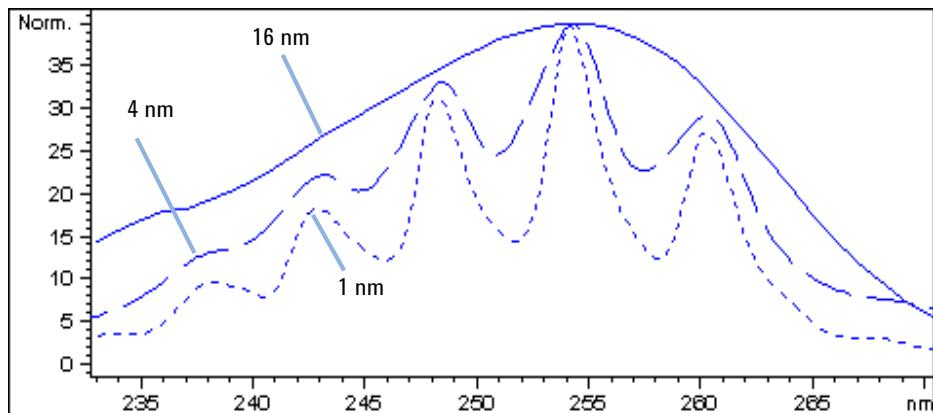


Figure 41 Benzene at 1, 4 and 16 nm slit width (principle)

A wide slit uses more of the light shining through the flow cell. This gives lower baseline noise as shown in **Figure 42** on page 128.

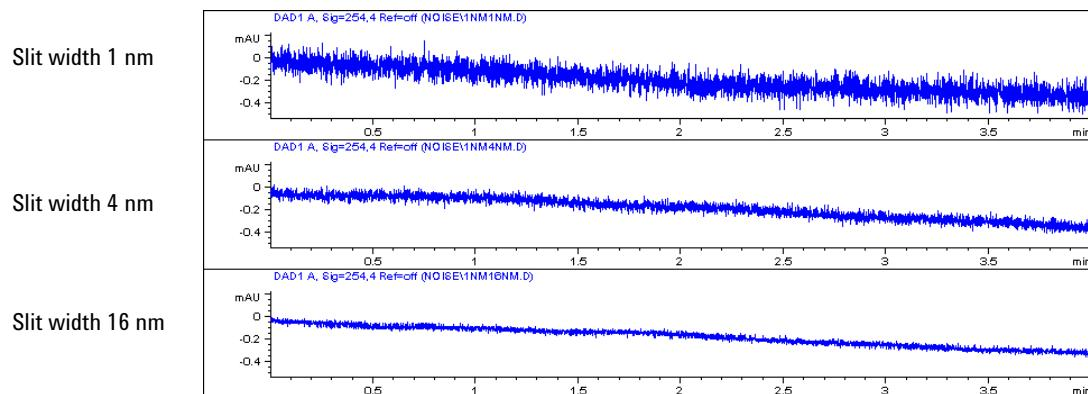


Figure 42 Influence of the Slit Width on Baseline Noise

However, with a wider slit, the spectrograph's optical resolution (its ability to distinguish between different wavelengths) diminishes. Any photodiode receives light within a range of wavelength determined by the slit width. This explains why the fine spectral structure of benzene disappears when using a 16-nm wide slit.

Furthermore, the absorbance is no longer strictly linear with concentration for wavelengths at a steep slope of a compound's spectrum.

Substances with fine structures and steep slopes like benzene are very rare.

In most cases the width of absorbance bands in the spectrum is more like 30 nm as with anisic acid ([Figure 38](#) on page 125).

In most situations, a slit width of 4 nm will give the best results.

Use a narrow slit (1 or 2 nm) if you want to identify compounds with fine spectral structures or if you need to quantify at high concentrations (> 1000 mAU) with a wavelength at the slope of the spectrum. Signals with a wide bandwidth can be used to reduce baseline noise. Because (digital) bandwidth is computed as average of absorbance, there is no impact on linearity.

Use a wide (8 or 16 nm) slit when your sample contains very small concentrations. Always use signals with bandwidth at least as wide as the slit width.

Optimizing Spectral Acquisition (DAD only)

Storage of all spectra consumes a lot of disk space. It is very useful to have all spectra available during optimization of a method or when analyzing unique samples. However when running many samples of the same type, the large size of data files with all spectra may become a burden. The detector provides functions to reduce the amount of data, yet retaining the relevant spectral information.

For spectra options see [Table 23](#) on page 134.

Range

Only the wavelength range where the compounds in your sample absorb contains information that is useful for purity checks and library searches. Reducing the spectrum storage range saves disk space.

Step

Most substances have broad absorbance bands. Display of spectra, peak purity and library search works best if a spectrum contains 5 to 10 data points per width of the absorbance bands. For anisic acid (the example used before) a

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

step of 4 nm would be sufficient. However a step of 2 nm gives a more optimal display of the spectrum.

Threshold

Sets the peak detector. Only spectra from peaks higher than threshold will be stored when a peak-controlled storage mode is selected.

Margin for Negative Absorbance

The detector adjusts its gain during *balance* such that the baseline may drift slightly negative (about -100 mAU). In some special case, for example, when gradient with absorbing solvents are used, the baseline may drift to more negative values.

Only for such cases, increase the margin for negative absorbance to avoid overflow of the analog-to-digital converter.

Optimizing Selectivity

Quantifying Coeluting Peaks by Peak Suppression

In chromatography, two compounds may often elute together. A conventional dual-signal detector can only detect and quantify both compounds independently from each other if their spectra do not overlap. However, in most cases this is highly unlikely.

With a dual-channel detector based on diode-array technology, quantifying two compounds is possible even when both compounds absorb over the whole wavelength range. The procedure is called peak suppression or signal subtraction. As an example, the analysis of hydrochlorothiazide in the presence of caffeine is described. If hydrochlorothiazide is analyzed in biological samples, there is always a risk that caffeine is present which might interfere chromatographically with hydrochlorothiazide. As the spectra in [Figure 43](#) on page 131 shows, hydrochlorothiazide is best detected at 222 nm, where caffeine also shows significant absorbance. It would therefore be impossible, with a conventional variable wavelength detector, to detect hydrochlorothiazide quantitatively when caffeine is present.

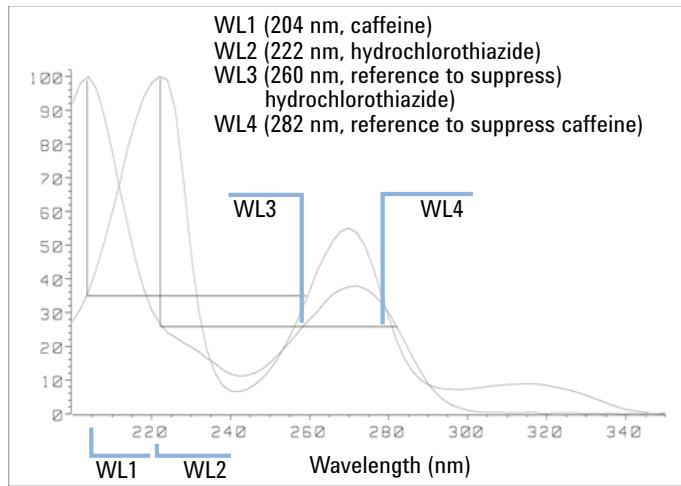


Figure 43 Wavelength Selection for Peak Suppression

With a UV-visible detector based on a diode array and the correct choice of a reference wavelength setting, quantitative detection is possible. To suppress caffeine, the reference wavelength must be set to 282 nm. At this wavelength, caffeine shows exactly the same absorbance as at 222 nm. When the absorbance values are subtracted from each other, any indication of the presence of caffeine is eliminated. In the same way, hydrochlorothiazide can be suppressed if caffeine is to be quantified. In this case the wavelength is set to 204 nm and the reference wavelength to 260 nm. [Figure 44](#) on page 132 shows the chromatographic results of the peak suppression technique.

The trade-off for this procedure is a loss in sensitivity. The sample signal decreases by the absorbance at the reference wavelength relative to the signal wavelength. Sensitivity may be decreased by as much as 10–30 %.

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

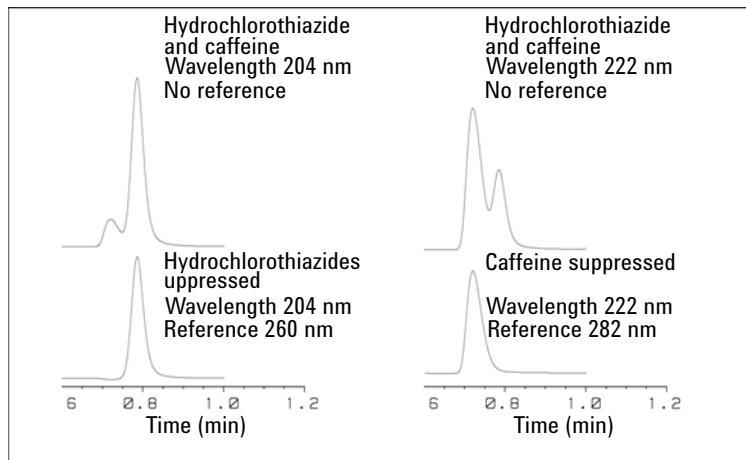
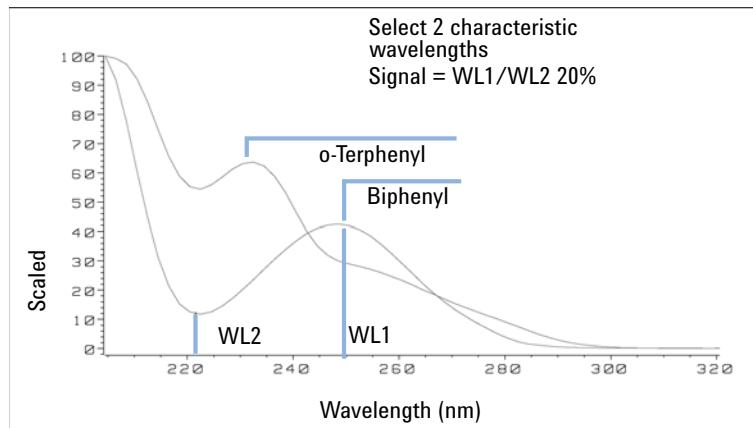
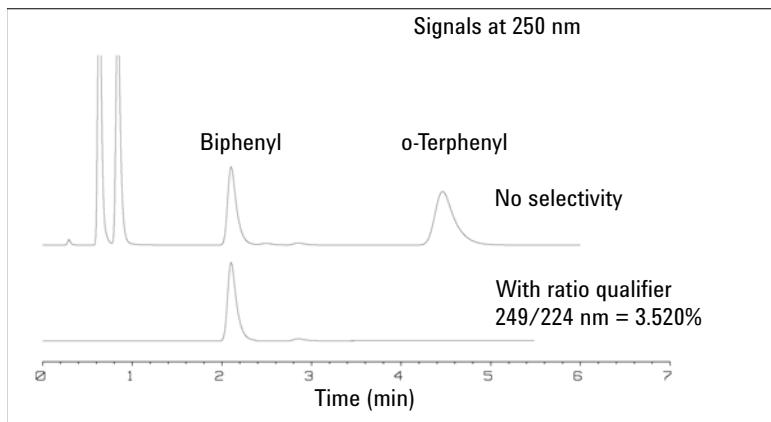


Figure 44 Peak Suppression Using Reference Wavelength

Ratio Qualifiers for Selective Detection of Compound Classes

Ratio qualifiers can be used where, in a complex sample, only one particular class needs to be analyzed – a parent drug and its metabolites in a biological sample, for example. Another example is the selective analysis of derivatives after pre- or post-column derivatization. Specifying a signal ratio that is typical for the sample class is one way of selectively plotting only those peaks that are of interest. The signal output remains at zero so long as the ratio is out of the user-specified ratio range. When the ratio falls within the range, the signal output corresponds to the normal absorbance, giving single, clear peaks on a flat baseline. An example is shown in [Figure 45](#) on page 133 and [Figure 46](#) on page 133.

**Figure 45** Wavelength Selection for Ratio Qualifiers**Figure 46** Selectivity by Ratio Qualifiers

In a four-component mixture, only biphenyl was recorded. The other three peaks were suppressed because they did not meet the ratio-qualifier criterion and therefore the output was set to zero. The characteristic wavelengths 249 nm (λ_1) and 224 nm (λ_2) were found from the spectra shown in [Figure 45](#) on page 133. The ratio range was set at 2 – 2.4 (2.2 ± 10%). Only when the ratio between 249 and 224 nm was within this range, is the signal plotted. Of all four peaks, only the third fulfilled the criterion ([Figure 46](#) on page 133). The others were not plotted.

8 Detector Description

Agilent 1220 Infinity II LC Diode Array Detector (DAD)

Spectrum Settings (DAD only)

To change the Spectra settings open.

- 1 To change the Spectra settings select **Setup Detector Signals**.
- 2 In the section Spectrum click on the drop-down list and chose a parameter. [Table 23](#) on page 134 shows the possible parameters.
- 3 Change the Range, Step width and Threshold according to your needs.

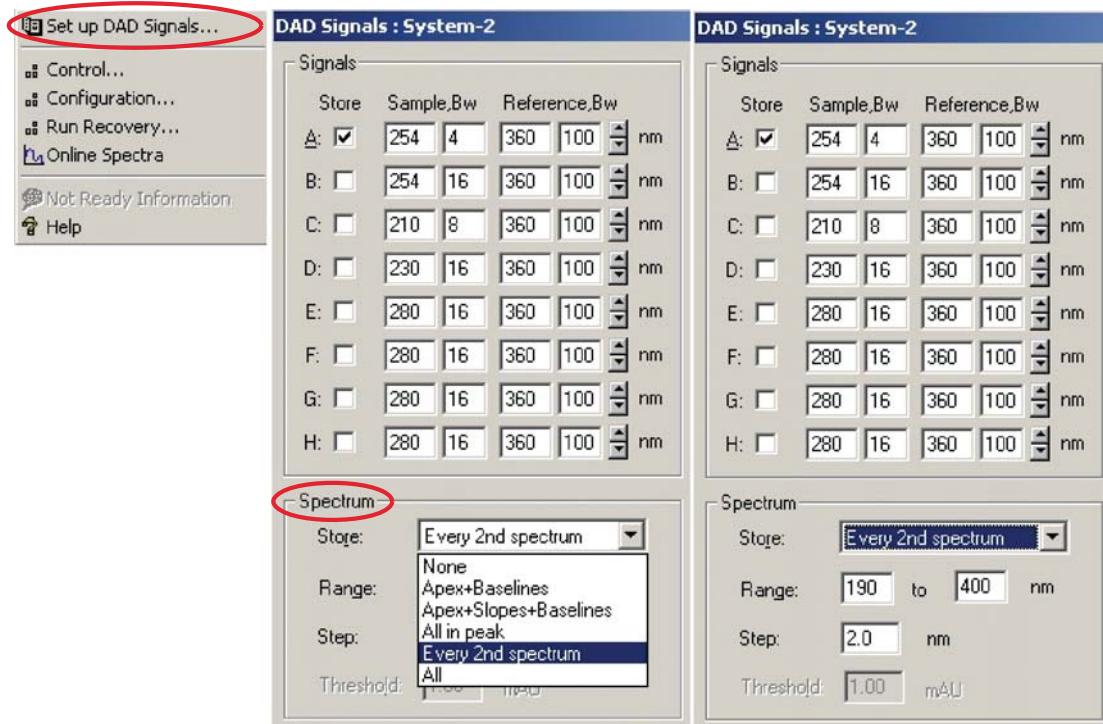


Figure 47 Spectra Settings

Table 23 Spectrum Settings

Store	Defines at which points on "signal A" spectra will be taken and saved. Signal A is used to control the "peak controlled spectra acquisition"; the other signals have no influence on spectra acquisition.
None	No spectra are taken.

Table 23 Spectrum Settings

Apex + Baselines	Spectra are taken at the apex and baselines of the peak.
Apex + Slopes + Baselines	Spectra are taken at the apex, baselines, upslope, and downslope of the peak.
All in Peak	All spectra within the peak are taken.
NOTE	The three spectra acquisition types mentioned above are also referred to as peak-controlled spectra acquisition. The peak detection is done by the detector firmware based on the threshold and peakwidth parameters you set for the DAD. If you want to use peak-controlled spectra storage, make sure that you set these parameters to recognize all the peaks of interest. The integration algorithm also includes peak detection based on the threshold and peakwidth parameters set in the integration events.
Every 2nd spectrum	Spectra are taken continuously as for All, but only every second spectrum is stored; other spectra are discarded. This reduces the amount of data storage necessary.
All	Spectra are taken continuously depending on the setting of the Peakwidth. Eight spectra are acquired per Peakwidth. The acquisition time for one spectrum is slightly less than the Peakwidth divided by 8, that is, greater than or equal to 0.01s and less than or equal to 2.55s.
NOTE	If there are no peaks in Signal A, there are no spectra. You cannot process spectra present in other signals.
Range	Range defines the wavelength range for spectral storage. Limits: 190 to 950 nm in steps of 1 nm for both low and high values. The high value must be greater than the low value by at least 2 nm.
Step	Step defines the wavelength resolution for spectral storage. Limits: 0.10 to 100.00 nm in steps of 0.1 nm.
Threshold	The threshold is the height in mAU of the smallest expected peak. The peak detector ignores any peaks which are lower than the threshold value and does not save spectra. Limits: 0.001 to 1000.00 mAU in steps of 0.001 mAU. Usable for modes Apex + Baselines, Apex + Slopes + Baselines and All in Peak

8 Detector Description

Match the Flow Cell to the Column

Match the Flow Cell to the Column

Figure 48 on page 136 shows recommendations for flow cells that match the column used. If more than one selection is appropriate, use the larger flow cell to get the best detection limit. Use the smaller flow cell for best peak resolution.

Choosing a Flow Cell for the VWD

Column length	Typical peak width	Recommended flow cell			
<= 5 cm	0.025 min				
10 cm	0.05 min		Semi-micro flow cell		
20 cm	0.1 min			Standard flow cell	
>= 40 cm	0.2 min				
	Typical flow rate	0.2 ml/min	0.2 - 0.4 ml/min	0.4 - 0.8 ml/min	1 - 5 ml/min
	Internal column diameter	1.0 mm	2.1 mm	3.0 mm	4.6 mm

Figure 48 Choosing a flow cell

Choosing a Flow Cell for the DAD

Typical column length	Typical peak width	Recommended flow cell					
T <= 5 cm	0.025 min	Micro or Semi-nano					
10 cm	0.05 min		Semi-micro flow cell				
20 cm	0.1 min			Standard flow cell			
>= 40 cm	0.2 min						
	Typical flow rate	0.01 ... 0.2 ml/min	0.2 ... 0.4 ml/min	0.4 ... 0.4 ml/min	1 ... 5 ml/min	0.01 ... 5 ml/min	
	Internal column diameter	0.5 ... 1 mm	2.1 mm	3.0 mm	4.6 mm		

Figure 49 Choosing a Flow Cell in HPLC

Flow Cell Path Length

Lambert-Beer's law shows a linear relationship between the flow cell path length and absorbance.

$$\text{Absorbance} = -\log T = \log \frac{I_0}{I} = \epsilon \cdot C \cdot d$$

where

- T** is the transmission, defined as the quotient of the intensity of the transmitted light I divided by the intensity of the incident light, I_0 ,
- e** is the extinction coefficient, which is a characteristic of a given substance under a precisely-defined set of conditions of wavelength, solvent, temperature and other parameters,
- C** is the concentration of the absorbing species (usually in g/l or mg/l), and
- d** is the path length of the cell used for the measurement.

8 Detector Description

Match the Flow Cell to the Column

Therefore, flow cells with longer path lengths yield higher signals. Although noise usually increases a little with increasing path length, there is a gain in signal-to-noise ratio. For example, the noise increases by less than 10 %, but a 70 % increase in signal intensity is achieved by increasing the path length from 6 mm to 10 mm.

When increasing the path length, the cell volume usually increases – in our example, from 5 μL to 13 μL . Typically, this causes more peak dispersion. As [Figure 50](#) on page 138 demonstrates, this does not affect the resolution in the gradient separation in our example.

As a rule-of-thumb, the flow cell volume should be about 1/3 of the peak volume at half height. To determine the volume of your peaks, take the peak width as reported in the integration results, multiply it by the flow rate and divide it by 3.

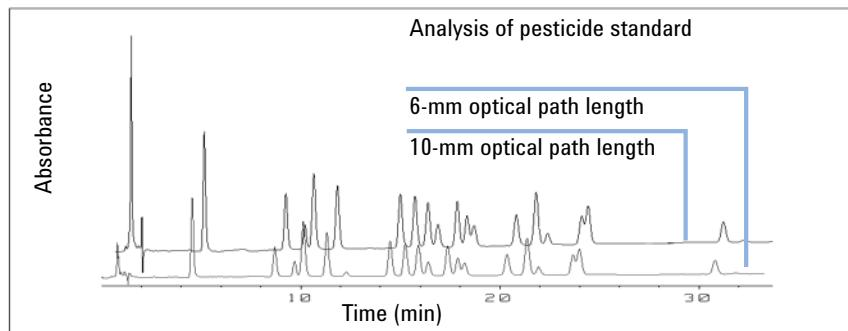


Figure 50 Influence of cell path length on signal height

Traditionally, LC analysis with UV detectors is based on comparing measurements with internal or external standards. To check photometric accuracy of the detector, it is necessary to have more precise information on path lengths of the flow cells.

The correct response is:

expected response * correction factor

Details of the flow cells are shown in [Table 24](#) on page 139 and [Table 25](#) on page 140.

Correction factors for VWD flow cells

Table 24 Correction factors for Agilent VWD flow cells

Flow cell type	Cell volume	Part number	Path length (nominal)	Path length (actual)	Correction factor
Standard flow cell	14 µL	Standard flow cell, 10 mm, 14 µL, 40 bar (G1314-60086)	10 mm	10.15 ± 0.19 mm	10/10.15
Semi-micro flow cell	5 µL	Semi-micro flow cell, 6 mm, 5 µL, 40 bar (G1314-60083)	6 mm	6.10 ± 0.19 mm	6/6.10
Micro flow cell	2 µL	Micro flow cell, without I.D. tag, 3 mm, 2 µL, 120 bar (G1314-60087)	3 mm	2.80 ± 0.19 mm	3/2.8
High Pressure flow cell	14 µL	High pressure flow cell, 10 mm, 14 µL, 400 bar (G1314-60082)	10 mm	10.00 ± 0.19 mm	6/5.75

NOTE

Be aware that there are additional tolerance of gasket thickness and its compression ratio, which are considered to be very small in comparison with the machining tolerance.

8 Detector Description

Match the Flow Cell to the Column

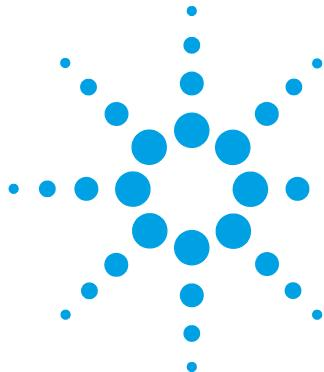
Correction factors for DAD flow cells

Table 25 Correction factors for flow cells

Flow cell	Path length (actual)	Correction factor
Standard flow cell, 10 mm, 13 µL, 120 bar (12 MPa) (G1315-60022)	9.80±0.07 mm	10/9.8
Semi-micro flow cell, 6 mm, 5 µL, 120 bar (12 MPa) (G1315-60025)	5.80±0.07 mm	6/5.8
Micro flow cell, 3 mm, 2 µL, 120 bar (12 MPa) (G1315-60024)	3.00+0.05 mm/-0.07 mm	3/3
Semi-nano flow cell kit, 10 mm, 500 nL, 5 MPa (G1315-68724)	10.00±0.02 mm	10/10
Nano flow cell kit, 6 mm, 80 nL, 5 MPa (G1315-68716)	6.00±0.02 mm	6/6
Standard flow cell bio-inert, 10 mm, 13 µL, 120 bar (12 MPa) for MWD/DAD, includes Capillary Kit Flow Cells BIO (p/n G5615-68755) (G5615-60022)	9.80±0.07 mm	10/9.8

NOTE

Be aware that there are additional tolerance of gasket thickness and its compression ratio, which are considered to be very small in comparison with the machining tolerance.



9

Test Functions and Calibration

1220 Infinity II LC System	143
Installation Check	143
Module Info	144
Diagnostic Buffers	144
State Info	144
Solvent Delivery System	145
Pump Leak Test Description	145
Running the Leak Test	147
Evaluating the Leak Test Results	148
Pressure Too High Check	151
Pressure Too High Check Evaluation	151
Purge Pump	152
Degasser Exchange	153
Autosampler	154
Maintenance Positions	154
Injector Steps	155
Alignment Teaching	156
Gripper Verification	159
Column Oven	160
Oven Test	160
Oven Calibration	161
Variable Wavelength Detector (VWD)	162
Cell Test	162
Dark Current Test	163
Holmium Oxide Test	164
Intensity Test	165
Filter/Grating Motor Test	167
Detector Calibration	168



9 Test Functions and Calibration

Match the Flow Cell to the Column

Test Chromatogram	169
Spectral Scan	170
Diode Array Detector (DAD)	171
Self-test	171
Filter Test	172
Slit Test	174
Dark-Current Test	175
Intensity Test	177
Holmium Oxide Test	180
Spectral flatness test	182
ASTM Noise Test	184
Cell Test	184
Using the Built-in Test Chromatogram	186
Wavelength Verification and Recalibration	188
Diagnosis Information on Agilent ChemStation	189
D/A Converter (DAC) Test	191

This chapter describes the tests, calibrations and tools that are available with the Instrument Utilities software or the Lab Advisor (Service and Diagnostics Section).

1220 Infinity II LC System

Installation Check

The **Installation Check** switches on all available modules, purges the system for five minutes at 1 mL/min, tests the flow path by applying a pressure up to 200 bar and switches on the oven (if available) and detector.

NOTE

The pump and detector are mandatory for this check; the oven and autosampler are optional.

The **Installation Check** passes if the following conditions are met:

- All modules switch on successfully within the timeout period (120 s).
- The pump achieves 200 bar after 5 min.
- The oven reaches 2 K above its actual temperature.
- The detector lamp ignites and the detector reaches a **Ready** state.

Module Info

The **Module Info** tool collects diagnostic information from a module and writes the results to a file. You can view the results in three tabs:

General

The **General** tab shows information about the module's firmware and options in a two-column table.

Tables

The **Tables** tab allows you to display the contents of all available diagnostic tables for the module. You click the **[+]** sign to open a table, or the **[-]** sign to close an open table.

Signals

The **Signals** tab shows the plots of the available diagnostic signals from the module. The signal plots that are available are module-dependent; where available, both short-term and long-term plots are displayed for a signal.

Diagnostic Buffers

Each module contains a set of diagnostic buffers that can be queried.

Lab Advisor allows you to query all diagnostic buffers for a selected module and write the results to a text file. The text file is written in English.

State Info

The **State Info** tool displays the current status of all Agilent 1220 Infinity II LC modules; the status is continuously updated. Unless aborted, the tool runs continuously for 60 min. Click **Stop Test** to stop the tool.

Solvent Delivery System

Pump Leak Test Description

The leak test is a built-in troubleshooting test designed to verify the tightness of pump components. The test should be used when problems with the pump are suspected. The test involves monitoring the pressure increase at very low flow rates while different plungers are delivering solvent. At these very low flow rates, very small leaks can be detected by evaluating the pressure profile as the pump runs through a predefined pumping sequence. The test requires blocking the pump with a blank nut, then running the test with isopropanol (IPA), while monitoring the pressure profile.

NOTE

Make absolutely sure that all parts of the flow path that are included in the test are very thoroughly flushed with IPA before starting to pressurize the system! Any trace of other solvents or the smallest air bubble inside the flow path will definitely cause the test to fail!

Ramp 1

After initialization, plunger 2 is at the top of its stroke. The test begins with plunger 1 delivering with a stroke length of 100 μL and a flow of 153 $\mu\text{L}/\text{min}$. The plunger sequence during the pressure ramp is 1-2-1-2. The pressure increase during this phase should be linear. Pressure disturbances during this phase indicate larger leaks or defective pump components.

Plateau 1

Plunger 2 continues to pump with a flow rate of 2 $\mu\text{L}/\text{min}$ for approximately one minute. The pressure during the plateau should remain constant or increase slightly. A falling pressure indicates a leak of $>2 \mu\text{L}/\text{min}$.

Ramp 2

The flow is changed to 153 $\mu\text{L}/\text{min}$, and plunger 2 continues to deliver for the rest of its stroke. Then plunger 1 continues to pump to complete the second half of the ramp.

9 Test Functions and Calibration

Solvent Delivery System

Plateau 2

The flow is reduced to 2 µL/min for approximately one minute (plunger 1 still delivering). The pressure during the plateau should remain constant or increase slightly. A falling pressure indicates a leak of >2 µL/min.

Ramp 3

The flow increases to 220 µL/min and the stroke is changed to 100 µL. Plunger 1 completes its stroke. Next, the flow is changed to 510 µL/min. The ramp reaches 390 bar with the plunger sequence 2-1-2-1.

Plateau 3

When the system pressure reaches 390 bar, the flow is reduced to zero, and the pressure stabilizes just below 400 bar.

One minute after reaching the maximum pressure, the pressure drop should not exceed 2 bar/min.

Running the Leak Test

When If problems with the pump are suspected

Tools required **Description**
Wrench, 1/4 inch (supplied in HPLC Tool-Kit)

Parts required # p/n Description
1 G1313-87305 Restriction Capillary
1 01080-83202 Blank nut
1 500 mL Isopropanol

Preparations • Place a bottle of LC-grade isopropyl alcohol in the solvent cabinet and connect its solvent tube to the active inlet valve of the pump.

NOTE Make absolutely sure that all parts of the flow path that are part of the test are very thoroughly flushed with IPA before starting to pressurize the system. Any trace of other solvents, or the smallest air bubble inside the flow path, will definitely cause the test to fail.

Running the test from the Agilent Lab Advisor

- 1 Select the **Leak Test** from the **Test Selection** menu.
- 2 Start the test and follow the instructions.

NOTE Make sure to release the pressure by slowly opening the purge valve when the test has finished.

NOTE Evaluating the Results describes the evaluation and interpretation of the leak test results ([“Evaluating the Leak Test Results”](#) on page 148).

NOTE For detailed instructions refer to the Agilent Lab Advisor tool.

9 Test Functions and Calibration

Solvent Delivery System

Evaluating the Leak Test Results

Defective or leaky components in the pump head lead to changes in the leak-test pressure plot. Typical failure modes are described below.

NOTE

Please notice the difference between an **error** in the test and a **failure** of the test! An **error** means that during the operation of the test there was an abnormal termination. If a test **failed**, this means that the results of the test were not within the specified limits.

NOTE

Often it is only a damaged blank nut itself (poorly shaped from overtightening) that causes a failure of the test. Before investigating on any other possible sources of failure make sure that the blank nut you are using is in good condition and properly tightened!

Table 26 No Pressure increase at Ramp 1

Potential Cause	Corrective Action
Pump not running.	Check the logbook for error messages.
Purge valve open.	Close the purge valve, and restart the test.
Loose or leaky fittings.	Ensure all fittings are tight, or exchange capillary.
Wrong solvent-line connections.	Ensure the solvent lines from the degasser are connected correctly.
Contaminated purge valve.	Open and close purge valve to flush out contamination. Exchange the valve if still leaky.
Large leaks (visible) at the pump seals.	Exchange the pump seals.
Large leaks (visible) at active inlet valve, outlet valve, or purge valve.	Ensure the leaky components are installed tightly. Exchange the component if required.

Table 27 Pressure limit not reached but plateaus horizontal or positive

Potential Cause	Corrective Action
Degasser and pump not flushed sufficiently (air in the pump head).	Purge the degasser and pump thoroughly with isopropanol under pressure (use the restriction capillary).
Wrong solvent.	Install isopropanol. Purge the degasser and pump thoroughly.

Table 28 All plateaus negative

Potential Cause	Corrective Action
Loose or leaky fittings.	Ensure all fittings are tight, or exchange capillary.
Loose purge valve.	Tighten the purge valve (14 mm wrench).
Contaminated purge valve.	Open and close purge valve to flush out contamination. Exchange the valve if still leaky.
Loose pump head screws.	Ensure the pump head screws are tight.
Leaking seals or scratched plungers.	Exchange the pump seals. Check the plungers for scratches. Exchange if scratched.
Leaking outlet valve.	Exchange the outlet valve.
Leaky damper.	Exchange damper.

Table 29 First plateau positive, second and third plateau negative

Potential Cause	Corrective Action
Air in pump or new seals not yet seated.	Flush pump thoroughly with isopropanol under pressure (use restriction capillary).
Loose active inlet valve.	Tighten the active inlet valve (14 mm wrench). Do not overtighten!
Loose pump head screws.	Ensure the pump head screws are tight.
Loose outlet valve.	Ensure the sieve in the outlet valve is installed correctly. Tighten the outlet valve.

9 Test Functions and Calibration

Solvent Delivery System

Table 29 First plateau positive, second and third plateau negative

Leaking seal or scratched plunger.	Exchange the pump seals. Check the plungers for scratches. Exchange if scratched.
Defective active inlet valve.	Exchange the active inlet valve.

Table 30 First plateau negative, second plateau positive

Potential Cause	Corrective Action
Leaking outlet valve.	Clean the outlet valve. Ensure the sieve in the outlet valves are installed correctly. Tighten the outlet valve.
Loose pump head screws.	Ensure the pump head screws are tight.
Leaking seals or scratched plungers.	Exchange the pump seals. Check the plunger for scratches. Exchange if scratched.

Table 31 Ramp 3 does not reach limit

Potential Cause	Corrective Action
Pump stopped due to error.	Check the logbook for error messages.
Large leaks (visible) at the pump seals.	Exchange the pump seals.
Large leaks (visible) at active inlet valve, outlet valve, or purge valve.	Ensure the leaky components are installed tightly. Exchange the component if required.

Table 32 Third plateau negative (pressure drop > 2 bar/min)

Potential Cause	Corrective Action
Loose or leaky fittings.	Ensure all fittings are tight, or exchange capillary.
Loose purge valve.	Tighten the purge valve (14 mm wrench).
Contaminated purge valve.	Open and close purge valve to flush out contamination. Exchange the valve if still leaky.
Loose pump head screws.	Ensure the pump head screws are tight.

Table 32 Third plateau negative (pressure drop > 2 bar/min)

Leaking seals or scratched plungers.	Exchange the pump seals. Check the plungers for scratches. Exchange if scratched.
Leaking outlet valve.	Exchange the outlet valve.
Leaky damper.	Exchange damper.

Pressure Too High Check

This test checks the flow path for a blockage, and tries to identify the module that is causing the blockage. If the blockage is in the autosampler, the test can identify whether the blockage occurs in the needle or needle seat.

The pump and autosampler are necessary to run the **Pressure Too High Check**.

Pressure Too High Check Evaluation

Start Conditions

The pump and autosampler are brought to the READY state, and an operating pressure of 200 bar is applied to the system.

The pump ripple is measured, and the start of the test is delayed until the ripple is within the defined limits (typically 1 % of operating pressure).

Test Part 1

Part 1 of the test tries to determine in which part of the system the pressure problem lies.

After the system has achieved the start conditions, the autosampler valve is switched from mainpass to bypass, and the pressure slice is tested against a limit.

If the limit is exceeded, the pressure problem lies in the autosampler; otherwise, it lies somewhere in the rest of the flow path.

9 Test Functions and Calibration

Solvent Delivery System

Test Part 2

In Part 2 of the test, an empty vial is driven to the needle, and the valve is switched from bypass to mainpass, so that the pressure should drop dramatically. The pressure drop is checked against a limit.

If the autosampler is identified as the source of the problem, and the limit is not reached, the problem lies in the needle, needle loop or metering drive; otherwise, the problem lies in the needle seat or needle seat capillary.

If the autosampler is not the source of the problem, the problem lies either in the pump (typically the filter or frit), or after the autosampler (heater capillary or column). If the pressure drop limit is not achieved, the problem lies in the pump; otherwise, the problem occurs after the autosampler.

Purge Pump

Purge Pump Description

The **Purge Pump** tool enables you to purge the pump with solvent at a specified flow rate for a specified time. For multi-channel pumps, and pumps with solvent selection valve (SSV), you select the channels to purge; each channel can be purged with different conditions.

You can select a flow rate between 1 and 5 mL/min in steps of 1 mL/min.

You can select a time from 1, 2, 3, 5, 7, 10 and 15 minutes.

NOTE

The G1361A Prep Pump has an automatic purge cycle; there are no user-configurable options.

Purging the Pump

To purge the pump

- 1 Prepare each channel with the appropriate purge solvents.
- 2 Select **Purge Pump** from the **Tool Selection** screen.
- 3 In the **Purge Configuration** dialog box,
 - a If necessary, select the channel(s) that you want to purge.

- b For each selected channel, select a **Flow** and a purge **Time**.
 - c Click **OK** to close the **Purge Configuration** dialog box.
 - 4 When the request to open the purge valve appears, open the purge valve on the pump, then click **OK** to close the message box.

During purging, the **General** tab shows the current channel that is being purged, and the remaining purge time. The **Signals** tab shows a plot of pressure against time for the complete purge cycle.

 - 5 When the purge time has elapsed and the request to close the purge valve appears, close the purge valve on the pump, then click **OK** to close the message box.

The pump purge process is complete.

Degasser Exchange

The **Degasser Exchange Tool** allows you to set a new calibration value after exchanging the degasser unit.

When you start the **Degasser Exchange Tool**, a dialog box is displayed asking you to enter the calibration value written on the degasser label. Type the calibration value in the field and click **OK** to upload it to the degasser.

NOTE

With the introduction of Lab Advisor B.02.04[093], the **Degasser Exchange Tool** function has been moved to **Instrument Control** tab.

To use the tool, do as follows:

- 1 Go to **Instrument Control** tab
- 2 Select **G4281B** pump module
- 3 Expand the pump function by clicking the small right arrow
- 4 In **special commands** you will find the tool

9 Test Functions and Calibration

Autosampler

Autosampler

Maintenance Positions

Change Needle

The **Change Needle** function moves the safety flap out of position, and positions the needle for easy exchange and alignment.

Start	moves the safety flap away from the needle, and positions the needle approximately 15 mm above the needle seat.
Up	moves the needle arm up stepwise.
Down	moves the needle arm down stepwise. The lowest position is used to align the needle at the correct position in the needle arm.
End	repositions the safety flap around the needle.

Change Piston

The **Change Piston** function draws the piston away from the home position, relieving the tension on the spring. In this position, the analytical head assembly can be removed and reinstalled easily after maintenance.

Start	draws the piston away from the home position, relieving the tension on the spring.
End	repositions the piston at the home position.

Change Gripper

The **Change Gripper** function moves the gripper to the front of the autosampler enabling easy access to the gripper release mechanism.

Start	moves the gripper to the front of the sample-tray area.
End	repositions the gripper at the home position.

Arm Position

Move Arm Home	Moves the gripper arm to its home position for better access and exchange of trays.
Park Arm	Secures the gripper arm to the park position behind the sampling unit. Before parking the gripper arm, ensure there is no vial in the gripper.

Injector Steps

Each movement of the sampling sequence can be done under manual control. This is useful during troubleshooting, where close observation of each of the sampling steps is required to confirm a specific failure mode or verify successful completion of a repair.

Each injector step command actually consists of a series of individual commands that move the autosampler components to predefined positions, enabling the specific step to be done.

Table 33 Injector step commands

Step	Action	Comments
Valve Bypass	Switches injection valve to the bypass position.	
Piston Home	Moves the piston to the home position.	
Needle Up	Lifts the needle arm to the upper position.	Command also switches the valve to bypass if it is not already in that position.
Vial to Seat	Moves the selected vial to the seat position.	Command also lifts the needle to the upper position.

9 Test Functions and Calibration

Autosampler

Table 33 Injector step commands

Step	Action	Comments
Needle into Sample	Lowers the needle into the sample.	Command also positions the vial at the seat, and lifts the needle to the upper position.
Draw	Metering device draws the defined injection volume.	Command also positions the vial at the seat, lifts the needle, and lowers the needle into vial. Command can be done more than once (maximum draw volume of 100 µL cannot be exceeded). Use Plunger Home to reset the metering device.
Needle Up	Lifts the needle out of the vial.	Command also switches the valve to bypass if it is not already in that position.
Vial to Tray	Returns the selected vial to the tray position.	Command also lifts the needle to the upper position.
Needle into Seat	Lowers the needle arm into the seat.	Command also returns the vial to the tray position.
Valve Mainpass	Switches the injection valve to the mainpass position.	
Reset	Resets the injector.	

Alignment Teaching

ALS Alignment Teaching Description

The Alignment Teaching tool is required to compensate for small deviations in positioning of the autosampler gripper that may occur after the module has been disassembled for repair. It requires that a 100-vial tray be inserted into the autosampler.

The alignment procedure uses two tray positions as reference points; because the tray is rectangular, a two-point alignment is sufficient to correct all other

vial positions on the tray. When the correction calculation is complete, the values for both X and theta are rounded to one decimal place. On completion of the alignment procedure, the corrected gripper positions are stored in the module's firmware.

NOTE

To ensure correct operation of the autosampler, the alignment procedure must be carried out in the correct sequence and in full (that is, without skipping any part).

ALS Alignment Controls

Button	Description	Keyboard Shortcut
	Rotate the gripper by increasing theta.	Cursor Up
	Move the gripper horizontally to the left.	Cursor Left
	Move the gripper horizontally to the right.	Cursor Right
	Rotate the gripper by decreasing theta.	Cursor Down
Arm Up	Lifts the gripper arm.	Page Up
Arm Down	Lowers the gripper arm.	Page Down
Open Gripper	Opens the gripper.	
Close Gripper	Closes the gripper.	
Start >>	Starts the execution of the procedure. Shown only at the start.	Enter
Continue >>	Jumps to the next step of the procedure. Shown only during alignment.	Enter
Restart	Restarts the execution of the step.	

9 Test Functions and Calibration

Autosampler

Running the ALS Alignment Teaching

NOTE

To ensure correct operation of the autosampler, the alignment procedure must be carried out in the correct sequence and in full (that is, without skipping any part).

To align the ALS:

- 1 Insert a 100-vial tray into the autosampler.
- 2 Place capped vials into positions 15 and 95.
- 3 Click **Start >>**.

The gripper arm moves to a position above vial 15.

- 4 Answer **Yes** to reset the correction values to their factory defaults, or **No** to leave them as they are.
- 5 Use **Arm Down** to move the fingers of the gripper as close as possible to the top of the vial without touching.
- 6 Use **▲** and **▼** (for rotation), and **◀** and **▶** (for movement left and right) to adjust the gripper position in the horizontal plane.
- 7 Use **Open Gripper** to open the gripper fingers.
- 8 Use **Arm Down** to move the gripper arm down a further 5 mm until the vial cap and the rubber of the gripper fingers are the same height.
- 9 Check that the vial is in the center of the gripper fingers and readjust the position if necessary (step 6).
- 10 When you are satisfied that the gripper position is correct, click **Continue**.
The gripper arm moves to a position above vial 95.

- 11 Repeat steps 6 to 9 to align the gripper at position 95.

- 12 Click **Continue**.

On completion of the calculation, the values for both X and theta are rounded to one decimal place. The correction values are stored permanently in the non-volatile memory of the sampler, and the sampler is initialized.

Gripper Verification

ALS Gripper Verification Description

The verification procedure uses several vial positions as reference points to verify the gripper alignment is correct. If verification indicates one or more positions are out of alignment, the alignment procedure should be done.

Verifying the ALS Gripper Positions

Vial positions 1, 10, 55, 81, 91 (Lab Advisor B.02.07 and later) and 100 can be used for position verification.

- 1 Insert empty capped vials into the vial tray at the positions to be verified.
- 2 Select the first vial position in the vial-position menu.
- 3 Select **Go to selected position**.
- 4 If the fingers of the gripper arm are aligned centrally above the vial, select **Pick vial** to verify the gripper arm lifts the vial out of the tray correctly. If there is a misalignment, the gripper must be realigned.
- 5 Select **Put vial** to verify that the gripper replaces the vial correctly. If there is a misalignment, the gripper must be realigned.
- 6 Repeat the procedure for the next vial position.

9 Test Functions and Calibration

Column Oven

Column Oven

Oven Test

Oven Test Description

The Oven Test is used to evaluate the heating performance of the heating element.

The heating rate over a 10 K interval from the start temperature is determined. The start temperature must be between 30 °C and 50 °C, and is determined as follows:

- If the current oven temperature is below 30 °C, the oven tries to attain a temperature of 30 °C. 30 °C is used as the start temperature.
- If the current oven temperature is above 30 °C but below 50 °C, the current oven temperature is used as the start temperature.
- If the current oven temperature is above 50 °C, an error message is displayed. The oven must then be allowed to cool to below 50 °C before the test can be run.

Oven Test Evaluation

At the end of the **Oven Test**, the slope of the temperature rise of the left and right channels are evaluated. The test passes if both slopes are $\geq 4^{\circ}\text{C}/\text{min}$.

Oven Calibration

Oven Calibration Description

The oven calibration procedure enables the oven temperature to be measured against an external, calibrated measuring device.

Normally, temperature calibration is not required during the lifetime of the instrument; however, in order to comply with local regulatory requirements, the 2-point calibration and verification procedure may be performed.

Running the Oven Calibration

NOTE

For the measurement and calibration procedure, we recommend a measuring device that provides the necessary resolution and precision, for example, Heraeus Quat340 quartz surface-temperature measurement sensor. Contact your local Agilent Technologies support representative for ordering information.

- 1 Install the calibrated temperature measuring device.
- 2 Select the **Oven Calibration** in the user interface.
- 3 Wait for the oven to reach the first set point (40 °C).
- 4 Measure the temperature of the heat exchanger and enter the value in the field.
- 5 Wait for the oven to reach the second set point (50 °C).
- 6 Measure the temperature of the heat exchanger and enter the value in the field.
- 7 Click **OK** to save the calibration values to the oven, or **Cancel** to abort the calibration process.

9 Test Functions and Calibration

Variable Wavelength Detector (VWD)

Variable Wavelength Detector (VWD)

Cell Test

VWD Cell Test Description

The **Cell Test** compares the intensity of the deuterium lamp measured by the sample and reference diodes (unfiltered and not logarithmized) when the grating is in the zero-order position. The resulting intensity ratio (sample:reference) is a measure of the amount of light absorbed by the flow cell. The test can be used to check for dirty or contaminated flow cell windows. When the test is started, the gain is set to -1. To eliminate effects due to absorbing solvents, the test should be done with water in the flow cell.

NOTE

The test should not be performed using the micro flow cell, since the reduction in light intensity will cause the test to fail.

Evaluating the VWD Cell Test Results

The intensity ratio is dependent on the degree of contamination of the flow cell windows, and on the type of flow cell used. The lower the ratio, the more light is absorbed by the flowcell.

Table 34 Probable causes of excessive flowcell absorbance

Cause	Corrective action
Absorbing solvent or air bubble in flow cell.	Ensure the flow cell is filled with water, and free from air bubbles.
Dirty or contaminated flow cell.	Exchange the flow cell windows.

Dark Current Test

VWD Dark Current Test Description

The dark-current test measures the leakage current from the sample and reference circuits. The test is used to check for defective sample or reference diodes or ADC circuits which may cause non-linearity or excessive baseline noise. During the test, the lamp is switched off. Next, the leakage current from both diodes is measured. The test evaluates the results automatically.

Evaluating the VWD Dark Current Test

Table 35 Limits

Max. Sample Intensity	790000 counts
Max. Reference Intensity	790000 counts

Probable causes of test failure

Table 36 Sample circuit noise exceeds limit:

Cause	Corrective action
Defective sample diode	Exchange the sample diode.
Defective sample ADC board.	Exchange the sample ADC board.

Table 37 Reference circuit noise exceeds limit:

Cause	Corrective action
Defective reference diode.	Exchange the reference diode.
Defective reference ADC board.	Exchange the reference ADC board.

9 Test Functions and Calibration

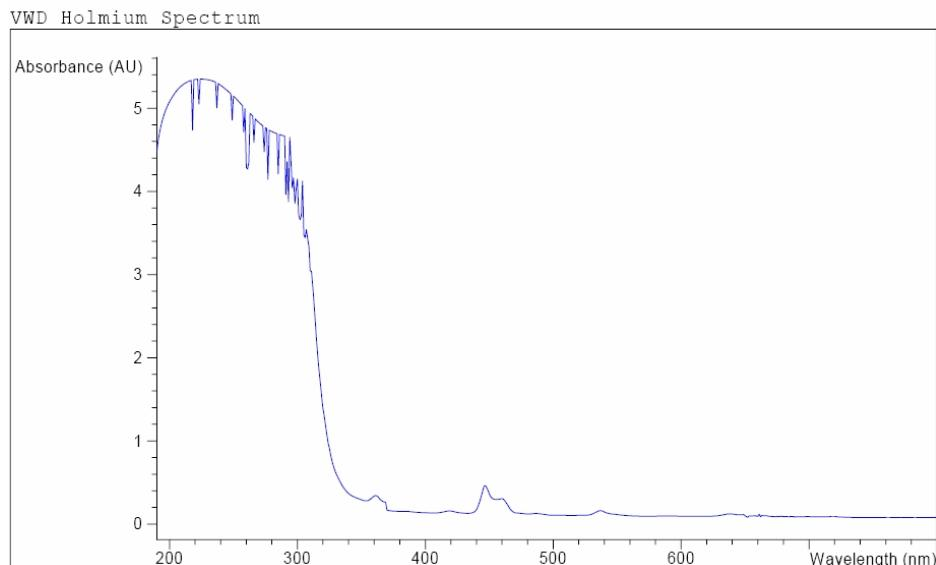
Variable Wavelength Detector (VWD)

Holmium Oxide Test

VWD Holmium Oxide Test Description

The **Holmium Oxide Test** uses three characteristic absorbance maxima of the built-in holmium oxide filter to verify wavelength accuracy (see also **Wavelength Calibration**). The test evaluates the results automatically, and provides a spectrum of the holmium oxide filter. To eliminate effects due to absorbing solvents, the test should be done with water in the flow cell. On completion of the test, the results are displayed automatically.

Holmium Oxide Test Report



VWD Holmium Test Results

	Specification	Measured	Result
<hr/>			
Deviation from wavelength 1:	360.8 nm	-1..1 nm	0.0 nm Passed
Deviation from wavelength 2:	418.5 nm	-1..1 nm	0.1 nm Passed
Deviation from wavelength 3:	536.4 nm	-1..1 nm	0.0 nm Passed

Evaluating the VWD Holmium Oxide Test

The test is evaluated by the instrument, and the measured maxima are displayed automatically. The test fails if one or more of the maxima lies outside the limits.

Table 38 Limits

Absorbance Maxima	Limits
360.8 nm	-1 to +1 nm
418.5 nm	-1 to +1 nm
536.4 nm	-1 to +1 nm

Intensity Test

VWD Intensity Test Description

The **Intensity Test** measures the intensity of the UV lamp over the full VWD wavelength range (190 – 800 nm). The test evaluates the results automatically, and provides an intensity spectrum. The test evaluates the highest intensity, average intensity, and lowest intensity across the full wavelength range. The test is used to determine the performance of the lamp and optics (see also cell test). To eliminate effects due to absorbing solvents, the test should be done with water in the flow cell. The shape of the intensity spectrum is primarily dependent on the lamp and grating. Therefore, intensity spectra will differ slightly between instruments. On completion of the test, the intensity spectrum and intensity values are displayed. (“[VWD Cell Test Description](#)” on page 162)

NOTE

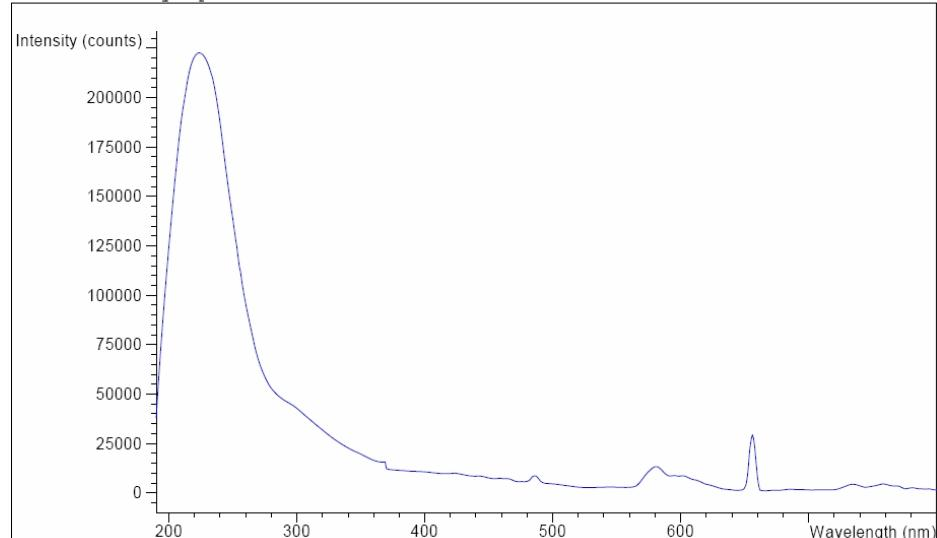
The test should not be performed using the micro flow cell, since the reduction in light intensity will cause the test to fail.

9 Test Functions and Calibration

Variable Wavelength Detector (VWD)

Intensity Test Report

VWD Intensity Spectrum



VWD Intensity Test Results

	Specification	Measured	Result
Accumulated lamp on time		94.35 h	
Highest intensity	> 320000 cts	7123680 cts	Passed
Average intensity	> 160000 cts	951488 cts	Passed
Lowest intensity	> 6400 cts	36384 cts	Passed

Evaluating the VWD Intensity Test

Table 39 Limits

Intensity	Limits (counts)
Highest	>320000
Average	>160000
Lowest	>6400

Table 40 Probable causes of test failure

Cause	Corrective action
Lamp off.	Switch on the lamp.
Old lamp.	Exchange the lamp.
Absorbing solvent or air bubble in flow cell.	Ensure the flow cell is filled with water, and free from air bubbles.
Dirty or contaminated flow cell.	Run the Cell Test. If the test fails, exchange the flow cell windows ("VWD Cell Test Description" on page 162).

Filter/Grating Motor Test

VWD Filter/Grating Test Description

The actual position of the filter motor and grating motor is defined as the number of steps from the reference (sensor) positions. The **VWD Filter/Grating Test** counts the number of motor steps required to move the filter motor and grating motor back to the reference (sensor) position. If the number of steps required to reach the reference positions are the same as the expected step number, the test is passed. If a motor fails to move, or loses motor steps, the test fails. The test evaluates the results automatically.

VWD Filter/Grating Test Results

Probable causes of test failure:

Table 41 Filter Motor Test

Cause	Corrective action
Defective filter motor assembly.	Exchange the filter motor assembly.
Defective VWM board.	Exchange the VWM board.

Table 42 Grating Motor Test

Cause	Corrective action
Defective filter motor assembly.	Exchange the filter motor assembly.
Defective VWM board.	Exchange the VWM board.

9 Test Functions and Calibration

Variable Wavelength Detector (VWD)

Detector Calibration

Wavelength Verification/Calibration

Wavelength Calibration of the detector is done using the zero-order position and 656 nm (alpha-emission line) and beta-emission line at 486 nm emission-line positions of the deuterium lamp. The calibration procedure involves three steps. First the grating is calibrated on the zero-order position. The stepper-motor step position where the zero-order maximum is detected is stored in the detector. Next, the grating is calibrated against the deuterium emission-line at 656 nm, and the motor position at which the maximum occurs is stored in the detector. Finally, the grating is calibrated against the deuterium emission-line at 486 nm, and the motor position at which the maximum occurs is stored in the detector.

NOTE

The wavelength verification/calibration takes about 2.5 min and is disabled within the first 10 min after ignition of the lamp because initial drift may distort the measurement.

When the lamp is turned ON, the 656 nm emission line position of the deuterium lamp is checked automatically.

When to Calibrate the Detector

The detector is calibrated at the factory, and under normal operating conditions should not require recalibration. However, it is advisable to recalibrate:

- after maintenance (flow cell or lamp),
- after repair of components in the optical unit,
- after exchange of the optical unit or VWM board,
- at a regular interval, at least once per year (for example, prior to an Operational Qualification/Performance Verification procedure), and
- when chromatographic results indicate the detector may require recalibration.

Test Chromatogram

A built-in pre-defined test chromatogram on the VWM board is processed through ADC like normal signals from the diodes and can be used to check the ADC and the data handling system. The signal is available at the analog output and on the GPIB.

NOTE

The run time of the chromatogram is depending on the setting for response time (peak width). If no stop time is set the chromatogram will repeat continuously.

Response Time	Stop Time
0.06 sec	0.8 min
0.12 sec	0.8 min
0.25 sec	0.8 min
0.50 sec	0.8 min
1.00 sec	1.6 min
2.00 sec	3.2 min (Default settings)
4.00 sec	6.4 min
8.00 sec	12.8 min

The test chromatogram has four main peaks with the following absorbances:

Peak	Absorbance (approx.)
1	38 mAU
2	100 mAU
3	290 mAU
4	20 mAU

9 Test Functions and Calibration

Variable Wavelength Detector (VWD)

Spectral Scan

The **Spectral Scan** tool is available for diode-array and variable wavelength detectors (DAD/MWD and VWD). It allows you to scan a spectrum over a specified wavelength range and export the data to a csv (comma-separated values) file that can be used in other applications (for example, Microsoft Excel).

Scan Parameters

UV Lamp On	Switches on the UV lamp.
Blank Scan (VWD only)	Scans a blank spectrum (solvent only) over the specified wavelength range at the specified resolution. You specify the wavelength range in the from and to fields, and the resolution in the step field.
Sample Scan	Scans the sample spectrum over the specified wavelength range at the specified resolution. You specify the wavelength range in the from and to fields, and the resolution in the step field.
Export Data	Exports the selected data in csv format for use in other applications.

Diode Array Detector (DAD)

This chapter describes the detector's built in test functions.

Self-test

The DAD self-test (see [Figure 51](#) on page 172) runs a series of individual tests, and evaluates the results automatically. The following tests are run:

- Filter Test
- Slit Test
- Dark Current Test
- Intensity Test
- Wavelength Calibration Test
- Holmium Test
- Spectral Flatness Test
- ASTM Noise Test (optional)

The self-test can be run once or repetitively. When set up to run repetitively, the tests run in series continually until stopped by the user. Running the test repetitively is useful when troubleshooting problems which occur intermittently.

The ASTM noise test determines the detector baseline noise (254 nm) while pumping water at 1 mL/min. The test requires approximately 20 minutes to run, and can be included or excluded from the self-test sequence as required.

Setup of the self test is done in the Self Test dialog box. Select either Single Test or Repetitive Tests. Check the ASTM Noise Test checkbox to include the noise test in the self test.

9 Test Functions and Calibration

Diode Array Detector (DAD)

Agilent G1315 Self Test Results			
	Limits	Measured	Result
Filter test	0.005..0.5 AU	0.27 AU	Passed
Slit test	0.7..1.3	1.08	Passed
Dark current test	0..12000 cts	7977..8026 cts	Passed
Min. intensity (190nm - 220nm)	> 2000 cts	18042 cts	Passed
Min. intensity (221nm - 350nm)	> 5000 cts	28451 cts	Passed
Min. intensity (351nm - 500nm)	> 2000 cts	16886 cts	Passed
Min. intensity (501nm - 950nm)	> 2000 cts	14683 cts	Passed
Max. intensity (190nm - 350nm)	< 450000 cts	89461 cts	Passed
Max. intensity (700nm - 950nm)	< 300000 cts	63609 cts	Passed
Max. intensity (D2 alpha line)	< 1200000 cts	169947 cts	Passed
Wavelength at 486.0nm	485.5..486.5 nm	486.31 nm	Passed
Wavelength at 656.1nm	655.6..656.6 nm	656.25 nm	Passed
Holmium test	-1..1 nm	0.30 nm	Passed
Spectral flatness	< 0.002 AU	0.0000 AU	Passed
ASTM noise (20 min. at 254nm)	<= 0.02 mAU	0.0096 mAU	Passed

Figure 51 Self-test Results (report)

For details refer to the individual tests on the following pages.

Filter Test

The filter test checks the correct operation of the filter assembly. When the test is started, the holmium oxide filter is moved into position. During filter movement, the absorbance signal is monitored. As the edge of the filter passes through the light path, an absorbance maximum is seen. Once the filter is in position, the absorbance maximum (of holmium oxide) is determined. Finally, the filter is moved out of the light path. During movement, an additional absorbance maximum is expected as the edge of the filter passes through the light path. The test passes successfully, if the two maxima resulting from the edge of the filter assembly (during filter movement) are seen, and the absorbance maximum of holmium oxide is within the limits.

```

Instrument: G1315C
Serial Number: PP00000024
Operator: Wolfgang
Date: 25.02.2005
Time: 14:55:42
File: C:\CHEM32\2\DIAGNOSE\DAD_FILTER3.DGR

```

	Result	Status
DAD Filter Test		
Date: 25.02.2005; Time: 14:55:42		
Expected total time: approx. 45s.		
Test Procedure:		
1. Measuring and evaluating filter test result	0.16	done

Figure 52 Filter Test Results (report)

Filter Test Evaluation

Filter Test Failed

Test Failed

Probable cause	Suggested actions
1 Filter assembly (lever and filter) not installed.	Install the filter assembly.
2 Defective filter motor.	Please contact your Agilent service representative.

Holmium Oxide Maximum out of Limits

Probable cause	Suggested actions
1 Holmium oxide filter not installed.	Install the holmium oxide filter.
2 Dirty or contaminated filter.	Exchange the holmium oxide filter.

Slit Test

The slit test verifies correct operation of the micromechanical slit. During the test, the slit is moved through all slit positions while the detector monitors the lamp intensity change. When the slit position is changed, the intensity drop (move to smaller slit) or intensity increase (move to larger slit) must be within a defined range. If the intensity changes are outside the expected range, the test fails.

Slit Test Evaluation

Limit: 0.7 - 1.3 (for standard flow cell)

Test Failed

Probable cause	Suggested actions
1 Flow cell still installed.	Remove the flow cell.
2 Old or non-Agilent lamp.	Run the “ Intensity Test ” on page 177. Exchange the lamp if old or defective.
3 Defective slit assembly (stray light).	Exchange the slit assembly.
4 Defective detector main board.	Exchange the detector main board.
5 Defective PDA/optical unit.	Exchange the optical unit.

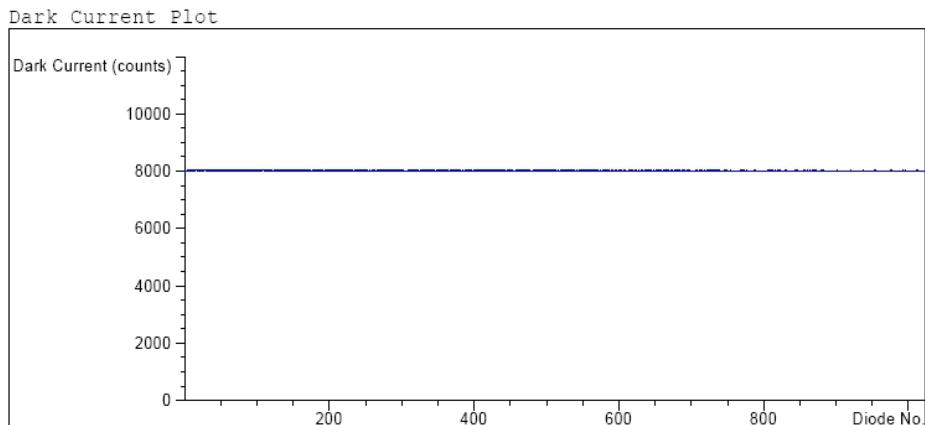
Dark-Current Test

The dark-current test measures the leakage current from each diode. The test is used to check for leaking diodes which may cause non-linearity at specific wavelengths. During the test, the slit assembly moves to the dark position, cutting off all light falling onto the diode array. Next, the leakage current from each diode is measured, and displayed graphically (see [Figure 53](#) on page 176). The leakage current (represented in *counts*) for each diode should fall within the limits (red bands) shown in the plot (see [Figure 53](#) on page 176).

Dark-Current Test Evaluation

Limit: 0... 12000 counts

Instrument: G1315C
Serial Number: FP00000024
Operator: Wolfgang
Date: 25.02.2005
Time: 14:50:12
File: C:\CHEM32\2\DIAGNOSE\DAD_DARK2.DGR



Dark Current Test Results

	Specification	Measured	Result
Dark current maximum value	<= 12000 cts	8056 cts	Passed
Dark current minimum value	> 0 cts	7994 cts	Passed

Figure 53 Dark-Current Test Results (report)

9 Test Functions and Calibration

Diode Array Detector (DAD)

Test Failed

Probable cause	Suggested actions
1 Defective slit assembly (stray light).	Run the “ Slit Test ” on page 174 (part of the “ Self-test ” on page 171).
2 Defective detector main board.	Please contact your Agilent service representative.
3 Defective PDA/optical unit.	Please contact your Agilent service representative.

Intensity Test

NOTE

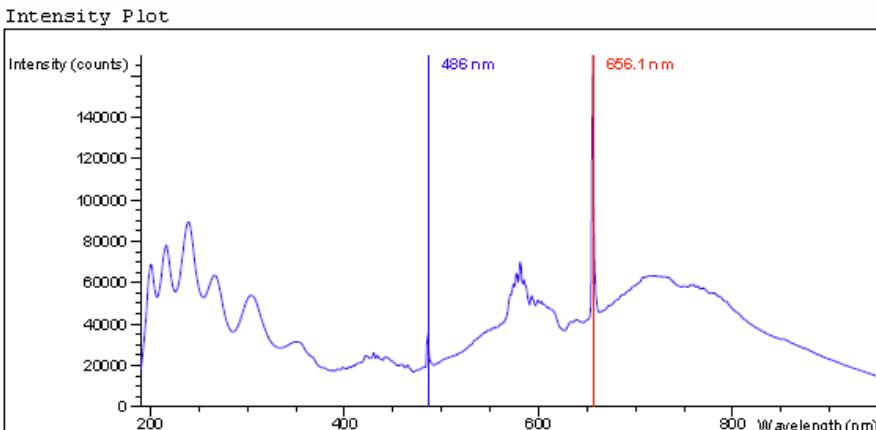
The test is for the standard flow cells (10 mm and 6 mm pathlength) only. The nano-flow cells (80 nL and 500 nL) cannot be run with this test due to its low volume.

The intensity test measures the intensity of the deuterium and tungsten lamps over the full wavelength range (190 – 950 nm). Four spectral ranges are used to evaluate the intensity spectrum. The test is used to determine the performance of the lamps and optics (see also “[Cell Test](#)” on page 184). When the test is started, the 1-nm slit is moved into the light path automatically, and the gain is set to zero. To eliminate effects due to absorbing solvents, the test should be done with water in the flow cell. The shape of the intensity spectrum is primarily dependent on the lamp, grating, and diode array characteristics. Therefore, intensity spectra will differ slightly between instruments. [Figure 54](#) on page 178 shows a typical intensity test spectrum.

Intensity Test Evaluation

The Agilent LabAdvisor, ChemStation and Instant Pilot evaluate four spectral ranges automatically, and display the limits for each range, the measured intensity counts, and *passed* or *failed* for each spectral range (see [Figure 54](#) on page 178).

Instrument: G1315C
Serial Number: DE60755000
Operator:
Date: 3/17/2009
Time: 1:24:38 PM



Intensity Test Results

	Specification	Measured	Result
Accumulated UV lamp on time		68.08 h	
Lowest intensity in range 190nm - 220nm:	> 2000 cts	18069 cts	Passed
Lowest intensity in range 221nm - 350nm:	> 5000 cts	28464 cts	Passed
Lowest intensity in range 351nm - 500nm:	> 2000 cts	16889 cts	Passed
Lowest intensity in range 501nm - 950nm:	> 2000 cts	14665 cts	Passed
Highest intensity in range 190nm - 350nm:	< 450000 cts	89478 cts	Passed
Highest intensity in range 700nm - 950nm:	< 300000 cts	63598 cts	Passed
Highest intensity for the D2 alpha line:	< 1200000 cts	169933 cts	Passed

Figure 54 Intensitiy Test Results (report)

In case of low counts in one or more ranges, start the testing with the comparison of values with flow cell vs. flow cell removed.

Contaminations of the cell windows and/or the lenses (there are 3 between vis-lamp and flow cell), will reduce the light throughput.

If the detector fails in the range 501 nm - 950 nm, check

- is the VIS-lamp ON? If not, turn it on.
- is VIS-lamp glass bulb blackended or broken? If yes, replace VIS-lamp.

9 Test Functions and Calibration

Diode Array Detector (DAD)

- does the UV-lamp show a reflective coating towards the VIS-lamp? If yes, replace UV-lamp.

Example (measured without flow cell):

VIS-LAMP OFF or defect:	110 counts
VIS-LAMP ON and OK:	13613 counts

Test Failed

Probable cause	Suggested actions
1 Absorbing solvent or air bubble in flow cell.	Ensure the flow cell is filled with water, and free from air bubbles.
2 Dirty or contaminated flow cell.	Run the cell test (see “ Cell Test ” on page 184). If the test fails, exchange the flow cell windows.
3 Dirty or contaminated optical components (achromat, windows).	Clean optical components with alcohol and lint-free cloth or replace the parts.
4 Old or non-Agilent lamp.	Exchange the lamp.

NOTE

If the lamp fails in a single range there might be no reason to change the lamp if the application is not run in that specific range.

Redo the test with removed flow cell. If the counts increase drastically (more than a factor of 2, then flow cell components are contaminated and may require maintenance/service.

If the intervals of lamp replacements are getting shorter, the Agilent service should check the optical unit for contaminated components in the light path (coupling lens, source lens, cell support assembly and flow cell windows).

Holmium Oxide Test

The holmium oxide test uses characteristic absorbance maxima of the built-in holmium oxide filter to verify wavelength accuracy (see also “[Wavelength Verification and Recalibration](#)” on page 188). When the test is started, the 1-nm slit is moved into the light path automatically. To eliminate effects due to absorbing solvents, the test should be done with water in the flow cell or with removed flow cell.

NOTE

See also “[Declaration of Conformity for HOX2 Filter](#)” on page 389.

Holmium Oxide Test Evaluation

Limits:

361.0 nm 360.0 - 362.0 nm ($\pm 1\text{nm}$)

453.7 nm 452.7 - 454.7 nm ($\pm 1\text{nm}$)

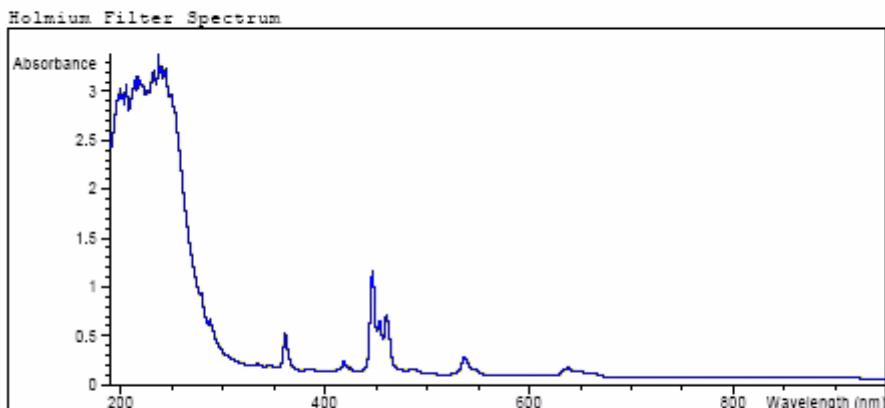
536.7 nm 535.7 - 537.7 nm ($\pm 1\text{nm}$)

The test is evaluated by the instrument, and the measured maxima are displayed automatically. The test fails if one or more of the maxima lies outside of the limits (see [Figure 55](#) on page 181).

9 Test Functions and Calibration

Diode Array Detector (DAD)

Instrument: G1315C
Serial Number: PP00000024
Operator: Wolfgang
Date: 25.02.2005
Time: 14:30:08
File: C:\CHEM32\2\DIAGNOSE\DA_D_FILTER2.DGR



Holmium Filter Test Results

	Specification	Measured	Result
Wavelength 1: 361.0 nm	360..362 nm	360.9 nm	Passed
Wavelength 2: 452.7 nm	452.7..454.7 nm	453.4 nm	Passed
Wavelength 3: 536.7 nm	535.7..537.7 nm	536.8 nm	Passed

Figure 55 Holmium Oxide Test Results (report)

Test Failed

Probable cause

- 1 Absorbing solvent or air bubble in flow cell. Ensure the flow cell is filled with water.
- 2 Incorrect calibration Recalibrate (see “[Wavelength Verification and Recalibration](#)” on page 188) and repeat the test.
- 3 Dirty or contaminated flow cell. Run the cell test (see “[Cell Test](#)” on page 184). If the test fails, exchange the flow cell windows.
- 4 Dirty or contaminated optical components (achromat, windows). Clean optical components with alcohol and lint-free cloth or replace the parts (see “[Intensity Test](#)” on page 177).
- 5 Old or non-Agilent lamp. Exchange the UV lamp.

Suggested actions

Spectral flatness test

The spectral flatness test determines the maximum noise (in mAU) on the spectrum. The test is run with the flowcell removed to eliminate effects due to absorbing solvent or a dirty flowcell. First, a detector balance is done. Next, a spectrum (no flowcell) is taken.

Theoretically, the spectrum should be flat, however, in practice, noise is superimposed on the spectrum. The amplitude of the noise is a measure of the optical and electronic performance.

NOTE

This test is part of the detector self test only, see “[Self-test](#)” on page 171.

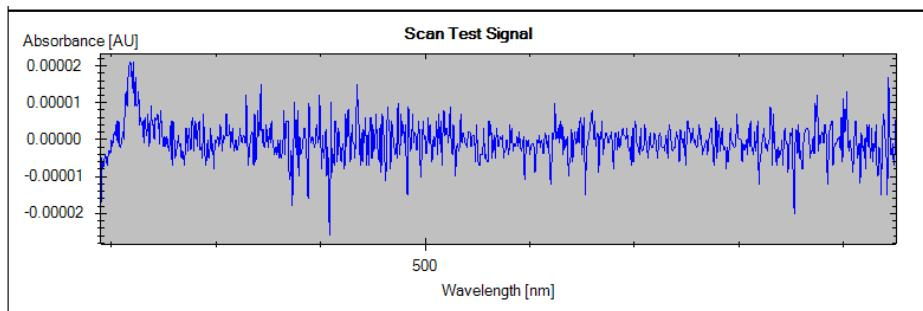


Figure 56 Spectral Flatness Test (Signals)

Spectral Flatness Evaluation

Test Evaluation

Limit is < 0.002 AU

9 Test Functions and Calibration

Diode Array Detector (DAD)

Test Failed

Probable cause	Suggested actions
1 Old or non-Agilent lamp.	<ul style="list-style-type: none">• Run the Intensity Test.• Exchange the lamp if old or defective.
2 Defective DAM board	Exchange the DAM board.
3 Defective PDA/optical unit.	Exchange the optical unit.

ASTM Noise Test

The ASTM noise test determines the detector noise over a period of 20 minutes. The test is done with the flowcell removed, so the test results are not influenced by solvent or pump effects. On completion of the test, the noise result is displayed automatically.

NOTE

This test is part of the detector self test only, see “[Self-test](#)” on page 171.

ASTM Noise Test Evaluation

Limit is ± 0.02 mAU

Test Failed

Probable cause	Suggested actions
1 Insufficient lamp warm-up time	Allow lamp to warm-up for at least 1 hour.
2 Old or non-Agilent lamp.	Exchange the lamp.

Cell Test

The cell test measures the intensity of the deuterium and tungsten lamps over the full wavelength range (190 – 950 nm), once with the flow cell installed, and once with the flow cell removed. The resulting intensity ratio is a measure of the amount of light absorbed by the flow cell. The test can be used to check for dirty or contaminated flow cell windows. When the test is started, the 1-nm slit is moved into the light path automatically, and the gain is set to zero. To eliminate effects due to absorbing solvents, the test should be done with water in the flow cell.

NOTE

This test should be performed initially with a new detector/flow cell. The values should be kept for later reference/comparison.

9 Test Functions and Calibration

Diode Array Detector (DAD)

Cell Test Evaluation

The Agilent ChemStation calculates the intensity ratio automatically. The intensity ratio (typically between 0.5 and 0.7 for new standard flow cells and 0.1 to 0.3 for new mico- and high pressure cells) is dependent on the degree of contamination of the flow cell windows, and on the type of flow cell used.

```
Instrument: G1315C
Serial Number: PP00000024
Operator: Wolfgang
Date: 25.02.2005
Time: 14:54:22
File: C:\CHEM32\2\DIAGNOSE\DAD_CELL2.DGR
```

	Result	Status
Detector Cell Test		
Date: 25.02.2005; Time: 14:54:22		
Expected total time: approx. 45 s.		
Test Procedure:		
1. If cell not in place, install it		done
2. Measuring intensity with cell	42053	done
3. Remove cell		done
4. Measuring intensity without cell	42034	done
5. Calculating intensity ratio	1.0	done

Figure 57 Cell Test Results (report) with no flow cell inserted

NOTE

This test can be used for the standard flow cells only. The nano flow cells will give very low values due to their design.

Test Failed (low ratio value)

Probable cause	Suggested actions
1 Absorbing solvent or air bubble in flow cell.	Ensure the flow cell is filled with water, and free from air bubbles.
2 Dirty or contaminated flow cell.	Exchange the flow cell windows.

Using the Built-in Test Chromatogram

This function is available from the Agilent ChemStation, Lab Advisor and Instant Pilot.

The built-in Test Chromatogram can be used to check the signal path from the detector to the data system and the data analysis or via the analog output to the integrator or data system. The chromatogram is continuously repeated until a stop is executed either by means of a stop time or manually.

NOTE

The peak height is always the same but the area and the retention time depend on the set peakwidth, see example below.

Procedure Using the Agilent LabAdvisor

This procedure works for all Agilent 1200 Infinity detectors (DAD, MWD, VWD, FLD and RID). The example figure is from the RID detector.

- 1 Assure that the default LC method is loaded via the control software.
- 2 Start the Agilent Lab Advisor software (B.01.03 SP4 or later) and open the detector's **Tools** selection.
- 3 Open the test chromatogram screen



- 4 Turn the **Test Chromatogram** on.
- 5 Change to the detector's **Module Service Center** and add the detector signal to the Signal Plot window.

9 Test Functions and Calibration

Diode Array Detector (DAD)

6 To start a test chromatogram enter in the command line: STRT

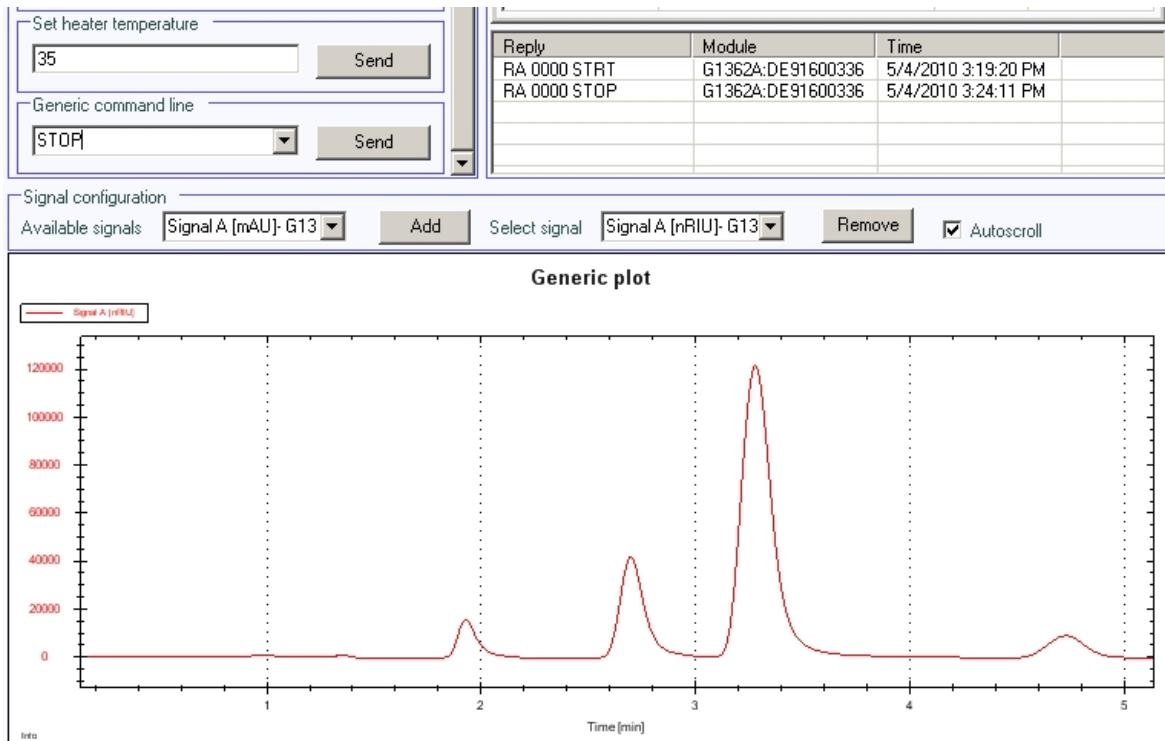


Figure 58 Test Chromatogram with Agilent Lab Advisor

7 To stop the test chromatogram enter in the command line: STOP

NOTE

The test chromatogram is switched off automatically at the end of a run.

Wavelength Verification and Recalibration

The detector uses the alpha (656.1 nm) and beta (486 nm) emission lines of the deuterium lamp for wavelength calibration. The sharp emission lines enable more accurate calibration than is possible with holmium oxide. When verification is started, the 1-nm slit is moved into the light path automatically, and the gain is set to zero. To eliminate effects due to absorbing solvents, the test should be done with bubble free degassed HPLC water in the flow cell.

If a deviation is found and displayed, it can be recalibrated by pressing Adjust. The deviations are tracked in the Calibration History (diagnosis buffer in the detector).

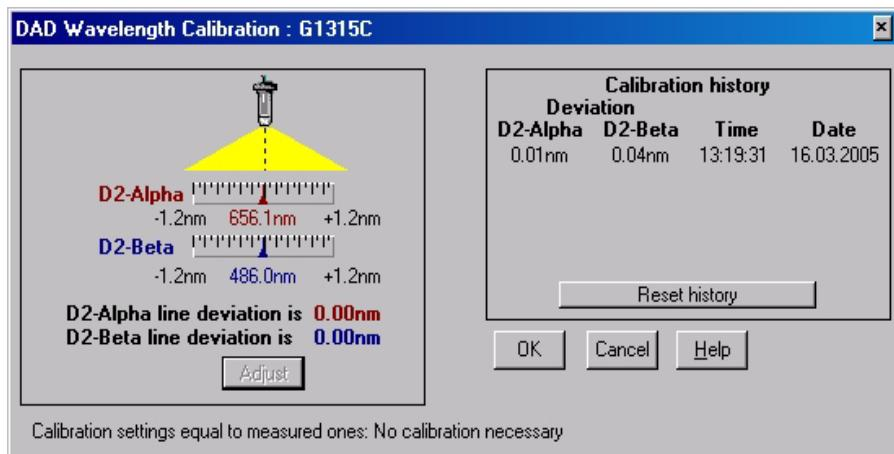


Figure 59 Wavelength Verification and Recalibration

Wavelength calibration should be done

- after maintenance of the flow cell,
- lamp exchange, or
- after a major repair, like processor board or optical unit exchange, see also “Replacing the Module’s Firmware” on page 325.

After calibration, the holmium oxide test (see [Figure 55](#) on page 181) provides verification of wavelength accuracy at three additional wavelengths.

Diagnosis Information on Agilent ChemStation

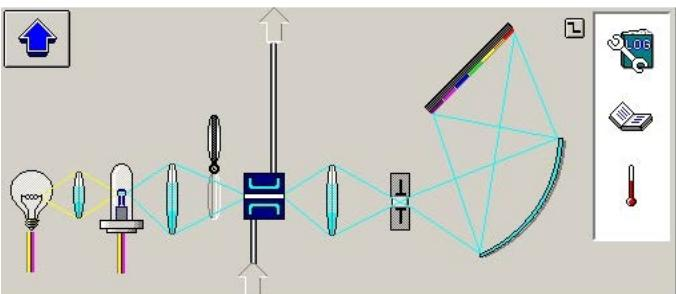
The Agilent ChemStation provides certain information on various assemblies in the detector.

- 1 Open the Diagnosis and change to detailed view.
- 2 Click on the assembly of interest and select “Update Variables Display”.
[Figure](#) on page 190 shows an example.

[Table 43](#) on page 189 lists the assemblies which provide detailed information.

Table 43 Diagnosis - detailed information

Assembly	Details
General	Product number, firmware revision, manufacturing date (of instrument or, if replaced, of main board), serial number, accumulated on-time, spectrometer serial number.
Vis lamp	Accumulated on-time, actual on-time, lamp switch on, accumulated on time switches, switch on voltage/current, lamp current, lamp voltage/current
UV lamp	Accumulated on-time, actual on-time, lamp ignitions, accumulated on time reset, ignition voltage/current, lamp voltage, lamp voltage/current.
	Available with ID-tag lamp only. product number, serial number, production date, last intensity test
Holmium Filter	Filter movements, filter moves reset log
Flow Cell	Available with ID-tag flow cells only. Product number, production date, pathlength, max pressure, serial number, volume, last cell test
Micro Slit	Slit movements, filter moves reset log



Diode Array Detector - UV Lamp

Accum. UV On Time	179.83 h	Accum. On Time Reset	Show
UV Lamp On Time	0.00 h	Ignition Voltage/Current	Show
Number of Ignitions	70	UV Lamp Voltage	Show
Product Number	2140-0820	UV Lamp Volt./Current	Show
Serial Number	213913	Last Intensity Test	02.01.2006
Production Date	08.10.2004		

Diode Array Detector - Cell

Product Number	G1315-60022	Serial Number	DE515H0876
Production Date	22.09.2005	Path Length	10 mm
Volume	13000 nl	Max Pressure	120 bar
Last Cell Test	22.09.2005		

Details of UV-lamp with RFID tag

Details of flow cell with RFID tag

Buttons: Copy to Memo Pad, Send Changes, Cancel Changes.

Figure 60 Diagnosis screen in Agilent ChemStation (detailed view)

D/A Converter (DAC) Test

The detector provides analog output of chromatographic signals for use with integrators, chart recorders or data systems. The analog signal is converted from the digital format by the digital-analog-converter (DAC).

The DAC test is used to verify correct operation of the digital-analog-converter by applying a digital test signal to the DAC.

The DAC outputs an analog signal of approximately 50 mV (if the zero offset of the analog output is set to the default value of 5 %) which can be plotted on an integrator. A continuous square wave with an amplitude of 10 µV and a frequency of approximately 1 cycle/24 seconds is applied to the signal.

The amplitude of the square wave and the peak-to-peak noise are used to evaluate the DAC test.

When If the analog detector signal is noisy or missing.

Preparations Lamp must be on for at least 10 minutes. Connect integrator, chart recorder or data system to the detector analog output.

Running the test with Agilent LabAdvisor

- 1 Run the **D/A Converter (DAC) Test** (for further information see Online-Help of user interface).

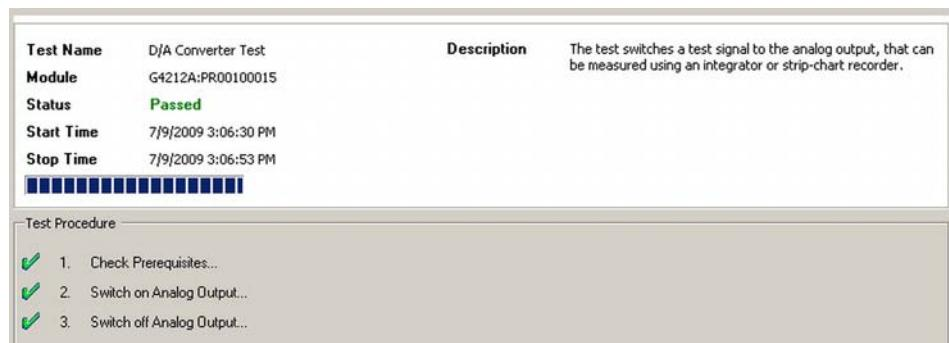


Figure 61 D/A Converter (DAC) Test – Results

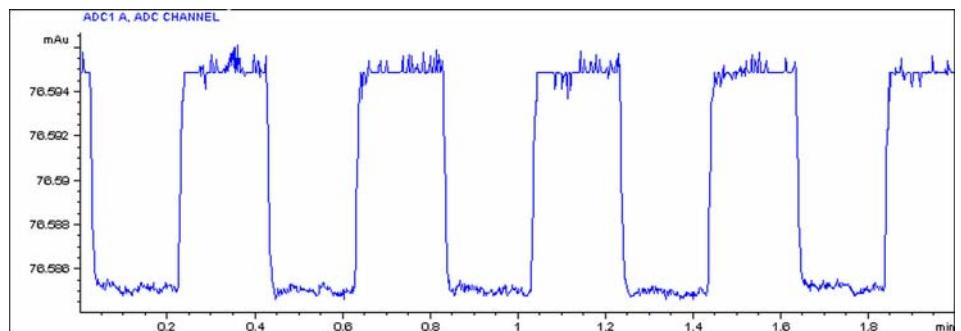


Figure 62 D/A Converter (DAC) Test – Example of Integrator Plot

Running the Test with Instant Pilot

The test can be started via the command line.

- 1 To start the test

TEST: DAC 1

Reply: RA 00000 TEST:DAC 1

- 2 To stop the test

TEST:DAC 0

Reply: RA 00000 TEST:DAC 0

Test Evaluation

The noise on the step should be less than 3 µV.

Probable cause	Suggested actions
----------------	-------------------

1 Bad cable or grounding problem between detector and external device.

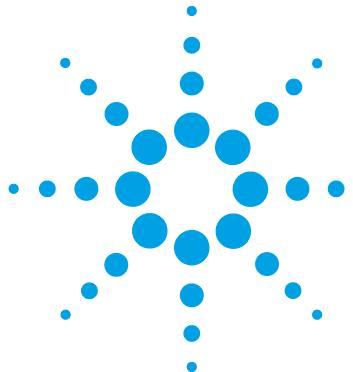
Check or replace the cable.

2 Defective detector main board.

Please contact your Agilent service representative.

9 Test Functions and Calibration

Diode Array Detector (DAD)



10 Error Information

What are Error Messages?	196
General Error Messages	197
Compensation Sensor Open	197
Compensation Sensor Short	197
Fan Failed	198
Leak	199
Leak Sensor Open	199
Leak Sensor Short	200
Remote Timeout	200
Shutdown	201
Lost CAN Partner	201
Timeout	202
Pump Error Messages	203
Encoder Missing	203
Index Adjustment	203
Index Limit	204
Index Missing	204
Initialization Failed	205
Missing Pressure Reading	205
Motor-Drive Power	206
Pressure Above Upper Limit	207
Pressure Below Lower Limit	207
Pressure Signal Missing	208
Pump Configuration	209
Pump Head Missing	209
Restart Without Cover	209
Servo Restart Failed	210
Stroke Length	211



10 Error Information

Diode Array Detector (DAD)

Temperature Limit Exceeded	211
Temperature Out of Range	212
Valve Failed (MCGV)	212
MCGV Fuse	213
Wait Timeout	213
Solvent Zero Counter	214
Autosampler Error Messages	216
Arm Movement Failed	216
Initialization Failed	217
Initialization with Vial	217
Invalid Vial Position	218
Metering Home Failed	219
Missing Vial	219
Missing Wash Vial	220
Motor Temperature	220
Needle Down Failed	221
Needle Up Failed	222
Safety Flap Missing	223
Valve to Bypass Failed	223
Valve to Mainpass Failed	224
Vial in Gripper	224
General Detector Error Messages	226
Heater at fan assembly failed	226
Heater Power At Limit	226
Illegal Temperature Value from Sensor on Main Board	227
Illegal Temperature Value from Sensor at Air Inlet	227
UV Lamp Current	227
UV Lamp Voltage	228
VWD Detector Error Messages	229
ADC Hardware Error	229
Wavelength calibration setting failed	229
Cutoff filter doesn't decrease the light intensity at 250 nm	230
Filter Missing	230
Grating or Filter Motor Errors	231
Grating Missing	232

No heater current	232
Wavelength holmium check failed	233
Ignition Failed	233
Wavelength test failed	234
DAD Detector Error Messages	235
Visible Lamp Current	235
Visible Lamp Voltage	235
Diode Current Leakage	236
UV Ignition Failed	236
UV Heater Current	237
Calibration Values Invalid	238
Holmium Oxide Test Failed	238
Wavelength Recalibration Lost	239
DSP Not Running	239
No Run Data Available In Device	239
Instrument Logbook	240

This chapter provides information on the error messages that might be displayed, and gives the possible causes and suggestions on their solutions.

10 Error Information

What are Error Messages?

What are Error Messages?

Error messages are displayed in the user interface when an electronic, mechanical, or hydraulic (flow path) failure occurs which requires attention before the analysis can be continued (for example, repair, or exchange of consumables is necessary). In the event of such a failure, the red status indicator at the front of the module is switched on, and an entry is written into the module logbook.

If an error occurs outside a method run, other modules will not be informed about this error. If it occurs within a method run, all connected modules will get a notification, all LEDs get red and the run will be stopped. Depending on the module type, this stop is implemented differently. For example, for a pump the flow will be stopped for safety reasons. For a detector, the lamp will stay on in order to avoid equilibration time. Depending on the error type, the next run can only be started, if the error has been resolved, for example liquid from a leak has been dried. Errors for presumably single time events can be recovered by switching on the system in the user interface.

Special handling is done in case of a leak. As a leak is a potential safety issue and may have occurred at a different module from where it has been observed, a leak always causes a shutdown of all modules, even outside a method run.

In all cases, error propagation is done via the CAN bus or via an APG/ERI remote cable (see documentation for the APG/ERI interface).

General Error Messages

General error messages are generic to all Agilent series HPLC modules and may show up on other modules as well.

Compensation Sensor Open

Error ID: 0081

The ambient-compensation sensor (NTC) on the power switch board in the module has failed (open circuit).

The resistance across the temperature compensation sensor (NTC) on the power switch board is dependent on ambient temperature. The change in resistance is used by the leak circuit to compensate for ambient temperature changes. If the resistance across the sensor increases above the upper limit, the error message is generated.

Probable cause	Suggested actions
1 Loose connection between the power switch board and the main board	Please contact your Agilent service representative.
2 Defective power switch assembly	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Compensation Sensor Short

Error ID: 0080

The ambient-compensation sensor (NTC) on the power switch board in the module has failed (open circuit).

10 Error Information

General Error Messages

The resistance across the temperature compensation sensor (NTC) on the power switch board is dependent on ambient temperature. The change in resistance is used by the leak circuit to compensate for ambient temperature changes. If the resistance across the sensor falls below the lower limit, the error message is generated.

Probable cause	Suggested actions
1 Defective power switch assembly	Please contact your Agilent service representative.
2 Loose connection between the power switch board and the main board	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Fan Failed

Error ID: 0068

The cooling fan in the module has failed.

The hall sensor on the fan shaft is used by the main board to monitor the fan speed. If the fan speed falls below a certain limit for a certain length of time, the error message is generated.

Depending on the module, assemblies (e.g. the lamp in the detector) are turned off to assure that the module does not overheat inside.

Probable cause	Suggested actions
1 Fan cable disconnected.	Please contact your Agilent service representative.
2 Defective fan.	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Leak

Error ID: 0064

A leak was detected in the module.

The signals from the two temperature sensors (leak sensor and board-mounted temperature-compensation sensor) are used by the leak algorithm to determine whether a leak is present. When a leak occurs, the leak sensor is cooled by the solvent. This changes the resistance of the leak sensor which is sensed by the leak-sensor circuit on the main board.

Probable cause	Suggested actions
1 Loose fittings.	Ensure all fittings are tight.
2 Broken capillary.	Exchange defective capillaries.

Leak Sensor Open

Error ID: 0083

The leak sensor in the module has failed (open circuit).

The current through the leak sensor is dependent on temperature. A leak is detected when solvent cools the leak sensor, causing the leak-sensor current to change within defined limits. If the current falls outside the lower limit, the error message is generated.

Probable cause	Suggested actions
1 Leak sensor not connected to the main board.	Please contact your Agilent service representative.
2 Defective leak sensor.	Please contact your Agilent service representative.
3 Leak sensor incorrectly routed, being pinched by a metal component.	Please contact your Agilent service representative.

10 Error Information

General Error Messages

Leak Sensor Short

Error ID: 0082

The leak sensor in the module has failed (short circuit).

The current through the leak sensor is dependent on temperature. A leak is detected when solvent cools the leak sensor, causing the leak sensor current to change within defined limits. If the current increases above the upper limit, the error message is generated.

Probable cause	Suggested actions
1 Defective leak sensor.	Please contact your Agilent service representative.
2 Leak sensor incorrectly routed, being pinched by a metal component.	Please contact your Agilent service representative.

Remote Timeout

Error ID: 0070

A not-ready condition is still present on the remote input. When an analysis is started, the system expects all not-ready conditions (for example, a not-ready condition during detector balance) to switch to run conditions within one minute of starting the analysis. If a not-ready condition is still present on the remote line after one minute the error message is generated.

Probable cause	Suggested actions
1 Not-ready condition in one of the instruments connected to the remote line.	Ensure the instrument showing the not-ready condition is installed correctly, and is set up correctly for analysis.
2 Defective remote cable.	Exchange the remote cable.
3 Defective components in the instrument showing the not-ready condition.	Check the instrument for defects (refer to the instrument's documentation).

Shutdown

Error ID: 0063

An external instrument has generated a shutdown signal on the remote line.

The module continually monitors the remote input connectors for status signals. A LOW signal input on pin 4 of the remote connector generates the error message.

Probable cause	Suggested actions
1 Leak detected in another module with a CAN connection to the system.	Fix the leak in the external instrument before restarting the module.
2 Leak detected in an external instrument with a remote connection to the system.	Fix the leak in the external instrument before restarting the module.
3 Shut-down in an external instrument with a remote connection to the system.	Check external instruments for a shut-down condition.
4 The degasser failed to generate sufficient vacuum for solvent degassing.	Check the vacuum degasser for an error condition. Refer to the <i>Service Manual</i> for the degasser or the pump that has the degasser built-in.

Lost CAN Partner

Error ID: 0071

During an analysis, the internal synchronization or communication between one or more of the modules in the system has failed.

The system processors continually monitor the system configuration. If one or more of the modules is no longer recognized as being connected to the system, the error message is generated.

10 Error Information

General Error Messages

Probable cause	Suggested actions
1 CAN cable disconnected.	<ul style="list-style-type: none">Ensure all the CAN cables are connected correctly.Ensure all CAN cables are installed correctly.
2 Defective CAN cable.	Exchange the CAN cable.
3 Defective main board in another module.	Switch off the system. Restart the system, and determine which module or modules are not recognized by the system.

Timeout

Error ID: 0062

The timeout threshold was exceeded.

Probable cause	Suggested actions
1 The analysis was completed successfully, and the timeout function switched off the module as requested.	Check the logbook for the occurrence and source of a not-ready condition. Restart the analysis where required.
2 A not-ready condition was present during a sequence or multiple-injection run for a period longer than the timeout threshold.	Check the logbook for the occurrence and source of a not-ready condition. Restart the analysis where required.

Pump Error Messages

These errors are specific to the pump.

Encoder Missing

Error ID: 2046, 2050, 2510

The optical encoder on the pump motor in the module is missing or defective.

The processor checks the presence of the pump encoder connector every 2 s. If the connector is not detected by the processor, the error message is generated.

Probable cause	Suggested actions
1 Defective or disconnected pump encoder connector.	Please contact your Agilent service representative.
2 Defective pump drive assembly.	Please contact your Agilent service representative.

Index Adjustment

Error ID: 2204, 2214

The encoder index position in the module is out of adjustment.

During initialization, the first piston is moved to the mechanical stop. After reaching the mechanical stop, the piston reverses direction until the encoder index position is reached. If the time to reach the index position is too long, the error message is generated.

10 Error Information

Pump Error Messages

Probable cause	Suggested actions
1 Irregular or sticking drive movement.	Remove the pump head, and examine the seals, pistons, and internal components for signs of wear, contamination or damage. Exchange components as required.
2 Defective pump drive assembly.	Please contact your Agilent service representative.

Index Limit

Error ID: 2203, 2213

The time required by the piston to reach the encoder index position was too short (pump).

During initialization, the first piston is moved to the mechanical stop. After reaching the mechanical stop, the piston reverses direction until the encoder index position is reached. If the index position is reached too fast, the error message is generated.

Probable cause	Suggested actions
1 Irregular or sticking drive movement.	Remove the pump head, and examine the seals, pistons, and internal components for signs of wear, contamination or damage. Exchange components as required.
2 Defective pump drive assembly.	Please contact your Agilent service representative.

Index Missing

Error ID: 2205, 2215, 2505

The encoder index position in the module was not found during initialization.

During initialization, the first piston is moved to the mechanical stop. After reaching the mechanical stop, the piston reverses direction until the encoder

index position is reached. If the index position is not recognized within a defined time, the error message is generated.

Probable cause	Suggested actions
1 Disconnected or defective encoder cable.	Please contact your Agilent service representative.
2 Defective pump drive assembly.	Please contact your Agilent service representative.

Initialization Failed

Error ID: 2207, 2217

The module failed to initialize successfully within the maximum time window. A maximum time is assigned for the complete pump-initialization cycle. If the time is exceeded before initialization is complete, the error message is generated.

Probable cause	Suggested actions
1 Blocked (passive or active) inlet valve.	Exchange the (passive or active) inlet valve.
2 Defective pump drive assembly.	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Missing Pressure Reading

Error ID: 2054

The pressure readings read by the pump ADC (analog-digital converter) are missing.

The ADC reads the pressure signal from the damper every 1ms. If the readings are missing for longer than 10 s, the error message is generated.

10 Error Information

Pump Error Messages

Probable cause	Suggested actions
1 Damper disconnected.	Please contact your Agilent service representative.
2 Defective damper.	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Motor-Drive Power

Error ID: 2041, 2042

The current drawn by the pump motor exceeded the maximum limit.

Blockages in the flow path are usually detected by the pressure sensor in the damper, which result in the pump switching off when the upper pressure limit is exceeded. If a blockage occurs before the damper, the pressure increase cannot be detected by the pressure sensor and the module will continue to pump. As pressure increases, the pump drive draws more current. When the current reaches the maximum limit, the module is switched off, and the error message is generated.

Probable cause	Suggested actions
1 Flow path blockage in front of the damper.	Ensure the capillaries and frits between the pump head and damper inlet are free from blockage.
2 Blocked (passive or active) inlet valve.	Exchange the (passive or active) inlet valve.
3 Blocked outlet valve.	Exchange the outlet valve.
4 High friction (partial mechanical blockage) in the pump drive assembly.	Remove the pump-head assembly. Ensure there is no mechanical blockage of the pump-head assembly or pump drive assembly.

Probable cause	Suggested actions
5 Defective pump drive assembly.	Please contact your Agilent service representative.
6 Defective main board.	Please contact your Agilent service representative.

Pressure Above Upper Limit

Error ID: 2014, 2500

The system pressure has exceeded the upper pressure limit.

Probable cause	Suggested actions
1 Upper pressure limit set too low.	Ensure the upper pressure limit is set to a value suitable for the analysis.
2 Blockage in the flowpath (after the damper).	Check for blockage in the flow path. The following components are particularly subject to blockage: inline filter frit, needle (autosampler), seat capillary (autosampler), sample loop (autosampler), column frits and capillaries with small internal diameters (e.g. 50 µm ID).
3 Defective damper.	Please contact your Agilent service representative.
4 Defective main board.	Please contact your Agilent service representative.

Pressure Below Lower Limit

Error ID: 2015, 2501

The system pressure has fallen below the lower pressure limit.

10 Error Information

Pump Error Messages

Probable cause	Suggested actions
1 Lower pressure limit set too high.	Ensure the lower pressure limit is set to a value suitable for the analysis.
2 Air bubbles in the mobile phase.	<ul style="list-style-type: none">Make sure that the degasser is in flow path and works correctly. Purge the module.Ensure solvent inlet filters are not blocked.
3 Leak.	<ul style="list-style-type: none">Inspect the pump head, capillaries and fittings for signs of a leak.Purge the module. Run a pressure test to determine whether the seals or other module components are defective.
4 Defective damper.	Please contact your Agilent service representative.
5 Defective main board.	Please contact your Agilent service representative.

Pressure Signal Missing

Error ID: 2016

The pressure signal of the damper is missing.

The pressure signal of the damper must be within a specific voltage range. If the pressure signal is missing, the processor detects a voltage of approximately -120 mV across the damper connector.

Probable cause	Suggested actions
1 Damper disconnected.	Please contact your Agilent service representative.
2 Defective damper.	Please contact your Agilent service representative.

Pump Configuration

Error ID: 2060

At switch-on, the gradient pump has recognized a new pump configuration.

The gradient pump is assigned its configuration at the factory. If the gradient valve is disconnected, and the gradient pump is rebooted, the error message is generated. However, the pump will function as an isocratic pump in this configuration. The error message reappears after each switch-on.

Probable cause	Suggested actions
1 Gradient valve disconnected.	Reconnect the gradient valve.

Pump Head Missing

Error ID: 2202, 2212

The pump-head end stop in the pump was not found.

When the pump restarts, the metering drive moves forward to the mechanical end stop. Normally, the end stop is reached within 20 s, indicated by an increase in motor current. If the end point is not found within 20 s, the error message is generated.

Probable cause	Suggested actions
1 Pump head not installed correctly (screws not secured, or pump head not seated correctly).	Install the pump head correctly. Ensure nothing (e.g. capillary) is trapped between the pump head and body.
2 Broken piston.	Exchange the piston.

Restart Without Cover

Error ID: 2502

The module was restarted with the top cover and foam open.

10 Error Information

Pump Error Messages

The sensor on the main board detects when the top foam is in place. If the module is restarted with the foam removed, the module switches off within 30 s, and the error message is generated.

Probable cause	Suggested actions
1 The module started with the top cover and foam removed.	Please contact your Agilent service representative.

Servo Restart Failed

Error ID: 2201, 2211

The pump motor in the module was unable to move into the correct position for restarting.

When the module is switched on, the first step is to switch on the C phase of the variable reluctance motor. The rotor should move to one of the C positions. The C position is required for the servo to be able to take control of the phase sequencing with the commutator. If the rotor is unable to move, or if the C position cannot be reached, the error message is generated.

Probable cause	Suggested actions
1 Mechanical blockage of the module.	Remove the pump-head assembly. Ensure there is no mechanical blockage of the pump-head assembly or pump drive assembly.
2 Disconnected or defective cable.	Please contact your Agilent service representative.
3 Blocked (passive or active) inlet valve.	Exchange the (passive or active) inlet valve.
4 Defective pump drive assembly.	Please contact your Agilent service representative.
5 Defective main board.	Please contact your Agilent service representative.

Stroke Length

Error ID: 2206, 2216

The distance between the lower piston position and the upper mechanical stop is out of limits (pump).

During initialization, the module monitors the drive current. If the piston reaches the upper mechanical stop position before expected, the motor current increases as the module attempts to drive the piston beyond the mechanical stop. This current increase causes the error message to be generated.

Probable cause	Suggested actions
1 Defective pump drive assembly.	Please contact your Agilent service representative.

Temperature Limit Exceeded

Error ID: 2517

The temperature of one of the motor-drive circuits is too high.

The processor continually monitors the temperature of the drive circuits on the main board. If excessive current is being drawn for long periods, the temperature of the circuits increases. If the temperature exceeds the upper limit, the error message is generated.

Probable cause	Suggested actions
1 High friction (partial mechanical blockage) in the pump drive assembly.	Remove the pump-head assembly. Ensure there is no mechanical blockage of the pump-head assembly or pump drive assembly.
2 Partial blockage of the flowpath in front of the damper.	Ensure the outlet valve is not blocked.

10 Error Information

Pump Error Messages

Probable cause	Suggested actions
3 Defective pump drive assembly.	Please contact your Agilent service representative.
4 Defective main board.	Please contact your Agilent service representative.

Temperature Out of Range

Error ID: 2517

The temperature sensor readings in the motor-drive circuit are out of range.

The values supplied to the ADC by the hybrid sensors must be between 0.5 V and 4.3 V. If the values are outside this range, the error message is generated.

Probable cause	Suggested actions
1 Defective main board.	Please contact your Agilent service representative.

Valve Failed (MCGV)

Error ID: 2040

Valve 0 Failed: valve A

Valve 1 Failed: valve B

Valve 2 Failed: valve C

Valve 3 Failed: valve D

One of the valves of the multi-channel gradient valve has failed to switch correctly.

The processor monitors the valve voltage before and after each switching cycle. If the voltages are outside expected limits, the error message is generated.

Probable cause	Suggested actions
1 Gradient valve disconnected.	Ensure the gradient valve is connected correctly.
2 Connection cable (inside instrument) not connected.	Please contact your Agilent service representative.
3 Connection cable (inside instrument) defective.	Please contact your Agilent service representative.
4 Gradient valve defective.	Exchange the gradient valve.

MCGV Fuse

Error ID: 2043

Valve Fuse 0: Channels A and B

Valve Fuse 1: Channels C and D

The gradient valve in the quaternary pump has drawn excessive current causing the electronic fuse to open.

Probable cause	Suggested actions
1 Defective gradient valve.	Restart the quaternary pump. If the error message appears again, exchange the gradient valve.
2 Defective connection cable (front panel to main board).	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Wait Timeout

Error ID: 2053

10 Error Information

Pump Error Messages

When running certain tests in the diagnostics mode or other special applications, the pump must wait for the pistons to reach a specific position, or must wait for a certain pressure or flow to be reached. Each action or state must be completed within the timeout period, otherwise the error message is generated.

Possible Reasons for a Wait Timeout:

- Pressure not reached.
- Pump channel A did not reach the delivery phase.
- Pump channel B did not reach the delivery phase.
- Pump channel A did not reach the take-in phase.
- Pump channel B did not reach the take-in phase.
- Solvent volume not delivered within the specified time.

Probable cause	Suggested actions
1 Purge valve open.	Ensure that purge valve is closed.
2 Leak at fittings, purge valve, active inlet valve, outlet valve or piston seals.	<ul style="list-style-type: none">· Ensure pump components are seated correctly. If there are still signs of a leak, exchange the appropriate seal (purge valve, active inlet valve, outlet valve, piston seal).· Exchange defective capillaries.
3 Flow changed after starting test.	Ensure correct operating condition for the special application in use.
4 Defective pump drive assembly.	Please contact your Agilent service representative.

Solvent Zero Counter

Error ID: 2055, 2524

Pump firmware version A.02.32 and higher allow to set solvent bottle fillings in the data system. If the volume level in the bottle falls below the specified value the error message appears when the feature is configured accordingly.

Probable cause	Suggested actions
1 Volume in bottle below specified volume.	Refill bottles and reset solvent counters.
2 Incorrect setting.	Make sure the limits are set correctly.

10 Error Information

Autosampler Error Messages

Autosampler Error Messages

These errors are specific to the autosampler.

Arm Movement Failed

Error ID: 4002

The transport assembly was unable to complete a movement in one of the axes.

The processor defines a certain time window for the successful completion of a movement in any particular axis. The movement and position of the transport assembly is monitored by the encoders on the stepper motors. If the processor does not receive the correct position information from the encoders within the time window, the error message is generated.

See figure [Figure 35](#) on page 109 for axes identification.

- **Arm Movement 0 Failed:** X-axis.
- **Arm Movement 1 Failed:** Z-axis.
- **Arm Movement 2 Failed:** Theta (gripper rotation).
- **Arm Movement 3 Failed:** Gripper (gripper fingers open/close).

Probable cause	Suggested actions
1 Mechanical obstruction.	Ensure unobstructed movement of the transport assembly.
2 High friction in the transport assembly.	Please contact your Agilent service representative.
3 Defective motor assembly.	Please contact your Agilent service representative.

Probable cause	Suggested actions
4 Defective sample transport assembly flex board.	Please contact your Agilent service representative.
5 Defective main board.	Please contact your Agilent service representative.

Initialization Failed

Error ID: 4020

The autosampler failed to complete initialization correctly.

The autosampler initialization procedure moves the needle arm and transport assembly to their home positions in a predefined sequence. During initialization, the processor monitors the position sensors and motor encoders to check for correct movement. If one or more of the movements is not successful, or is not detected, the error message is generated.

Probable cause	Suggested actions
1 Mechanical obstruction.	Ensure unobstructed movement of the transport assembly.
2 Defective sampling unit flex board.	Please contact your Agilent service representative.
3 Defective transport assembly flex board.	Please contact your Agilent service representative.
4 Defective sampling unit motor.	Please contact your Agilent service representative.
5 Defective main board.	Please contact your Agilent service representative.

Initialization with Vial

Error ID: 4028

10 Error Information

Autosampler Error Messages

The autosampler attempted to initialize with a vial still in the gripper.

During initialization, the autosampler checks correct operation of the gripper by closing and opening the gripper fingers while monitoring the motor encoder. If a vial is still in the gripper when initialization is started, the gripper fingers cannot close causing the error message to be generated.

Probable cause	Suggested actions
1 Vial still in gripper.	Remove the vial using the Release Vial function in the user interface. Reinitialize the autosampler.

Invalid Vial Position

Error ID: 4042

The vial position defined in the method or sequence does not exist.

The reflection sensors on the transport assembly flex board are used to automatically check which sample trays are installed (coding on tray). If the vial position does not exist in the current sample tray configuration, the error message is generated.

Probable cause	Suggested actions
1 Incorrect tray or trays installed.	Install the correct trays, or edit the method or sequence accordingly.
2 Incorrect vial positions defined in the method or sequence.	Check method settings vs. installed tray.
3 Tray recognition defective (dirty sample tray or defective transport assembly flex board).	Ensure the coding surfaces of the sample tray are clean (located at the rear of the sample tray). If this does not solve the problem, contact your Agilent service representative.

Metering Home Failed

Error ID: 4054, 4704

The metering piston has failed to move back to the home position.

The home position sensor on the sampling unit flex board monitors the home position of the piston. If the piston fails to move to the home position, or if the sensor fails to recognize the piston position, the error message is generated.

Probable cause	Suggested actions
1 Defective sensor or main board.	Please contact your Agilent service representative.
2 Broken piston.	Exchange the metering piston and seal.
3 Defective metering-drive motor.	Please contact your Agilent service representative.
4 Defective main board.	Please contact your Agilent service representative.

Missing Vial

Error ID: 4019, 4034, 4541, 4706

No vial was found in the position defined in the method or sequence.

When the gripper arm picks a vial out of the sample tray, the processor monitors the gripper motor encoder. If a vial is present, the closing of the gripper fingers is limited by the vial. However, if no vial is present, the gripper fingers close too far. This is sensed by the processor (encoder position), causing the error message to be generated.

Probable cause	Suggested actions
1 No vial in the position defined in the method or sequence.	Install the sample vial in the correct position, or edit the method or sequence accordingly.
2 Incorrect gripper alignment.	Align gripper.

10 Error Information

Autosampler Error Messages

Probable cause	Suggested actions
3 Defective gripper assembly (defective gripper fingers or belt).	Exchange the gripper assembly.
4 Defective transport assembly flex board.	Please contact your Agilent service representative.

Missing Wash Vial

Error ID: 4035, 4542, 4707

The wash vial programmed in the method was not found.

When the gripper arm picks a vial out of the sample tray, the processor monitors the gripper motor encoder. If a vial is present, the closing of the gripper fingers is limited by the vial. However, if no vial is present, the gripper fingers close too far. This is sensed by the processor (encoder position), causing the error message to be generated.

Probable cause	Suggested actions
1 No wash vial in the position defined in the method.	Install the wash vial in the correct position, or edit the method accordingly.

Motor Temperature

Error ID: 4027, 4040, 4261, 4451

One of the motors of the transport assembly has drawn excessive current, causing the motor to become too hot. The processor has switched OFF the motor to prevent damage to the motor.

See figure [Figure 35](#) on page 109 for motor identification.

- **Motor 0 temperature:** X-axis motor.
- **Motor 1 temperature:** Z-axis motor.
- **Motor 2 temperature:** Theta (gripper rotation) motor.
- **Motor 3 temperature:** Gripper motor (motor for gripper fingers).

The processor monitors the current drawn by each motor and the time the motor is drawing current. The current drawn by the motors is dependent on the load on each motor (friction, mass of components etc.). If the current drawn is too high, or the time the motor draws current is too long, the error message is generated.

Probable cause	Suggested actions
1 Mechanical obstruction.	Ensure unobstructed movement of the transport assembly.
2 High friction in the transport assembly.	Please contact your Agilent service representative.
3 Motor belt tension too high.	Switch OFF the autosampler at the power switch. Wait at least 10 min before switching on again.
4 Defective motor.	Please contact your Agilent service representative.
5 Defective transport assembly flex board.	Please contact your Agilent service representative.

Needle Down Failed

Error ID: 4018

The needle arm failed to move down into the needle seat.

The lower position of the needle arm is monitored by a position sensor on the sampling unit flex board. The sensor detects the successful completion of the needle movement to the needle seat position. If the needle fails to reach the end point, or if the sensor fails to recognize the needle arm movement, the error message is generated.

10 Error Information

Autosampler Error Messages

Probable cause	Suggested actions
1 Needle installed incorrectly, or wrong needle type (too long).	Ensure the correct needle type is used, and installed correctly.
2 Defective or dirty position sensor.	Please contact your Agilent service representative.
3 Defective motor.	Please contact your Agilent service representative.
4 Sticking spindle assembly.	Please contact your Agilent service representative.
5 Defective main board.	Please contact your Agilent service representative.

Needle Up Failed

Error ID: 4017

The needle arm failed to move successfully from the seat or out of the vial to the upper position.

The upper position of the needle arm is monitored by a position sensor on the sampling unit flex board. The sensor detects the successful completion of the needle movement to the upper position. If the needle fails to reach the end point, or if the sensor fails to recognize the needle arm movement, the error message is generated.

Probable cause	Suggested actions
1 Defective or dirty position sensor.	Please contact your Agilent service representative.
2 Defective motor.	Please contact your Agilent service representative.

Probable cause	Suggested actions
3 Sticking spindle assembly.	Please contact your Agilent service representative.
4 Defective main board.	Please contact your Agilent service representative.

Safety Flap Missing

Error ID: 4032

The safety flap was not detected.

Before the needle moves down into the needle seat to inject sample, the safety flap locks into position. Next, the gripper checks the safety flap by trying to move the safety flap away from the needle. If the gripper is able to move beyond the safety flap position (safety flap not in position), the error message is generated.

Probable cause	Suggested actions
1 Safety flap missing or broken.	Please contact your Agilent service representative.

Valve to Bypass Failed

Error ID: 4014, 4701

The injection valve failed to switch to the bypass position.

The switching of the injection valve is monitored by two microswitches on the valve assembly. The switches detect the successful completion of the valve movement. If the valve fails to reach the bypass position, or if the microswitch does not close, the error message is generated.

10 Error Information

Autosampler Error Messages

Probable cause	Suggested actions
1 Defective injection valve.	Please contact your Agilent service representative.
2 Defective main board.	Please contact your Agilent service representative.

Valve to Mainpass Failed

Error ID: 4015

The injection valve failed to switch to the mainpass position.

The switching of the injection valve is monitored by two microswitches on the valve assembly. The switches detect the successful completion of the valve movement. If the valve fails to reach the mainpass position, or if the microswitch does not close, the error message is generated.

Probable cause	Suggested actions
1 Defective injection valve.	Please contact your Agilent service representative.
2 Defective main board.	Please contact your Agilent service representative.

Vial in Gripper

Error ID: 4033

The gripper arm attempted to move with a vial still in the gripper.

During specific stages of the sampling sequence, no vial should be held by the gripper. The autosampler checks if a sample vial is stuck in the gripper by closing and opening the gripper fingers while monitoring the motor encoder. If the gripper fingers are unable to close, the error message is generated.

Probable cause	Suggested actions
1 Vial still in gripper.	Remove the vial using the Release Vial function in the user interface. Reinitialize the autosampler.

10 Error Information

General Detector Error Messages

General Detector Error Messages

These errors are specific to both detector types VWD and DAD.

Heater at fan assembly failed

Error ID: 1073

Every time the deuterium lamp or the tungsten lamp (DAD only) is switched on or off a heater self-test is performed. If the test fails an error event is created. As a result the temperature control is switched off.

Probable cause	Suggested actions
1 Defective connector or cable.	Please contact your Agilent service representative.
2 Defective heater.	Please contact your Agilent service representative.

Heater Power At Limit

Error ID: 1074

The available power of the heater reached either the upper or lower limit. This event is sent only once per run. The parameter determines which limit has been hit:

0 means upper power limit hit (excessive ambient temperature drop).

1 means lower power limit hit (excessive ambient temperature increase).

Probable cause	Suggested actions
1 Excessive ambient temperature change.	Wait until temperature control equilibrates.

Illegal Temperature Value from Sensor on Main Board

Error ID: 1071

This temperature sensor (located on the detector main board) delivered a value outside the allowed range. The parameter of this event equals the measured temperature in 1/100 centigrade. As a result the temperature control is switched off.

Probable cause	Suggested actions
1 Defective sensor or main board.	Please contact your Agilent service representative.
2 Detector is exposed to illegal ambient conditions.	Verify that the ambient conditions are within the allowed range.

Illegal Temperature Value from Sensor at Air Inlet

Error ID: 1072

This temperature sensor delivered a value outside the allowed range. The parameter of this event equals the measured temperature in 1/100 centigrade. As a result the temperature control is switched off.

Probable cause	Suggested actions
1 The temperature sensor is defect.	<ul style="list-style-type: none">Replace the cable to the main board.Please contact your Agilent service representative.
2 Detector is exposed to illegal ambient conditions.	Verify that the ambient conditions are within the allowed range.

UV Lamp Current

Error ID: 7450

The UV lamp current is missing.

10 Error Information

General Detector Error Messages

The processor continually monitors the anode current drawn by the lamp during operation. If the anode current falls below the lower current limit, the error message is generated.

Probable cause	Suggested actions
1 Lamp disconnected.	Ensure the UV lamp connector is seated firmly.
2 Defective UV lamp or non-Agilent lamp.	Exchange the UV lamp.
3 Defective detector main board.	Please contact your Agilent service representative.
4 Defective power supply.	Please contact your Agilent service representative.

UV Lamp Voltage

Error ID: 7451

The UV lamp anode voltage is missing.

The processor continually monitors the anode voltage across the lamp during operation. If the anode voltage falls below the lower limit, the error message is generated.

Probable cause	Suggested actions
1 Defective UV lamp or non-Agilent lamp.	Exchange the UV lamp.
2 Defective detector main board.	Please contact your Agilent service representative.
3 Defective power supply.	Please contact your Agilent service representative.

VWD Detector Error Messages

These errors are specific to the variable wavelength detector.

ADC Hardware Error

Error ID: 7830, 7831

A/D-Converter hardware is defective.

Probable cause	Suggested actions
1 A/D-Converter hardware is defective.	Please contact your Agilent service representative.

Wavelength calibration setting failed

Error ID: 7310

The intensity maximum was not found during wavelength calibration.

Calibration 0 Failed: Zero-order calibration failed.

Calibration 1 Failed: 656 nm calibration failed.

Probable cause	Suggested actions
1 Lamp is OFF.	Switch on the lamp.
2 Incorrect flow cell installation.	Ensure the flow cell is installed correctly.
3 Flow cell contamination or air bubbles.	Clean/replace flow cell windows or remove air bubbles.
4 Intensity too low.	Replace lamp.

10 Error Information

VWD Detector Error Messages

Probable cause	Suggested actions
5 Current step value too far from maximum.	<ul style="list-style-type: none">• Repeat the calibration.• Please contact your Agilent service representative.
6 Misaligned/defective grating assembly.	Please contact your Agilent service representative.
7 Defective main board.	Please contact your Agilent service representative.

Cutoff filter doesn't decrease the light intensity at 250 nm

Error ID: 7813

The automatic filter check after lamp ignition has failed. When the lamp is switched on, the detector moves the cutoff filter into the light path. If the filter is functioning correctly, a decrease in lamp intensity is seen. If the expected intensity decrease is not detected, the error message is generated.

Probable cause	Suggested actions
1 Motor is not connected.	Please contact your Agilent service representative.
2 Defective motor.	Please contact your Agilent service representative.
3 Defective/missing grating or filter.	Please contact your Agilent service representative.
4 Cable/connector defective.	Please contact your Agilent service representative.

Filter Missing

Error ID: 7816

The filter motor is not detected.

Probable cause	Suggested actions
1 Filter motor is not connected.	Please contact your Agilent service representative.
2 Cable/connector defective.	Please contact your Agilent service representative.

Grating or Filter Motor Errors

Error ID: Grating: 7800, 7801, 7802, 7803, 7804, 7805, 7806, 7808, 7809;
Filter: 7810, 7811, 7812, 7813, 7814, 7815, 7816

The motor test has failed.

Test 0 Failed:

Filter motor.

Test 1 Failed:

Grating motor.

During the motor tests, the detector moves the motor to the end position while monitoring the end-position sensor. If the end position is not found, the error message is generated.

Probable cause

1 Motor is not connected.

Suggested actions

Please contact your Agilent service representative.

2 Defective motor.

Please contact your Agilent service representative.

3 Defective/missing grating or filter.

Please contact your Agilent service representative.

4 Cable/connector defective.

Please contact your Agilent service representative.

10 Error Information

VWD Detector Error Messages

Grating Missing

Error ID: 7819

The grating motor is not detected.

Probable cause	Suggested actions
----------------	-------------------

- | | |
|-----------------------------------|---|
| 1 Grating motor is not connected. | Please contact your Agilent service representative. |
| 2 Cable/connector defective. | Please contact your Agilent service representative. |

No heater current

Error ID: 7453

The lamp heater current in the detector is missing. During lamp ignition, the processor monitors the heater current. If the current does not rise above the lower limit within 1, the error message is generated.

Probable cause	Suggested actions
----------------	-------------------

- | | |
|---|---|
| 1 Lamp disconnected. | Ensure the lamp is connected. |
| 2 Ignition started without the top foam in place. | Please contact your Agilent service representative. |
| 3 Fan not running (permitting lamp on). | Please contact your Agilent service representative. |
| 4 Defective main board. | Please contact your Agilent service representative. |
| 5 Defective or non-Agilent lamp. | Exchange the lamp. |
| 6 Defective power supply. | Please contact your Agilent service representative. |

Wavelength holmium check failed

Error ID: 7318

The holmium oxide test in the detector has failed. During the holmium test, the detector moves the holmium filter into the light path, and compares the measured absorbance maxima of the holmium oxide filter with expected maxima. If the measured maxima are outside the limits, the error message is generated.

Probable cause	Suggested actions
1 Misaligned/defective grating assembly.	<ul style="list-style-type: none">• Ensure the flow cell is inserted correctly, and is free from contamination (cell windows, buffers, and so on).• Run the filter-motor test to determine if the filter motor assembly is defective. If defective, please contact your Agilent service representative.• Run the grating-motor test to determine if the grating assembly is defective. If defective, please contact your Agilent service representative.

Ignition Failed

Error ID: 7452

The lamp failed to ignite. The processor monitors the lamp current during the ignition cycle. If the lamp current does not rise above the lower limit within 2 – 5 s, the error message is generated.

Probable cause	Suggested actions
1 Lamp disconnected.	Ensure the lamp is connected.
2 Defective or non-Agilent lamp.	Exchange the lamp.

10 Error Information

VWD Detector Error Messages

Probable cause	Suggested actions
3 Defective power supply.	Please contact your Agilent service representative.
4 Defective main board.	Please contact your Agilent service representative.

Wavelength test failed

Error ID: 7890

The automatic wavelength check after lamp ignition has failed. When the lamp is switched on, the detector waits 1 min to warm-up the lamp. Then a check of the deuterium emission line (656 nm) via the reference diode is performed. If the emission line is more than 3 nm away from 656 nm, the error message is generated.

Probable cause	Suggested actions
1 Calibration incorrect.	Recalibrate the detector.

DAD Detector Error Messages

These errors are specific to the diode array detector.

Visible Lamp Current

The visible lamp current is missing.

The processor continually monitors the lamp current during operation. If the current falls below the lower current limit, the error message is generated.

Probable cause	Suggested actions
1 Lamp disconnected.	Ensure the visible lamp connector is seated firmly.
2 Defective visible lamp.	Exchange the visible lamp.
3 Defective connector or cable.	Please contact your Agilent service representative.
4 Defective power supply.	Please contact your Agilent service representative.

Visible Lamp Voltage

The visible lamp voltage is missing.

The processor continually monitors the voltage across the lamp during operation. If the lamp voltage falls below the lower limit, the error message is generated.

10 Error Information

DAD Detector Error Messages

Probable cause	Suggested actions
1 Defective connector or cable.	Please contact your Agilent service representative.
2 Defective power supply.	Please contact your Agilent service representative.

Diode Current Leakage

Error ID: 1041

When the detector is switched on, the processor checks the leakage current of each of the optical diodes. If the leakage current exceeds the upper limit, the error message is generated.

Probable cause	Suggested actions
1 Defective PDA/optical unit.	Please contact your Agilent service representative.
2 Defective connector or cable.	Please contact your Agilent service representative.

UV Ignition Failed

Error ID: 7452

The UV lamp failed to ignite.

The processor monitors the UV lamp current during the ignition cycle. If the lamp current does not rise above the lower limit within 2 – 5 seconds, the error message is generated.

Probable cause	Suggested actions
1 Lamp too hot. Hot gas discharge lamps may not ignite as easily as cold lamps.	Switch off the lamp and allow it to cool down for at least 15 minutes.
2 Lamp disconnected.	Ensure the lamp is connected.
3 Defective UV lamp or non-Agilent lamp.	Exchange the UV lamp.
4 Defective detector main board.	Please contact your Agilent service representative.
5 Defective power supply.	Please contact your Agilent service representative.

UV Heater Current

Error ID: 7453

The UV lamp heater current is missing.

During UV lamp ignition, the processor monitors the heater current. If the current does not rise above the lower limit within one second, the error message is generated.

Probable cause	Suggested actions
1 Lamp disconnected.	Ensure the UV lamp is connected.
2 Ignition started without the top foam in place.	Please contact your Agilent service representative.
3 Defective UV lamp or non-Agilent lamp.	Exchange the UV lamp.
4 Defective detector main board.	Please contact your Agilent service representative.
5 Defective power supply.	Please contact your Agilent service representative.

10 Error Information

DAD Detector Error Messages

Calibration Values Invalid

Error ID: 1036

The calibration values read from the spectrometer ROM are invalid.

After recalibration, the calibration values are stored in ROM. The processor periodically checks if the calibration data are valid. If the data are invalid or cannot be read from the spectrometer ROM, the error message is generated.

Probable cause	Suggested actions
1 Defective connector or cable.	Please contact your Agilent service representative.
2 Defective PDA/optical unit.	Please contact your Agilent service representative.

Holmium Oxide Test Failed

Probable cause	Suggested actions
1 Lamps switched off.	Ensure the lamps are switched on.
2 Defective or dirty flow cell.	Ensure the flow cell is inserted correctly, and is free from contamination (cell windows, buffers etc.).
3 Defective filter assembly.	Please contact your Agilent service representative.
4 Defective achromat assembly.	Please contact your Agilent service representative.
5 Defective PDA/optical unit.	Please contact your Agilent service representative.

Wavelength Recalibration Lost

Error ID: 1037

The calibration information needed for your detector to operate correctly has been lost.

During calibration of the detector the calibration values are stored in ROM. If no data is available in the spectrometer ROM, the error message is generated.

Probable cause	Suggested actions
1 The detector is new.	Recalibrate the detector.
2 The detector has been repaired.	Please contact your Agilent service representative.

DSP Not Running

This error message comes up when the communication between the optical unit and the main board has a problem.

Probable cause	Suggested actions
1 Random communication error.	<ul style="list-style-type: none">Switch the detector off and on again at the power switch. If the error reoccurs:Please contact your Agilent service representative.
2 Defective detector main board.	Please contact your Agilent service representative.
3 Defective PDA/optical unit.	Please contact your Agilent service representative.

No Run Data Available In Device

In a very rare case the capacity of the CompactFlash Card is not sufficient. This could happen for example when the interrupt of LAN communication

10 Error Information

DAD Detector Error Messages

takes longer and the detector uses special settings (e.g full data rate at 80 Hz plus full spectra plus all signals) during data buffering.

Probable cause	Suggested actions
1 CompactFlash Card is full.	<ul style="list-style-type: none">Correct communication problem.Reduce data rate.

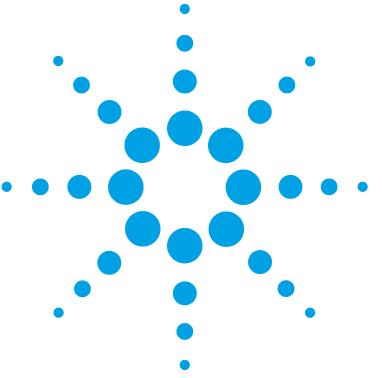
Instrument Logbook

Method	Instrument run started	09:44:46 11/20/05
1200 DAD	1 Power on	10:07:24 11/20/05
1200 DAD	1 UV-lamp on	10:07:24 11/20/05
1200 DAD	1 Vis-lamp on	10:07:24 11/20/05
1200 DAD	1 No Run data available in device!	10:07:24 11/20/05
CP Macro	Analyzing rawdata SHORT_02.D	10:07:25 11/20/05
Method	Instrument Error - Method/Sequence stopped	10:07:25 11/20/05
Method	Method aborted	10:09:52 11/20/05

Figure 63 Instrument Logbook

NOTE

The logbook does not indicate a communication loss (power fail). It just shows the recovering (Power on, Lamps on).



11 Maintenance and Repair

PM Scope of Work and Checklist	243
Cautions and Warnings	244
Solvent Delivery System	246
Introduction	246
Checking and Cleaning the Solvent Filter	248
Exchanging the Passive Inlet Valve	250
Exchanging the Outlet Valve	253
Exchanging the Purge Valve Frit or the Purge Valve	254
Removing the Pump Head Assembly	257
Maintenance of a Pump Head	259
Reinstalling the Pump Head Assembly	262
Exchanging the Plungers	263
Exchanging the Dual-Channel Gradient Valve (DCGV)	264
Manual Injector	267
Overview of Maintenance Procedures	267
Flushing the Manual Injector	267
Exchanging the Injection Valve Seal	268
Autosampler	271
Introduction	271
Exchanging the Needle Assembly	274
Exchanging the Needle Seat Assembly	278
Exchanging the Rotor Seal	280
Exchanging the Metering Seal	284
Exchanging the Gripper Arm	288
Variable Wavelength Detector (VWD)	290
Introduction	290
Exchanging the Deuterium Lamp	291
Replace the Flow Cell / Cuvette Holder	294



11 Maintenance and Repair

DAD Detector Error Messages

Repairing the Flow Cell	296
Using the Cuvette Holder	299
Correcting Leaks	301
Diode Array Detector (DAD)	302
Overview of Maintenance	302
Cleaning the Module	303
Exchanging a Lamp	304
Remove and Install a Flow Cell	307
Maintenance of the Flow Cell	310
Replacing Capillaries on a Standard Flow Cell	313
Cleaning or Exchanging the Holmium Oxide Filter	318
Correcting Leaks	321
Replacing Leak Handling System Parts	322
Algae Growth in HPLC Systems	323
Replacing the Module's Firmware	325

This chapter provides general information on maintenance and repair of the instrument.

PM Scope of Work and Checklist

Preventive Maintenance Scope of Work and Checklist

To perform a preventive maintenance (PM), follow the *PM Scope of Work* and *PM Checklist* step by step. The *PM Scope of Work* and *PM Checklist* documents can be found in the on the DVD for the Lab Advisor Software.

11 Maintenance and Repair

Cautions and Warnings

Cautions and Warnings

WARNING

Module is partially energized when switched off, as long as the power cord is plugged in.

Risk of stroke and other personal injury. Repair work at the module can lead to personal injuries, e. g. shock hazard, when the module cover is opened and the instrument is connected to power.

- Never perform any adjustment, maintenance or repair of the module with the top cover removed and with the power cord plugged in.
 - The security lever at the power input socket prevents that the module cover is taken off when line power is still connected. Never plug the power line back in when cover is removed.
-

WARNING

Sharp metal edges

Sharp-edged parts of the equipment may cause injuries.

- To prevent personal injury, be careful when getting in contact with sharp metal areas.
-

WARNING

Toxic, flammable and hazardous solvents, samples and reagents

The handling of solvents, samples and reagents can hold health and safety risks.

- When working with these substances observe appropriate safety procedures (for example by wearing goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the vendor, and follow good laboratory practice.
 - The volume of substances should be reduced to the minimum required for the analysis.
 - Do not operate the instrument in an explosive atmosphere.
-

CAUTION

Electronic boards and components are sensitive to electrostatic discharge (ESD). ESD can damage electronic boards and components.

- Be sure to hold the board by the edges, and do not touch the electrical components. Always use ESD protection (for example, an ESD wrist strap) when handling electronic boards and components.
-

WARNING**Eye damage by detector light**

Eye damage may result from directly viewing the UV-light produced by the lamp of the optical system used in this product.

- Always turn the lamp of the optical system off before removing it.
-

CAUTION**Safety standards for external equipment**

- If you connect external equipment to the instrument, make sure that you only use accessory units tested and approved according to the safety standards appropriate for the type of external equipment.
-

11 Maintenance and Repair

Solvent Delivery System

Solvent Delivery System

Introduction

The Agilent 1220 Infinity II LC pump is designed for easy repair. The procedures described in this section can be done with the pump in place in the rack.

The most frequent repairs, such as exchanging the plunger seals or purge valve seal, can be done from the front of the pump.

Table 44 Simple repair procedures - overview

Procedure	Typical frequency	Notes
Checking and cleaning the solvent filter	If solvent filter is blocked	Gradient performance problems, intermittent pressure fluctuations
Exchanging the Passive Inlet Valve	If internally leaking	Pressure ripple unstable, run Leak Test for verification
Exchanging the Outlet Ball Valve	If internally leaking	Pressure ripple unstable, run Leak Test for verification
Exchanging the Purge Valve Frit or the Purge Valve	If internally leaking	Solvent dripping out of waste outlet when valve closed
Exchanging the Purge Valve Frit or the Purge Valve	If the frit shows indication of contamination or blockage	A pressure drop of > 10 bar across the frit (5 mL/min H ₂ O with purge open) indicates blockage
Exchanging the Pump Seals	If pump performance indicates seal wear	Leaks at lower pump head side, unstable retention times, pressure ripple unstable — run Leak Test for verification
Seal Wear-in Procedure	After exchanging the pump seals	

Table 44 Simple repair procedures - overview

Procedure	Typical frequency	Notes
Removing the pump head assembly	Before exchanging the seals, or plungers	
Exchanging the Plungers	If scratched	Seal life time shorter than normally expected — check plungers while changing the seals

WARNING

Instrument is partially energized when switched off.

The power supply still uses some power, even if the switch on the front panel is turned off.

- To disconnect the Agilent 1220 Infinity II LC pump from line, unplug the power cord.

WARNING

Sharp metal edges

Sharp-edged parts of the equipment may cause injuries.

- To prevent personal injury, be careful when getting in contact with sharp metal areas.

WARNING

When opening capillary or tube fittings, solvents may leak out.

The handling of toxic and hazardous solvents and reagents can carry health risks.

- Observe appropriate safety procedures (for example, wear goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the solvent vendor, especially when toxic or hazardous solvents are used.

CAUTION

Electronic boards and components are sensitive to electrostatic discharge (ESD).

ESD can damage electronic boards and components.

- In order to prevent damage always use an ESD protection when handling electronic boards and components.

Checking and Cleaning the Solvent Filter

WARNING

Small particles can permanently block the capillaries and valves of the pump.

Damage to the Agilent 1220 Infinity II LC pump

- Always filter solvents.
 - Never use the pump without solvent inlet filter.
-

NOTE

If the filter is in good condition the solvent will freely drip out of the solvent tube (hydrostatic pressure). If the solvent filter is partly blocked only very little solvent will drip out of the solvent tube.

Cleaning the Solvent Filter

When If solvent filter is blocked

Parts required

Description
Concentrated nitric acid (65 %)
Bidistilled water
Beaker

Preparations Remove solvent inlet tube from the inlet valve.

- 1** Remove the blocked solvent filter from the bottle-head assembly and place it in a beaker with concentrated nitric acid (65 %) for one hour.
- 2** Thoroughly flush the filter with LC grade water (remove all nitric acid, some columns can be damaged by concentrated nitric acid; check with pH indicator).
- 3** Reinstall the filter.

Exchanging the Passive Inlet Valve

When If leaking internally (backflow)

Tools required

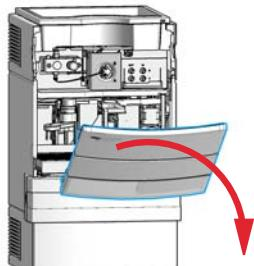
Description
Wrench, 14 mm
Pair of tweezers

Parts required

p/n	Description
G1312-60066	Passive inlet valve 1220/1260

Preparations Place the solvent bottles under the pump.

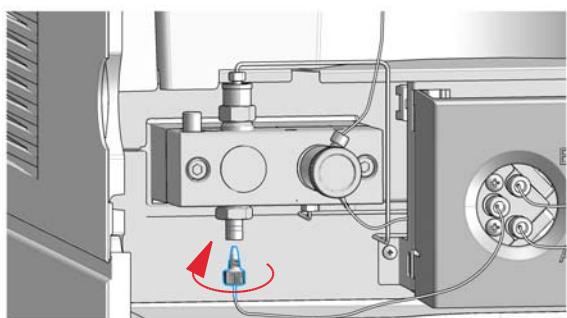
- 1** Remove the upper front cover.



- 2** Disconnect the solvent inlet tube from the inlet valve.

NOTE

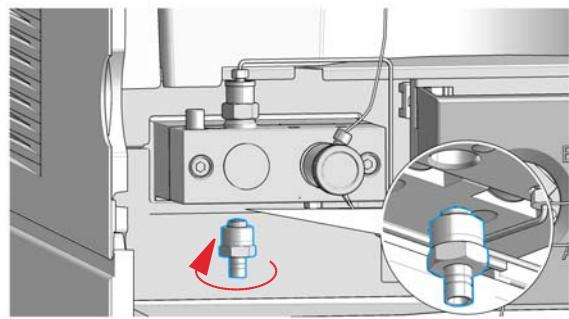
Beware of leaking solvents due to hydrostatic pressure.



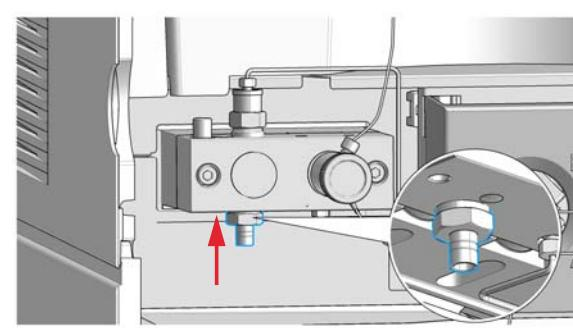
3 Unscrew the adapter from the inlet valve (optional).



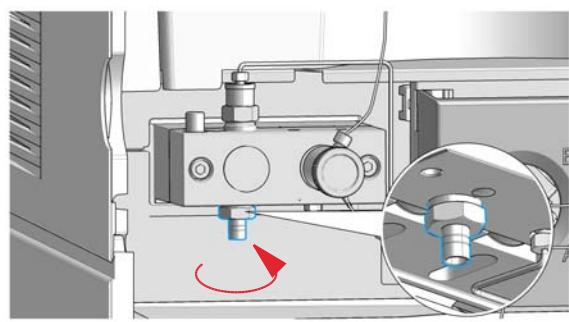
4 Using a 14 mm wrench, loosen the passive inlet valve and remove the valve from the pump head.



5 Insert the new valve into the pump head.



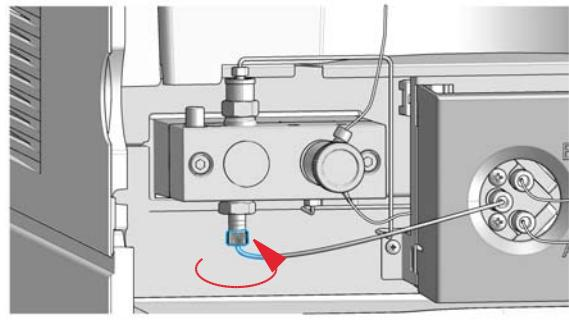
6 Using the 14 mm wrench, turn the nut until it is hand-tight.



7 Reconnect the adapter at the inlet valve (optional).



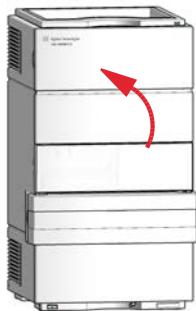
8 Reconnect the solvent inlet tube to the adapter.



11 Maintenance and Repair

Solvent Delivery System

9 Close the upper front cover.



Exchanging the Outlet Valve

When If leaking internally

Tools required

Description

Wrench, 14 mm

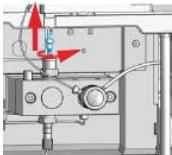
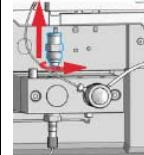
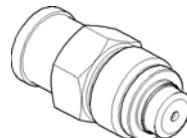
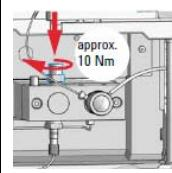
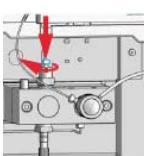
Parts required

p/n **Description**

G1312-60067 Outlet valve 1220/1260

Preparations Switch off pump at the main power switch

Remove the upper front cover

<p>1 Using a 1/4 inch wrench, disconnect the valve capillary from the outlet valve.</p> 	<p>2 Using the 14 mm wrench, loosen the valve and remove it from the pump body.</p> 
<p>3 Check that the new valve is assembled correctly and that the gold seal is present (if the gold seal is deformed, it should be replaced).</p> 	<p>4 Reinstall the outlet valve and tighten the valve.</p> 
<p>5 Reconnect the valve capillary.</p> 	

Exchanging the Purge Valve Frit or the Purge Valve

When

Frit: when plunger seals are exchanged or when contaminated or blocked (pressure drop of > 10 bar across the frit at a flow rate of 5 mL/min of H₂O with purge valve opened)

Purge valve: if internally leaking

Tools required

Wrench, 1/4 inch
Wrench, 14 mm
Pair of tweezers
OR
Toothpick

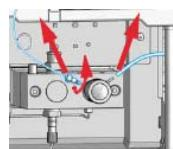
Parts required

#	p/n	Description
1	01018-22707	PTFE frits (pack of 5)
1	G4280-60061	Purge valve

Preparations

Switch off pump at the main power switch
Remove the upper front cover

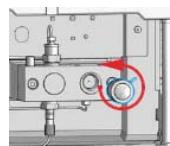
- 1 Using a 1/4 inch wrench, disconnect the pump outlet capillary at the purge valve.



NOTE

Beware of leaking solvents due to hydrostatic pressure.

- 2 Using the 14 mm wrench, unscrew the purge valve and remove it.



- 3 Remove the seal cap from the purge valve.

- Using a pair of tweezers or a toothpick, remove the frit.

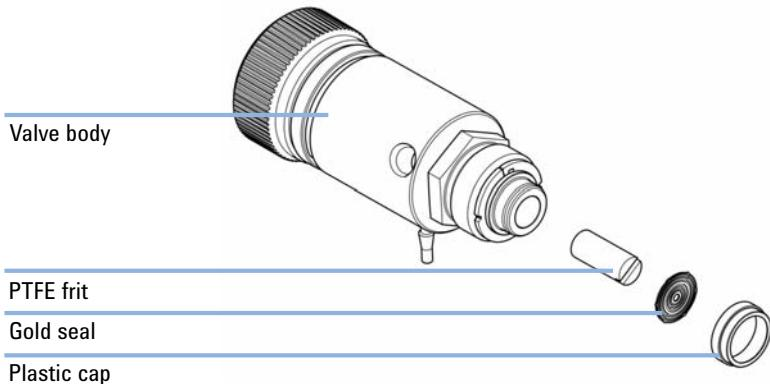


Figure 64 Purge Valve Parts

- Place a new frit into the purge valve, with the orientation of the frit as shown above.
- Reinstall the seal cap including the gold seal.

NOTE

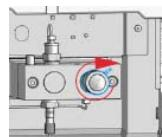
Before reinstallation always check the gold seal in the seal cap. A deformed seal cap should be exchanged.

CAUTION

Damage to the purge valve

- Do not lift the pump using the purge valve as a handle, it might get leaky.
- Do not try to turn the purge valve into the correct position when already fixed to the pump. The rubber o-ring might break.
- Anticipate the correct position of the connections before tightening the valve.

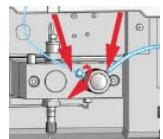
- Insert the purge valve into the pump head and fix it with a 14 mm wrench.



11 Maintenance and Repair

Solvent Delivery System

- 8 Reconnect the outlet capillary and the waste tubing.



Removing the Pump Head Assembly

When
Before exchanging the seals
Before exchanging the plungers

Tools required

p/n	Description
8710-0510	Wrench open 1/4 — 5/16 inch
8710-2392	Hexagonal key, 4.0 mm, 15 cm long, T-handle

Preparations
Switch off pump at the main power switch

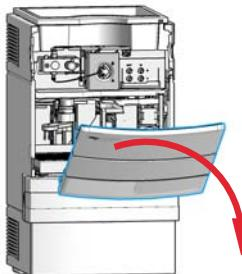
CAUTION

Damage of the pump drive

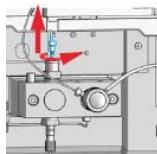
Starting the pump when the pump head is removed may damage the pump drive.

→ Never start the pump when the pump head is removed.

- 1** Remove the upper front cover.



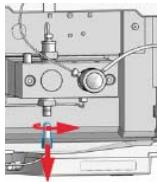
- 2** Using a 1/4 inch wrench disconnect the valve capillary from the outlet valve.



11 Maintenance and Repair

Solvent Delivery System

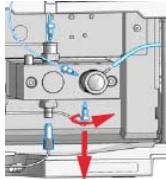
- 3** Disconnect the solvent inlet tube from the inlet valve.



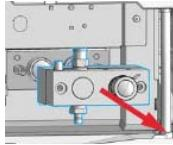
NOTE

Beware of leaking solvents due to hydrostatic pressure.

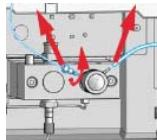
- 5** Remove the capillary at the bottom of the pump head.



- 7** Remove the pump head from the pump drive.



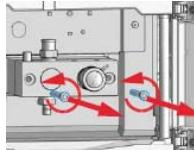
- 4** Using a 1/4 inch wrench disconnect the pump outlet capillary from the purge valve. Disconnect the waste tube.



NOTE

Beware of leaking solvents due to hydrostatic pressure.

- 6** Using a 4 mm hexagonal key, stepwise loosen the two pump head screws.



Maintenance of a Pump Head

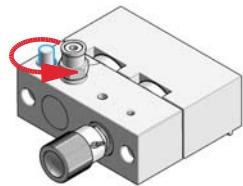
When Seal leaking, if indicated by the results of the leak test.

Tools required	p/n	Description
	8710-0510	Wrench open 1/4 — 5/16 inch
	8710-2392	Hexagonal key, 4.0 mm, 15 cm long, T-handle

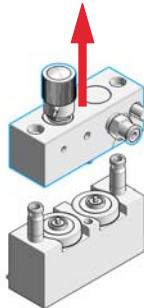
Parts required	p/n	Description
	5063-6589	Piston seal PTFE, carbon filled, black (pack of 2), default
	0905-1420	PE seals (pack of 2)
	5022-2159	Restriction capillary

Preparations Switch off the pump at the main power switch.
Remove the upper front cover.

- 1** Place the pump head on a flat surface. Loosen the lock screw (two turns).



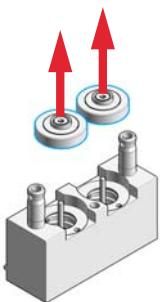
- 2** While holding the lower half of the assembly (piston housing), carefully pull the pump housing away from the piston housing.



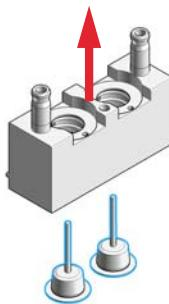
11 Maintenance and Repair

Solvent Delivery System

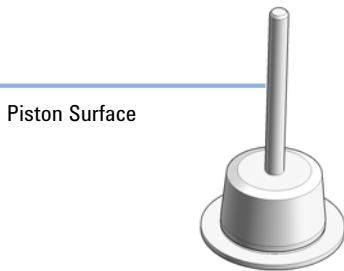
- 3** Remove the support rings from the piston housing.



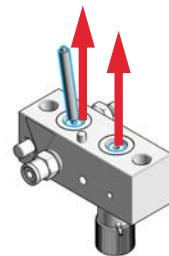
- 4** Lift the housing away from the pistons.



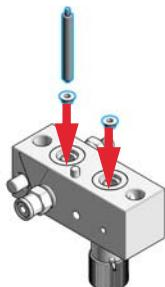
- 5** Check the piston surface and remove any deposits or layers: clean the piston surface with abrasive paper and rinse with 2-propanol. Replace piston if scratched.



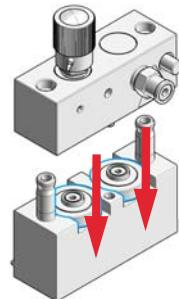
- 6** Using the steel side of the insert tool, carefully remove the seal from the pump housing.



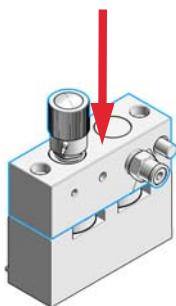
- 7** Using the plastic side of the insert tool, insert new seals into the pump head.



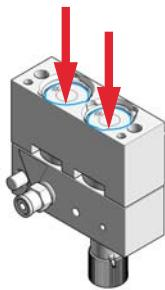
- 8** Place the support rings on the piston housing. Note the correct position of the pins.



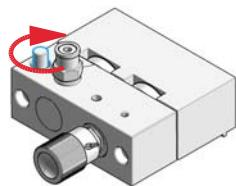
9 Place the pump housing onto the piston housing.



10 Insert the pistons and carefully press them into the seals.



11 Tighten the lock screw.



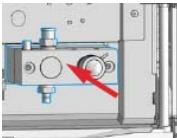
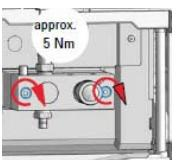
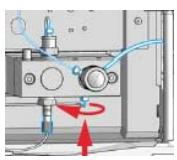
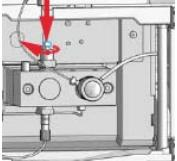
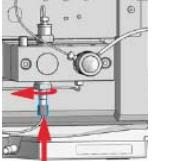
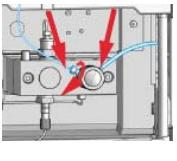
12 Install the pump head, see “[Reinstalling the Pump Head Assembly](#)” on page 262.

Reinstalling the Pump Head Assembly

When When reassembling the pump

Tools required **Description**

Hexagonal key, 4 mm

<p>1 If needed, apply a small amount of grease on the back of the screws. Normally, the grease added during manufacturing is sufficient for a long time.</p> 	<p>2 Slide the pump head assembly onto the pump drive.</p>
<p>3 Insert the pump head screws and tighten them stepwise with increasing torque (approx. 5 Nm) with a 4 mm hexagonal key.</p> 	<p>4 Install the capillary at the bottom of the pump head.</p> 
<p>5 Reconnect the valve capillary to the outlet valve.</p> 	<p>6 Reconnect the solvent inlet tube to the passive inlet valve.</p> 
<p>7 Reconnect the outlet capillary and the waste tubing.</p> 	<p>Next Steps:</p> <p>8 If installed, reconnect the active inlet valve cable to its connector.</p> <p>9 If a standard seal has been installed, run the seal wear-in procedure, which includes a replacement of the purge valve frit, see "Exchanging the Plungers" on page 263.</p> <p>10 For the normal phase seal, the purge valve frit should be replaced, see "Exchanging the Purge Valve Frit or the Purge Valve" on page 254.</p>

Exchanging the Plungers

CAUTION

This procedure is required for Standard seals (pack of 2) (5063-6589) only.

It will damage the PE seals (pack of 2) (0905-1420).

→ Never perform the seal wear-in procedure with normal-phase application seals.

- 1 Place a bottle with 100 mL of isopropanol in the solvent cabinet and attach a bottle-head assembly (including tubing) to the bottle.
- 2 Connect the inlet tube from the bottle head directly to the inlet valve.
- 3 Connect one end of the Restriction capillary (5022-2159) to the purge valve. Insert the other end into a waste container.
- 4 Open the purge valve and purge the system for 5 min with isopropanol at a flow rate of 2 mL/min.
- 5 Close the purge valve and set the flow to a rate adequate to achieve a pressure of 350 bar.
- 6 Pump for 15 min at this pressure to wear in the seals.
- 7 Turn off the pump and slowly open the purge valve to release the pressure from the system.
- 8 Disconnect the restriction capillary and reinstall the bottle containing the solvent for your application.
- 9 Rinse your system with the solvent used for your application.

Exchanging the Dual-Channel Gradient Valve (DCGV)

Tools required	p/n	Description
	8710-0899	Screwdriver, Pozidriv #1

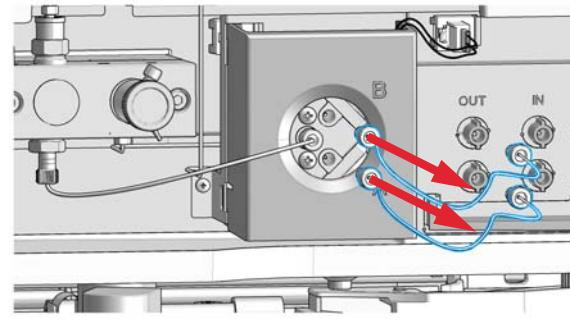
Parts required	#	p/n	Description
	1	G4280-60004	Dual-channel gradient valve

Preparations Switch off the pump at the power switch.
Remove the upper front cover to gain access to the pump mechanics.

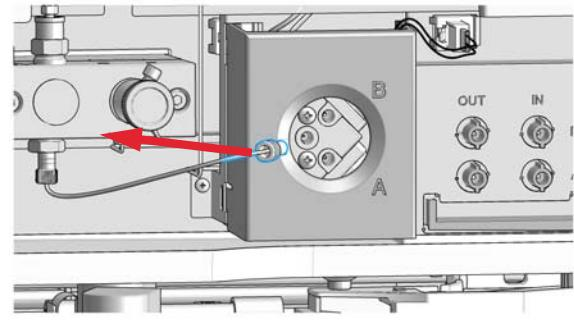
NOTE

The lifetime of the dual-channel gradient valve can be maintained by regularly flushing the valve, especially when using buffer solutions. If using buffer solutions, flush all channels of the valve with water to prevent precipitation of the buffer. Salt crystals can be forced into an unused channel and form plugs that may lead to leaks of that channel; such leaks will interfere with the general performance of the valve. When using buffer solutions and organic solvents in the Agilent 1220 Infinity II LC Pump, it is recommended to connect the buffer solution to the lower port of the gradient valve and the organic solvent to the upper port. It is best to have the organic channel directly above the salt solution channel (A: salt solution, B: organic solvent).

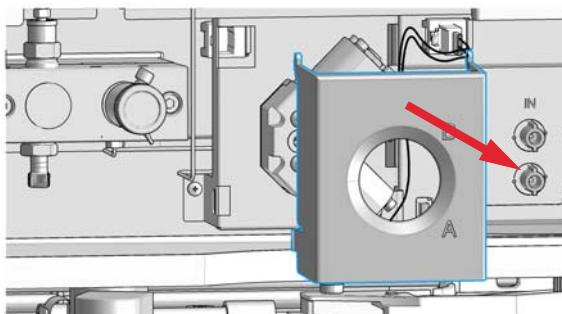
- 1** Disconnect the solvent tubes from the DCGV.



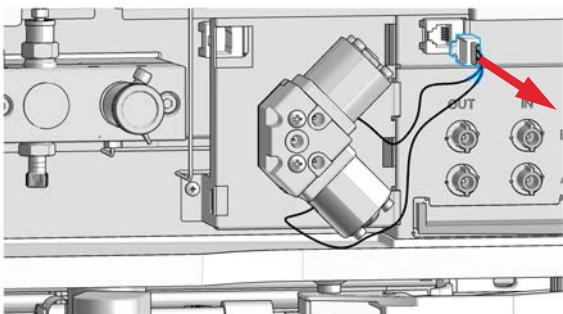
- 2** Disconnect the connecting tube, waste tube and the solvent tubes from the DCGV, unclip them from the tube clips and place them into the solvent cabinet to avoid flow by hydrostatic pressure.



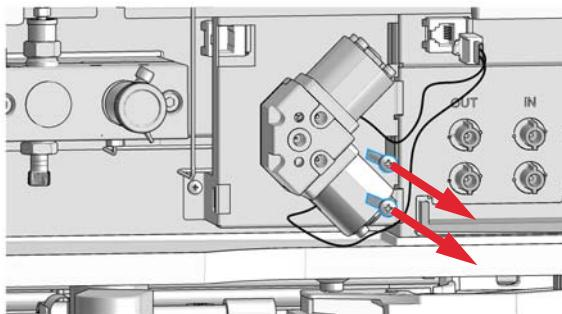
- 3** Press the lower sides of the cover to unclip it. Remove the cover.



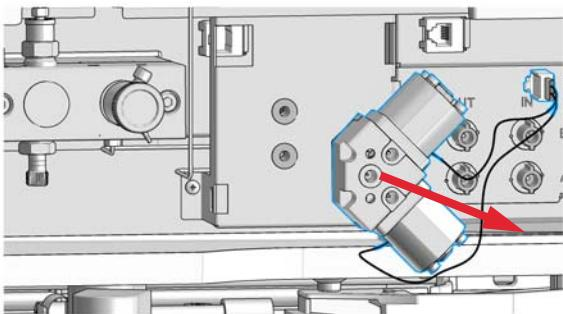
- 4** Disconnect the DCGV cable.



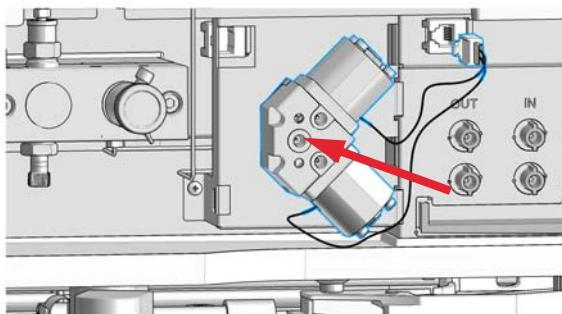
- 5** Unscrew the two holding screws.



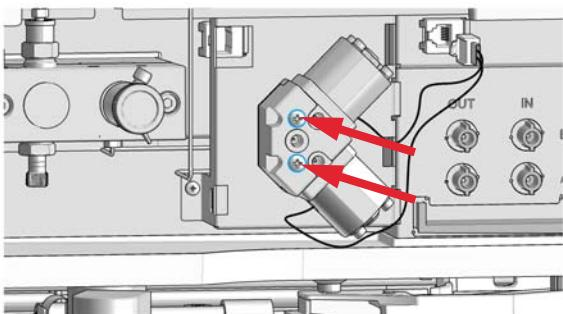
- 6** Remove the DCGV.



- 7** Place the new DCGV into position. Make sure that the valve is positioned with the A-channel at the bottom-right position. Tighten the two holding screws and connect the cable to its connector.



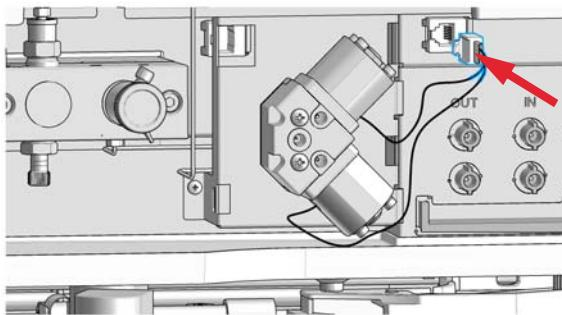
- 8** Tighten the two holding screws.



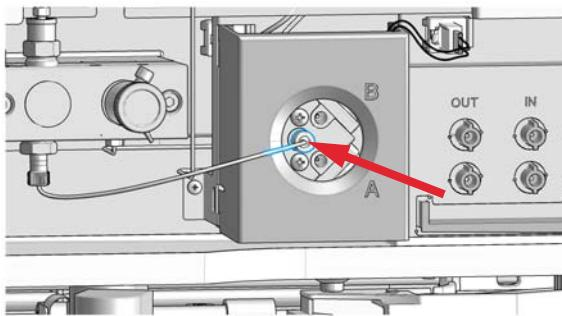
11 Maintenance and Repair

Solvent Delivery System

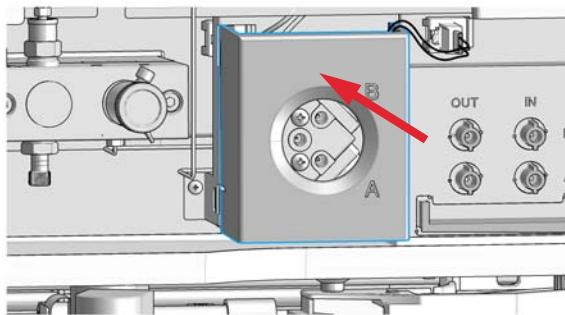
9 Connect the cable to its connector



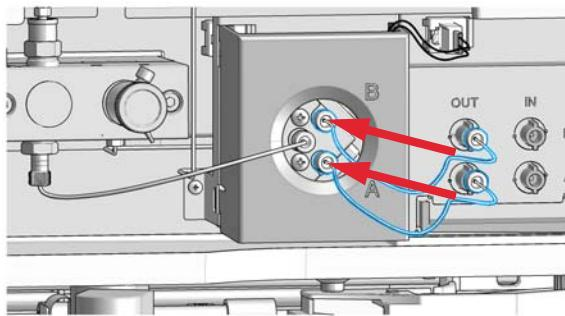
11 Reconnect the tube from the inlet valve to the middle position of the DCGV.



10 Replace the DCGV cover. Reconnect the waste funnel with the waste tube holder in the top cover. Insert waste tube in the holder in the waste pan and clip tube to the DCGV cover.



12 Connect the solvent tubes at channel A and B of the DCGV.



Manual Injector

Overview of Maintenance Procedures

Table 45 Overview of maintenance procedures

Procedure	Typical frequency	Time required
Flushing the injector	After using aqueous buffers or salt solutions	5 min
Exchanging the injection-valve seal	After approximately 10000 to 20000 injections, or when the valve performance shows indication of leakage or wear	10 min

Flushing the Manual Injector

CAUTION

The use of aqueous buffers or salt solutions can lead to crystal formation.

Crystal formation may cause scratches on the injection seal.

→ Always rinse the valve with water after using aqueous buffers or salt solutions.

- 1 Switch the valve to the INJECT position.
- 2 Use the pump to flush the sample loop and seal grooves.
- 3 Use the needle-port cleaner (supplied with the valve) and syringe to flush the needle port and vent capillary.

11 Maintenance and Repair

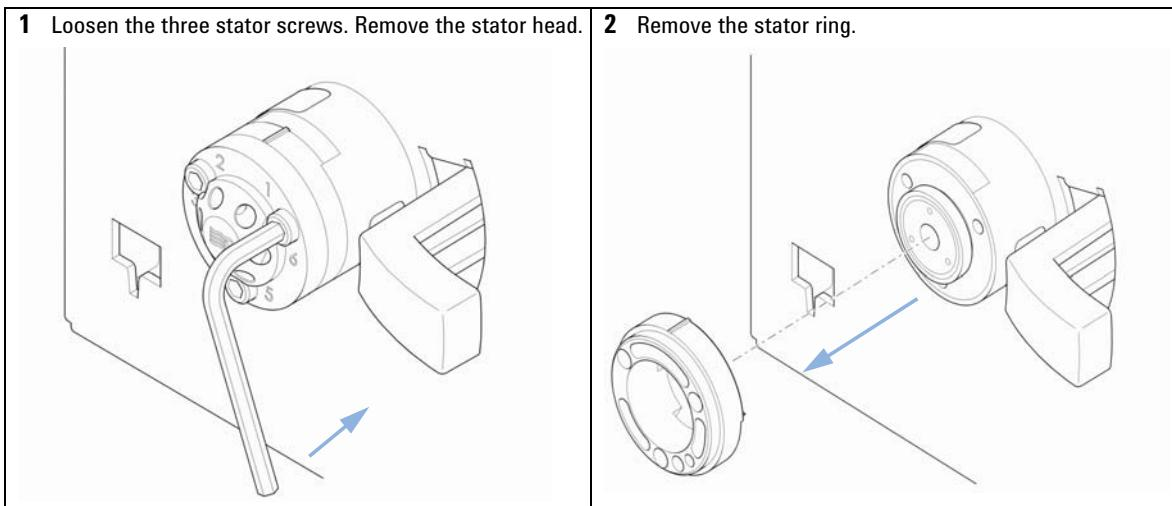
Manual Injector

Exchanging the Injection Valve Seal

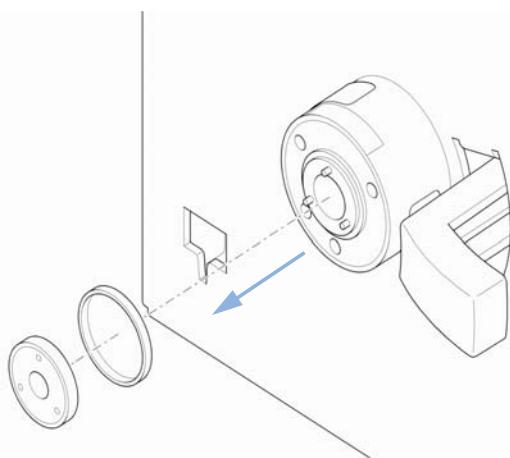
When Poor injection-volume reproducibility
Leaking injection valve

Tools required **Description**
Hex key, 9/64 inch
(supplied in the tool kit)

Parts required **p/n** **Description**
5068-0082 Rotor seal, PEEK

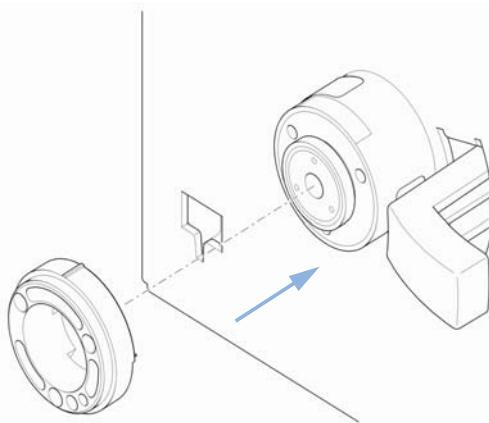


3 Remove the rotor seal.

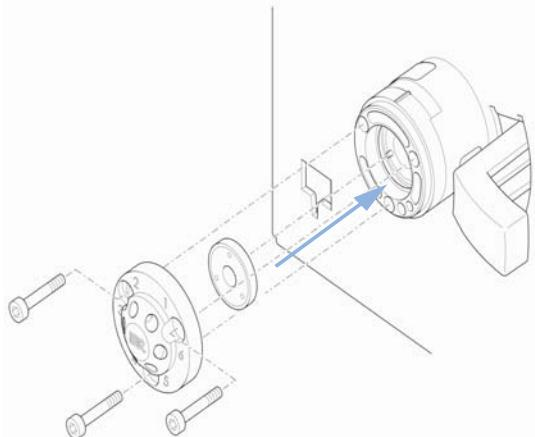


4 Install the new rotor seal.

5 Install the stator ring. Ensure the pin in the stator ring is aligned with the hole in the valve body.



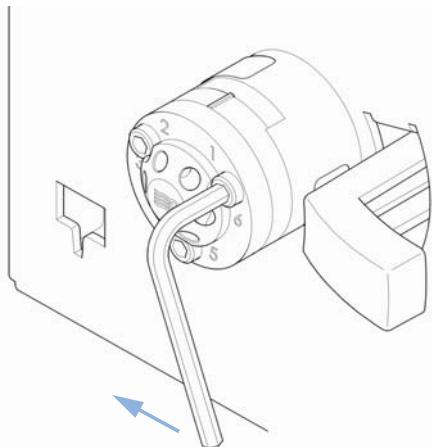
6 Install the stator head onto the valve.



11 Maintenance and Repair

Manual Injector

- 7 Secure stator head in place with the stator screws.
Tighten each screw alternately $\frac{1}{4}$ -turn until the stator head is secure.



Autosampler

Introduction

The autosampler is designed for easy repair. The most frequent repairs can be done from the front of the instrument with the instrument in place in the Agilent 1220 Infinity II LC. These repairs are described in following sections.

Overview of procedures

Procedure	Typical frequency	Time required
Exchanging the needle assembly	When needle shows indication of damage or blockage	15 min
Exchanging the seat assembly	When the seat shows indication of damage or blockage	10 min
Exchanging the rotor seal	After approximately 30000to 40000injections, or when the valve performance shows indication of leakage or wear	30 min
Exchanging the metering seal	When autosampler reproducibility indicates seal wear	30 min
Exchanging the gripper arm	When the gripper arm is defective	10 min

11 Maintenance and Repair

Autosampler

WARNING

The power supplies still use some power, even if the power switch on the front panel is turned off.

Repair work at the autosampler can lead to personal injuries, e.g. shock hazard, when the autosampler cover is opened and the instrument is connected to power.

- Make sure that it is always possible to access the power plug.
 - Remove the power cable from the instrument before opening the cover.
 - Do not connect the power cable to the Instrument while the covers are removed.
-

Exchanging Internal Parts

Some repairs may require exchange of defective internal parts. Exchange of these parts requires removing the autosampler unit from the Agilent 1220 Infinity II LC; these repairs have to be done by trained service personal only.

Safety flap, flex board

It is strongly recommended that the exchange of the safety flap, and flex board is done by Agilent-trained service personnel.

Transport assembly parts

The adjustment of the motors, and the tension on the drive belts are important for correct operation of the transport assembly. It is strongly recommended that exchange of drive belts, and the gripper assembly is done by Agilent-trained service personnel. There are no other field-replaceable parts in the transport assembly. If any other component is defective (flex board, spindles, plastic parts) the complete unit must be exchanged.

Cleaning the autosampler

WARNING

Electrical shock hazard

Liquid dripping into the autosampler could cause shock hazard and damage to the autosampler.

- Drain all solvent lines before opening any fittings.
-

The autosampler covers should be kept clean. Clean with a soft cloth slightly dampened with water or a solution of water and a mild detergent. Do not use an excessively damp cloth that liquid can drip into the autosampler.

Maintenance Functions

Certain maintenance procedures require the needle arm, metering device, and gripper assembly to be moved to specific positions to enable easy access to components. The maintenance functions move these assemblies into the appropriate maintenance position.

11 Maintenance and Repair

Autosampler

Exchanging the Needle Assembly

When When the needle is visibly damaged
When the needle is blocked

Tools required

	Description
	Wrench, 1/4 inch (supplied in HPLC Tool-Kit)
	Hexagonal key, 2.5 mm
	Pair of pliers

Parts required

#	p/n	Description
1	G1313-87201	Needle assembly

Preparations Select **Change Needle** in the **Tools** function in the Instrument Utilities or Lab Advisor Software and select **Start**.
When the needle is positioned approximately 15 mm above the needle seat, remove the upper front cover.

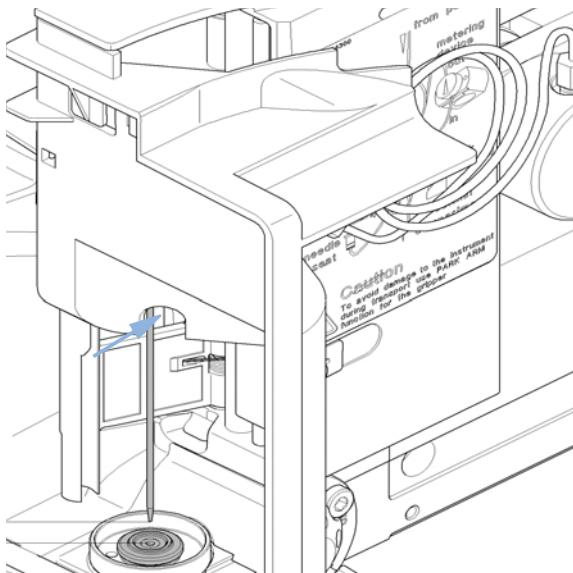
WARNING

Personal injury

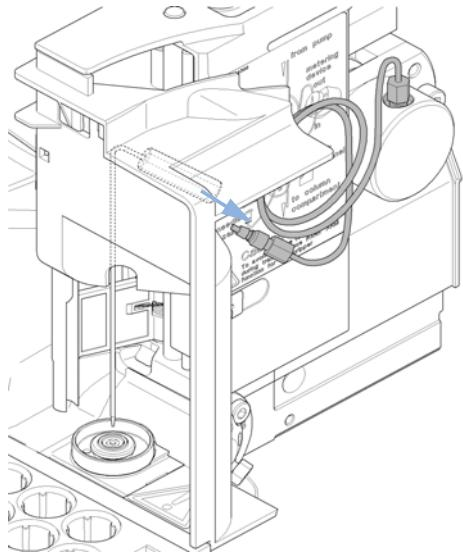
To avoid personal injury, keep fingers away from the needle area during autosampler operation.

- Do not bend the safety flap away from its position, or attempt to remove the safety cover.
- Do not attempt to insert or remove a vial from the gripper when the gripper is positioned below the needle.

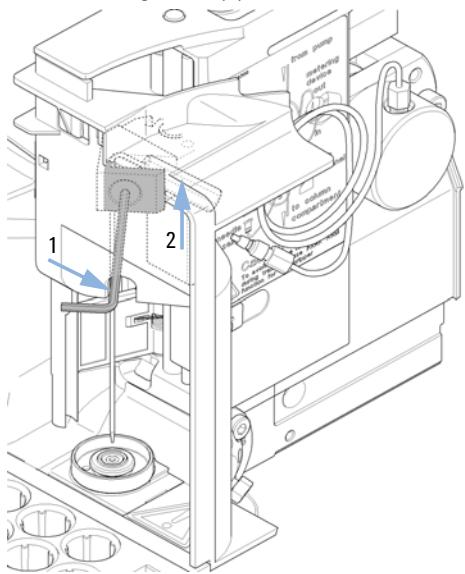
- 1** Select **Needle Down** until the needle screw is aligned with the hole in the safety cover.



- 2** Remove the sample-loop fitting from the needle fitting.



- 3** Loosen the fixing screw (1), and lift out the needle (2).



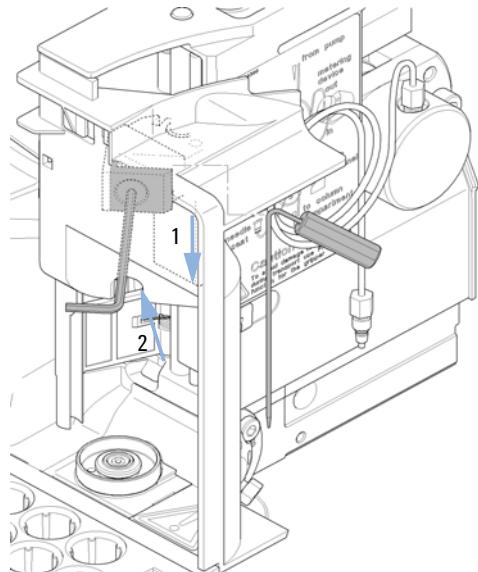
- 4** Select **Needle Down** to move the needle arm to its lowest position.

The needle arm must be in its lowest position before installing the new needle, otherwise leaks at the needle seat will occur due to incorrect needle installation.

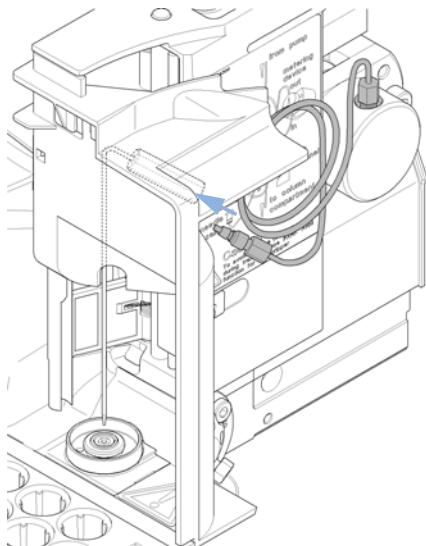
11 Maintenance and Repair

Autosampler

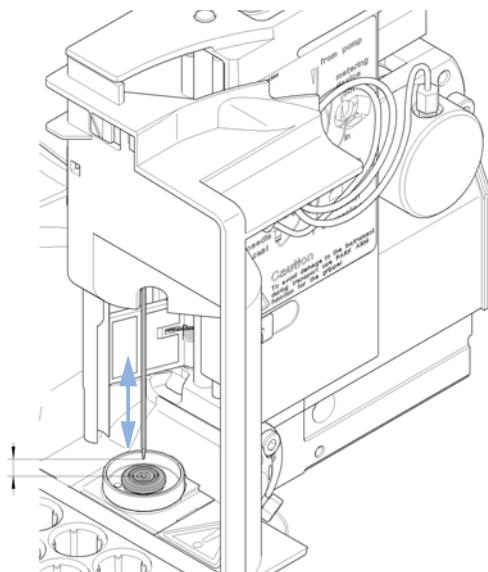
- 5** Insert the new needle (1). Align the needle in the seat, then tighten the screw firmly (2).



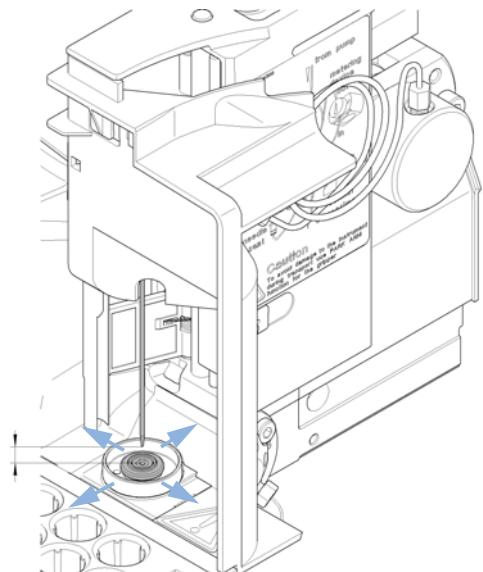
- 6** Reconnect the sample-loop fitting to the needle fitting.



- 7** Use **Needle Up** to lift the needle to a position approximately 2 mm above the seat.



- 8** Ensure the needle is aligned with the seat.



Next Steps:

- 9** On completion of this procedure: Install the front cover.
- 10** Select **End** in the **Tools** function **Change Needle**.

11 Maintenance and Repair

Autosampler

Exchanging the Needle Seat Assembly

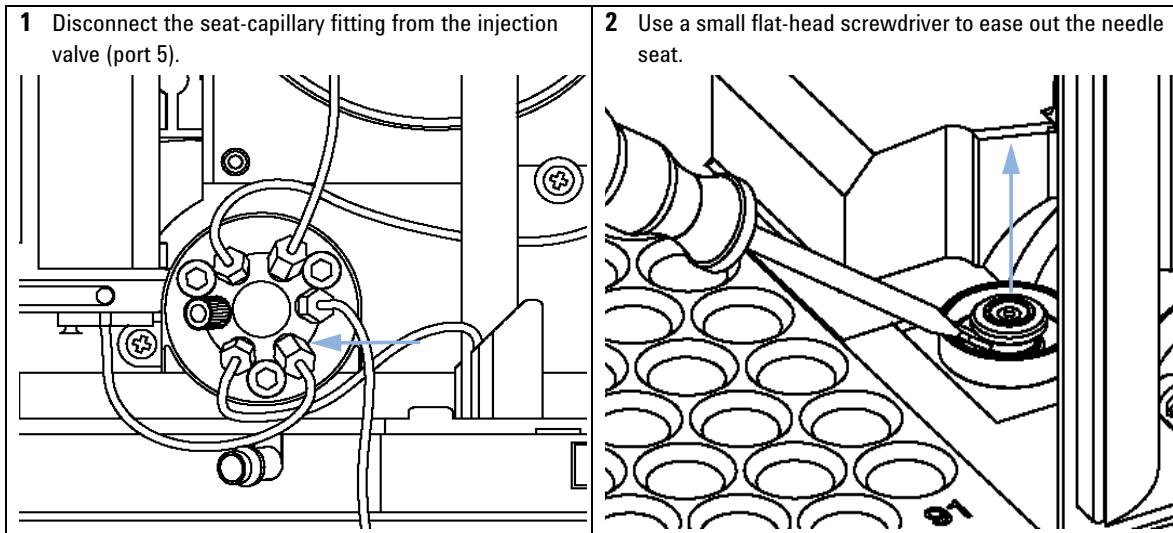
When When the seat is visibly damaged
When the seat capillary is blocked

Tools required **Description**
Wrench, 1/4 inch (supplied in HPLC Tool-Kit)
Screwdriver, flat-head

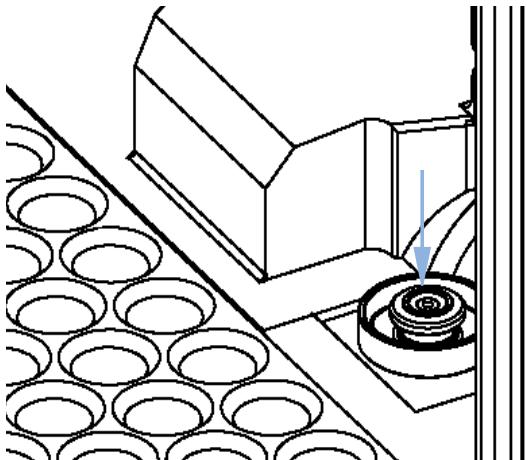
Parts required	#	p/n	Description
	1	G1313-87101	Needle-seat assy (0.17 mm i.d. 2.3 µL)

Preparations

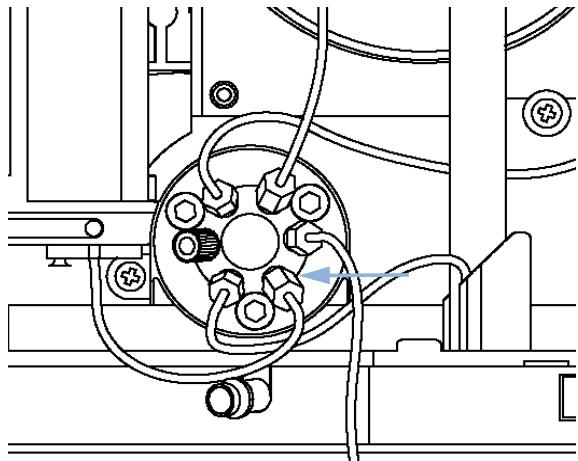
- Select **Start** in the **Tools** function in LMD Software **Change Needle**
- Remove the upper front cover.
- Use the **Needle Up** command in the **Change Needle** function to lift the needle an additional 1 cm.



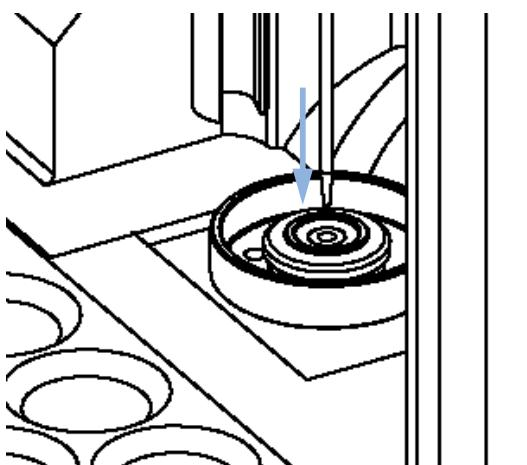
- 3** Insert the new needle-seat assembly. Press the seat firmly into position.



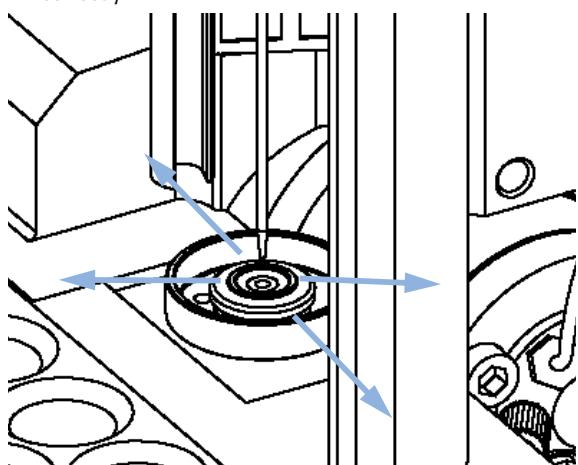
- 4** Connect the seat-capillary fitting to port 5 of the injection valve.



- 5** Use **Down** to position the needle approximately 2 mm above the seat.



- 6** Ensure the needle is aligned with the seat. If required, bend the needle slightly until the needle is aligned correctly.



Next Steps:

- 7** On completion of this procedure: Install the front cover.
8 Select **End** in the **Tools** function **Change Needle**.

11 Maintenance and Repair

Autosampler

Exchanging the Rotor Seal

When Poor injection-volume reproducibility
Leaking injection valve

Tools required **Description**
Wrench 1/4 inch
Hex key, 9/16 inch (supplied in the tool kit)

Parts required # p/n Description
1 0101-1416 Rotor seal (PEEK)

Preparations

- Remove upper front cover.
- Remove the leak tubing (if necessary).

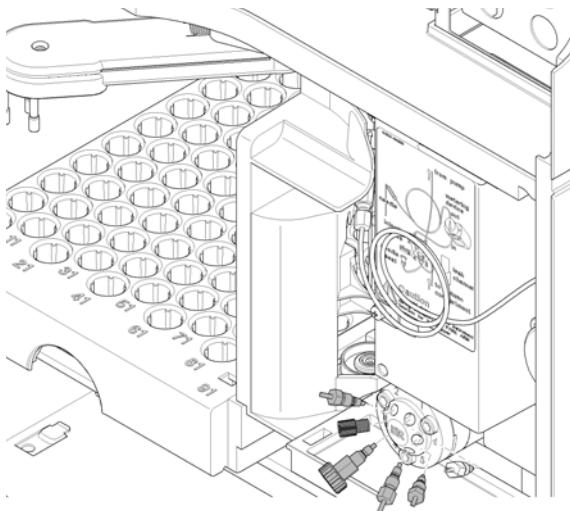
CAUTION

Removing the stator head

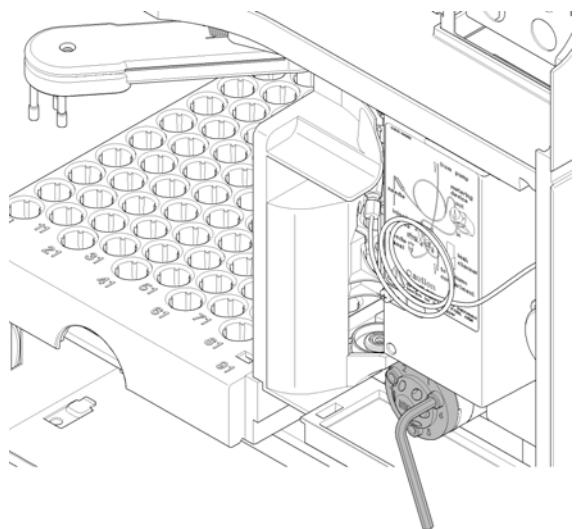
The stator face is held in place by the stator head. When you remove the stator head, the stator face can fall out of the valve.

→ Carefully handle the valve to prevent damage to the stator face

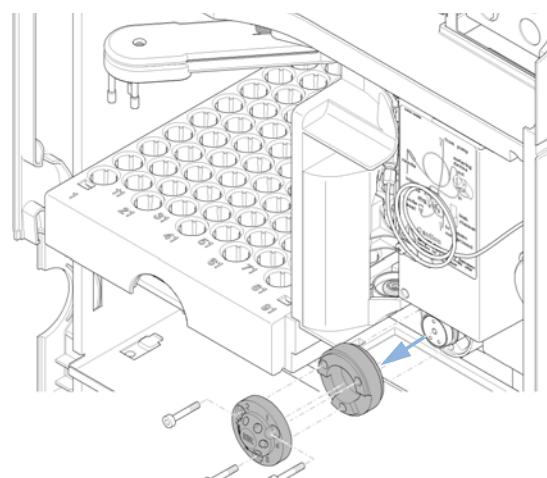
- 1** Remove all capillary fittings from the injection-valve ports.



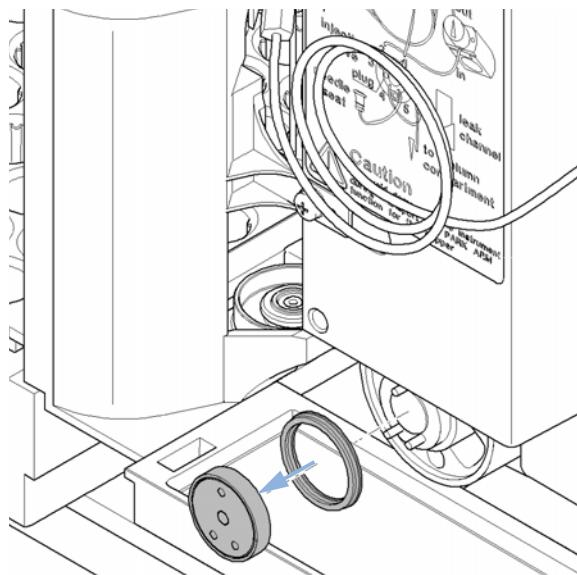
- 2** Loosen each fixing bolt two turns at a time. Remove the bolts from the head.



- 3** Remove the stator head and stator ring.



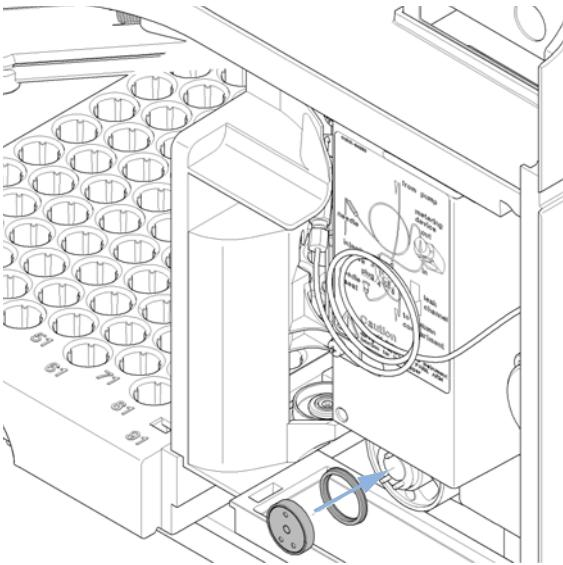
- 4** Remove the rotor seal and isolation seal.



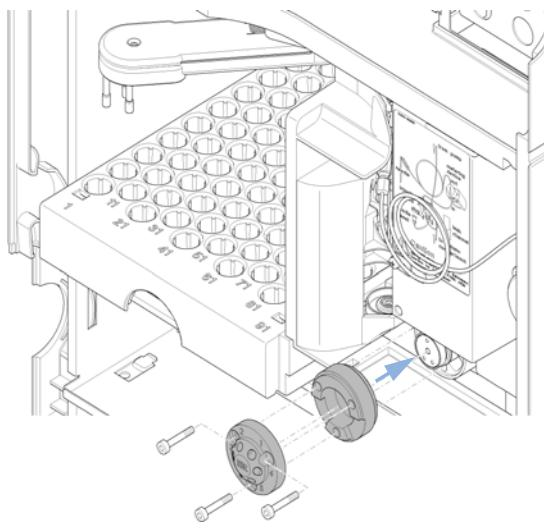
11 Maintenance and Repair

Autosampler

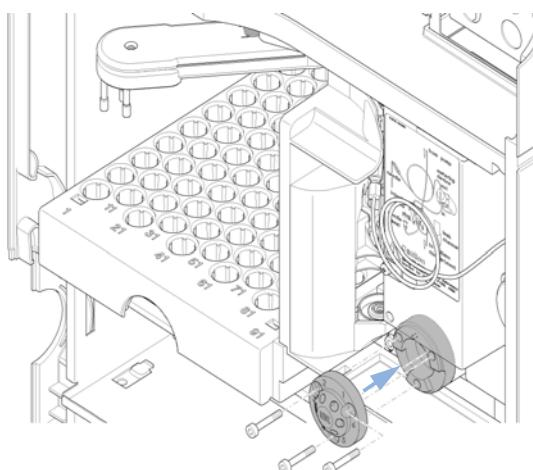
- 5** Install the new rotor seal and isolation seal. Ensure the metal spring inside the isolation seal faces towards the valve body.



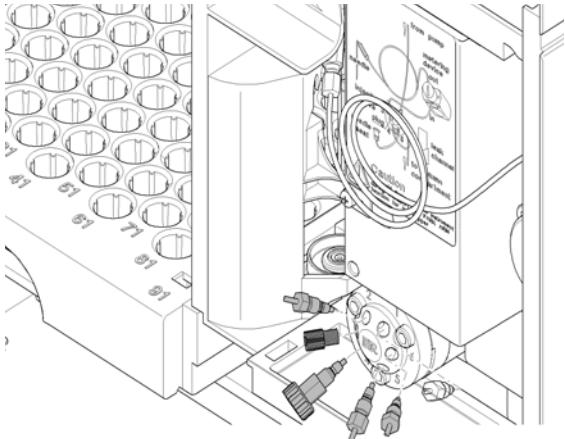
- 6** Install the stator ring with the short of the two pins facing towards you at the 12 o'clock position. Ensure the ring sits flat on the valve body.



- 7** Install stator head. Tighten the bolts alternately two turns at a time until the stator head is secure.



- 8** Reconnect the pump capillaries to the valve ports.



Next Steps:

- 9** Slide the waste tube into the waste holder in the leak tray.
- 10** On completion of this procedure: Install the front cover.

11 Maintenance and Repair

Autosampler

Exchanging the Metering Seal

When Poor injection-volume reproducibility
Leaking metering device

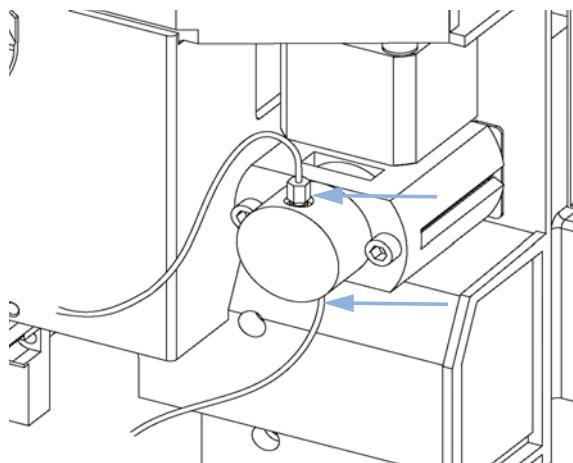
Tools required	p/n	Description
		Wrench, 1/4 inch (supplied in HPLC Tool-Kit)
		Hexagonal key, 4 mm (supplied in HPLC Tool-Kit)
	8710-2411	Hexagonal key, 3 mm (supplied in HPLC Tool-Kit)

Parts required	#	p/n	Description
	1	5063-6589	Metering seal (pack of 2) for 100 µL analytical head
	1	5063-6586	Sapphire piston

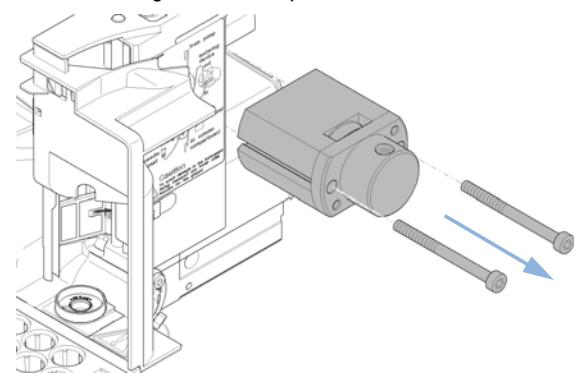
Preparations

- Select **Start** in the **Tools** function in the LMD software **Change piston**.
- Remove the upper front cover.

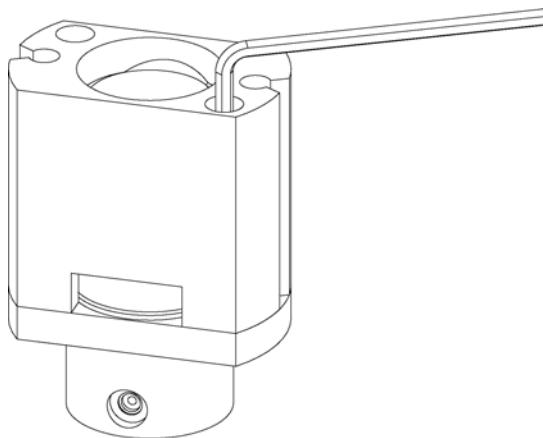
- 1** Remove the two capillaries from the metering-head assembly.



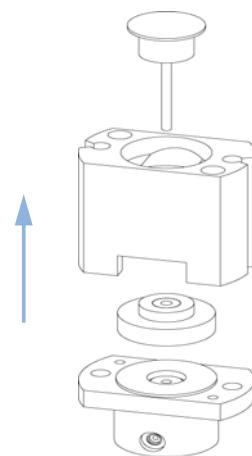
- 2** Remove the two fixing bolts, and pull the head assembly away from the sampler. Notice that the closed side of the metering head faces upwards.



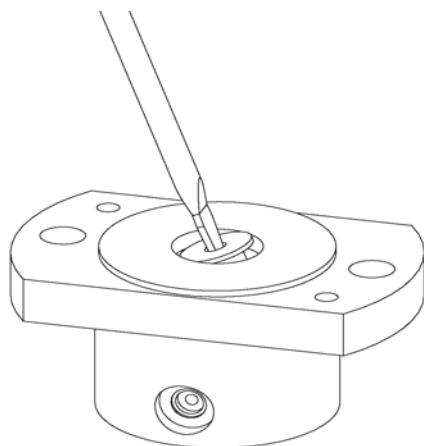
- 3** Remove the two fixing bolts from the base of the metering head assembly.



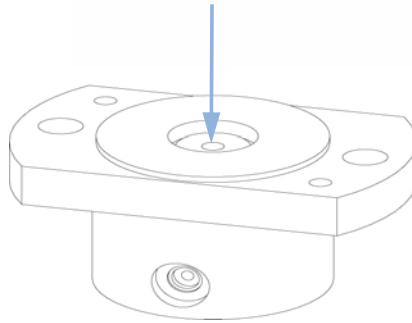
- 4** Disassemble the metering head assembly.



- 5** Use a small screwdriver to carefully remove the seal. Clean the chamber with lint-free cloth. Ensure all particular matter is removed.



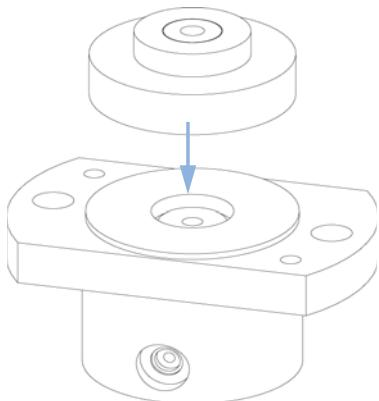
- 6** Install the new seal. Press the seal firmly into position.



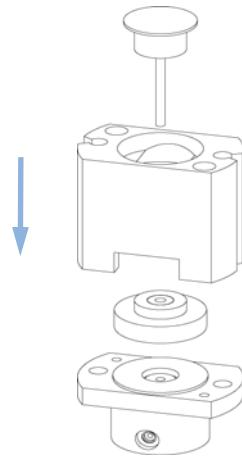
11 Maintenance and Repair

Autosampler

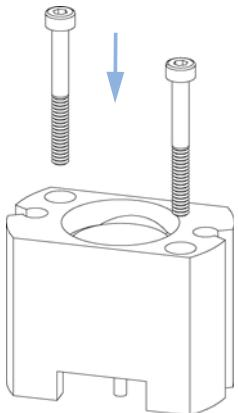
7 Place the piston guide on top of the seal.



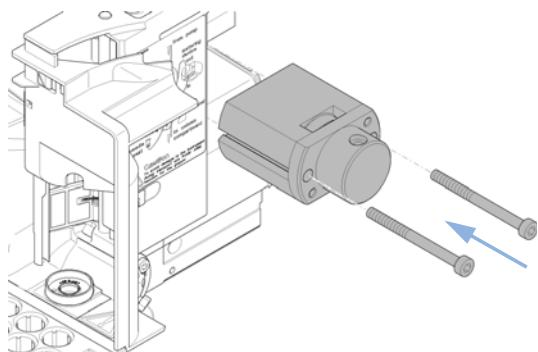
8 Reassemble the metering head assembly. Carefully insert the plunger into the base. The closed side of the metering head must be on the same side as the lower one of the two capillary drillings.

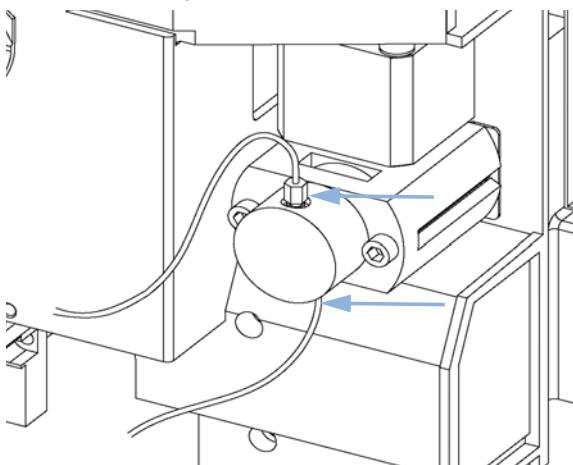


9 Install the fixing bolts. Tighten the bolts securely.



10 Install the metering head assembly in the autosampler. Ensure the large hole in the metering head is facing downwards.



11 Reinstall the capillaries.**Next Steps:**

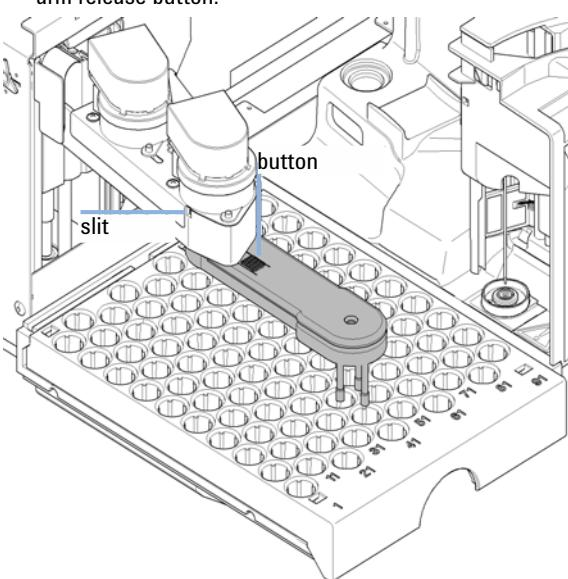
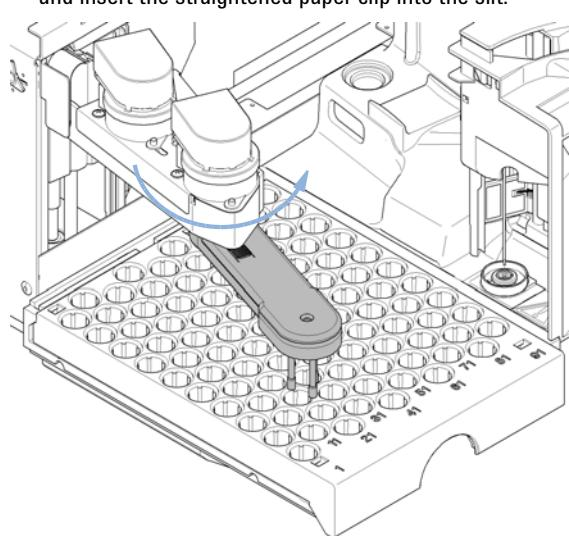
12 On completion of this procedure: Install the front cover.

13 Select **End** in the **Tools** function in the LMD software
Change piston.

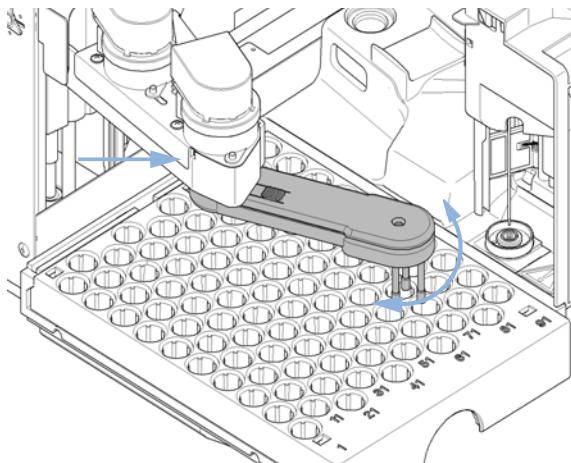
11 Maintenance and Repair

Autosampler

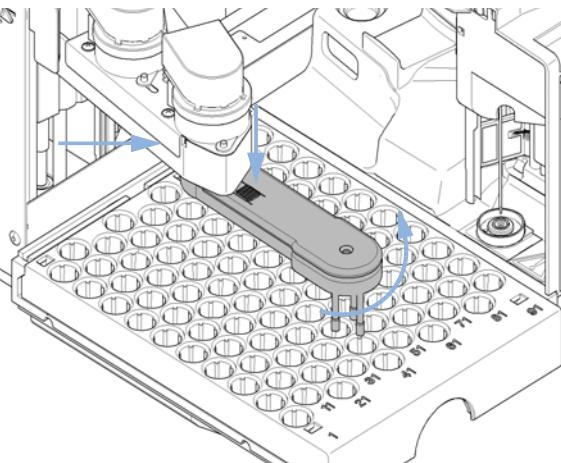
Exchanging the Gripper Arm

When	Defective gripper arm		
Tools required	Description Straightened paper clip.		
Parts required	#	p/n	Description
	1	G1313-60010	Gripper assembly
Preparations	<ul style="list-style-type: none">Select Start in the Tools function in the LMD software Change Gripper.Turn off the power to the instrument.Remove the upper front cover.		
1	Identify the slit below the gripper motor and the gripper arm release button.		
			
2	Rotate the arm approximately 2.5 cm (1 inch) to the left and insert the straightened paper clip into the slit.		
			

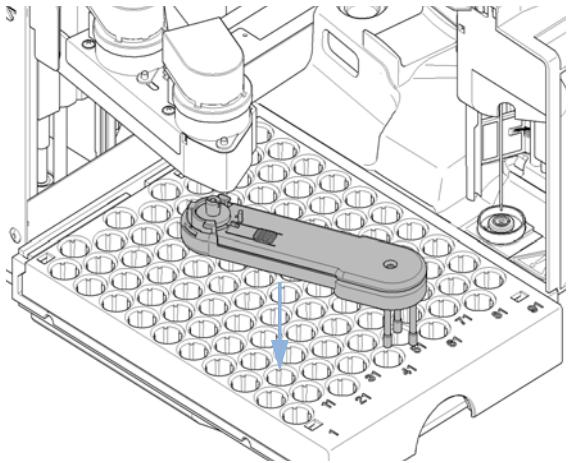
- 3** Rotate the gripper arm slowly from left to right and apply a gentle pressure to the paper clip. The clip will engage on an internal catch and the rotation of the arm will be blocked.



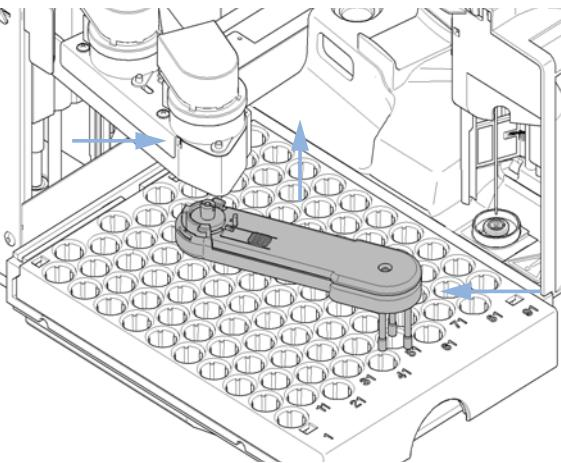
- 4** Hold the paper clip in place, press the gripper release button and rotate the gripper arm to the right.



- 5** The gripper arm will come off.



- 6** Replace the gripper arm by holding the paper clip in place, pushing the gripper arm into the holder and rotating the gripper arm to the left.



Next Steps:

- 7** On completion of this procedure: Install the front cover.
8 Turn the power to the Instrument ON.

11 Maintenance and Repair

Variable Wavelength Detector (VWD)

Variable Wavelength Detector (VWD)

Introduction

This section describes simple maintenance and repair procedures for the detector that can be carried out without opening the main cover.

Table 46 Detector maintenance and repair

Procedure	Typical Frequency	Notes
Exchanging the deuterium lamp	If noise and/or drift exceeds your application limits or lamp does not ignite.	A VWD test should be performed after replacement.
Exchanging the flow cell	If the application requires a different flow cell type.	A VWD test should be performed after replacement.
Repairing the flow cell	If leaking or if intensity drops due to contaminated flow cell windows.	A pressure tightness test should be done after repair.
Drying the leak sensor	If leak has occurred.	Check for leaks.

Exchanging the Deuterium Lamp

When If noise or drift exceeds application limits or lamp does not ignite.

Tools required **Description**
Screwdriver, Pozidriv #1 PT3

Parts required **#** **p/n** **Description**
1 G1314-60100 Deuterium lamp

Preparations Turn the lamp OFF.

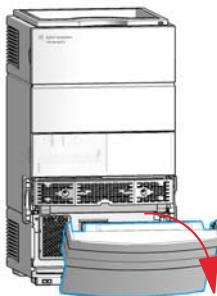
WARNING

Injury by touching hot lamp

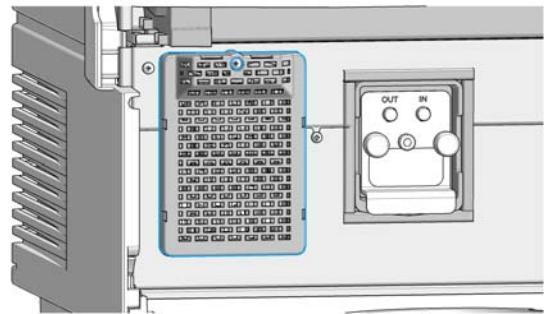
If the detector has been in use, the lamp may be hot.

→ If so, wait for lamp to cool down.

- 1** Remove the lower front cover.



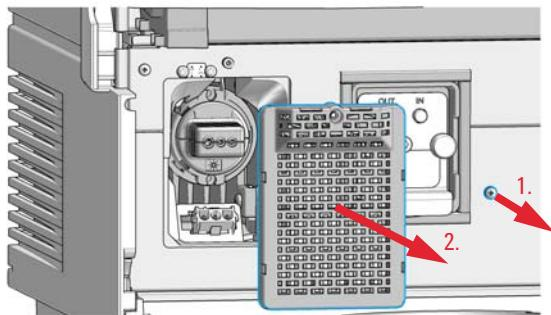
- 2** Locate the heater fan cover.



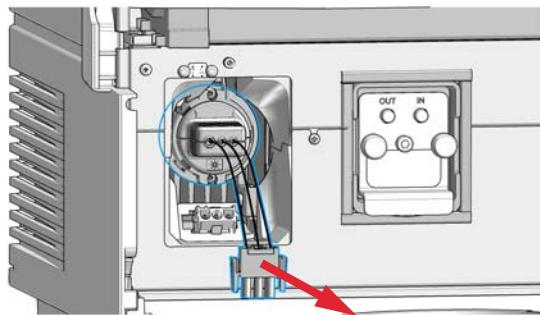
11 Maintenance and Repair

Variable Wavelength Detector (VWD)

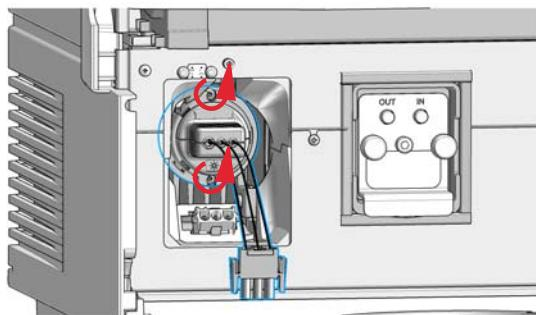
- 3 Unscrew the fan cover (1.) and remove it (2.).



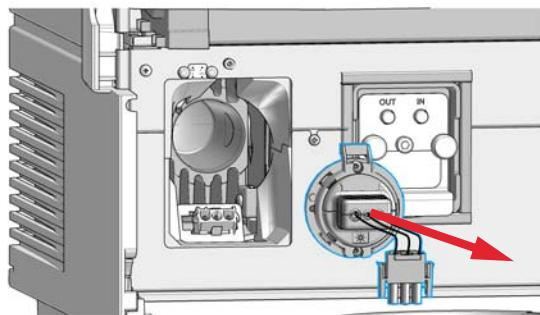
- 4 Unplug the lamp connector.



- 5 Unscrew the two lamp screws (Pozidriv).



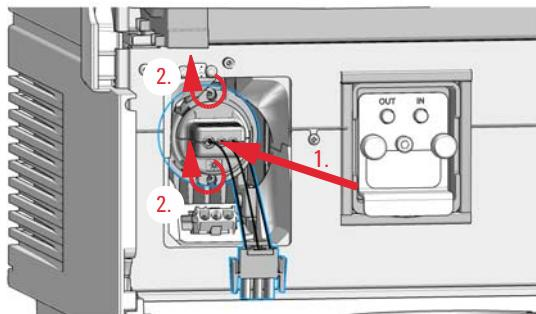
- 6 Remove the lamp and place it on a clean place.



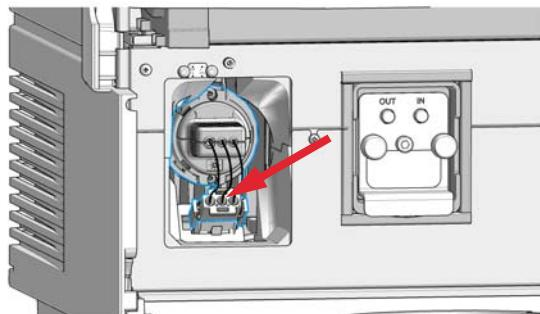
NOTE

Do not touch the glass bulb with your fingers. It may reduce the light output.

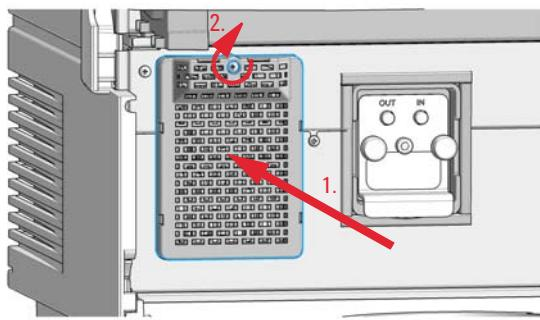
- 7 Insert the lamp (RFID tag on top) (1.) and fix the screws (2.).



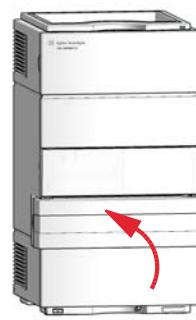
- 8 Reconnect the connector.



- 9** Replace the fan cover (1.) and fix its screw (2.).



- 10** Close the lower front cover.



Next Steps:

- 11** Reset the lamp counter as described in the Utilities software documentation.
- 12** Turn the lamp ON.
- 13** Give the lamp more than 10 min to warm-up.
- 14** Perform Wavelength Calibration to check the correct positioning of the lamp.

11 Maintenance and Repair

Variable Wavelength Detector (VWD)

Replace the Flow Cell / Cuvette Holder

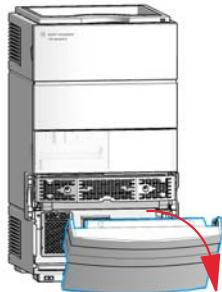
When If application needs a different type of flow cell or the flow cell needs repair.

Tools required **Description**
Wrench, 1/4 inch
for capillary connections

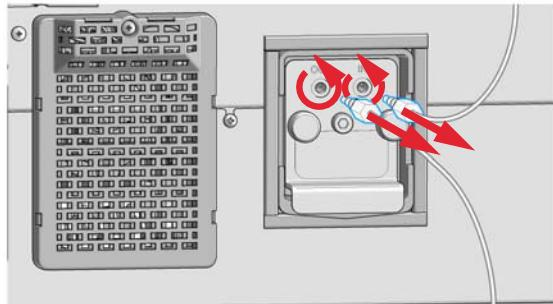
Parts required # **Description**
1 Flow cell

Preparations Turn the lamp OFF.

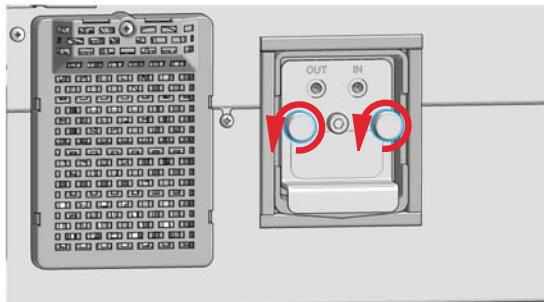
- 1** Remove the lower front cover.



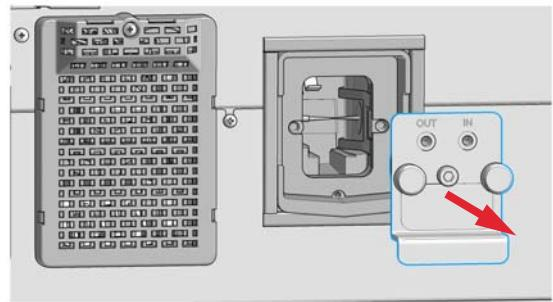
- 2** Disconnect the inlet capillary and outlet tubing from the flow cell.



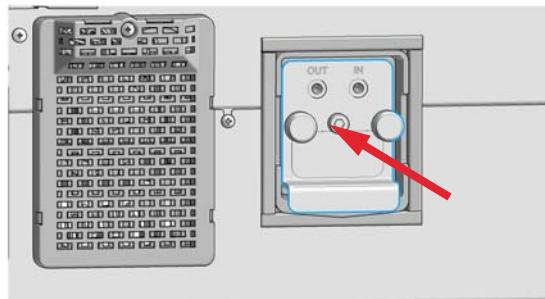
- 3** Unscrew the two thumb screws.



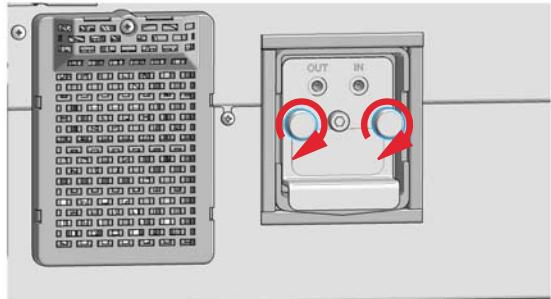
- 4** Pull the flow cell out of its location.



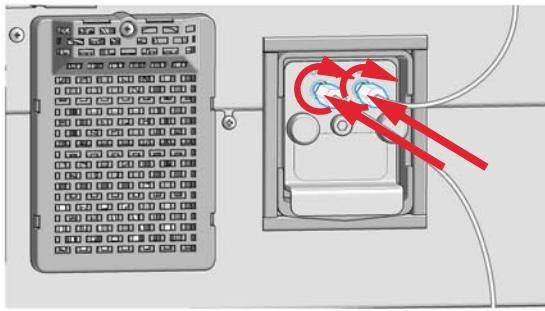
- 5 Insert the flow cell into its location.



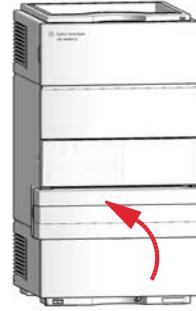
- 6 Fix the two thumb screws parallel.



- 7 Reconnect the inlet capillary.



- 8 Close the lower front cover.



Next Steps:

- 9 Configure the flow cell.

- 10 Perform a **Wavelength Verification-Calibration** to check the correct position of the flow cell/cuvette holder.

11 Maintenance and Repair

Variable Wavelength Detector (VWD)

Repairing the Flow Cell

Parts required	p/n	Description
	G1314-60086	Standard flow cell, 10 mm, 14 µL, 40 bar
	G1314-65061	Cell Repair Kit, includes 2x Gasket #1, 2x Gasket #2, 2x Window Quartz

CAUTION

Window surfaces scratched by tweezers

Window surfaces can easily be scratched by using tweezers to remove the windows.

→ Do not use tweezers to remove the windows

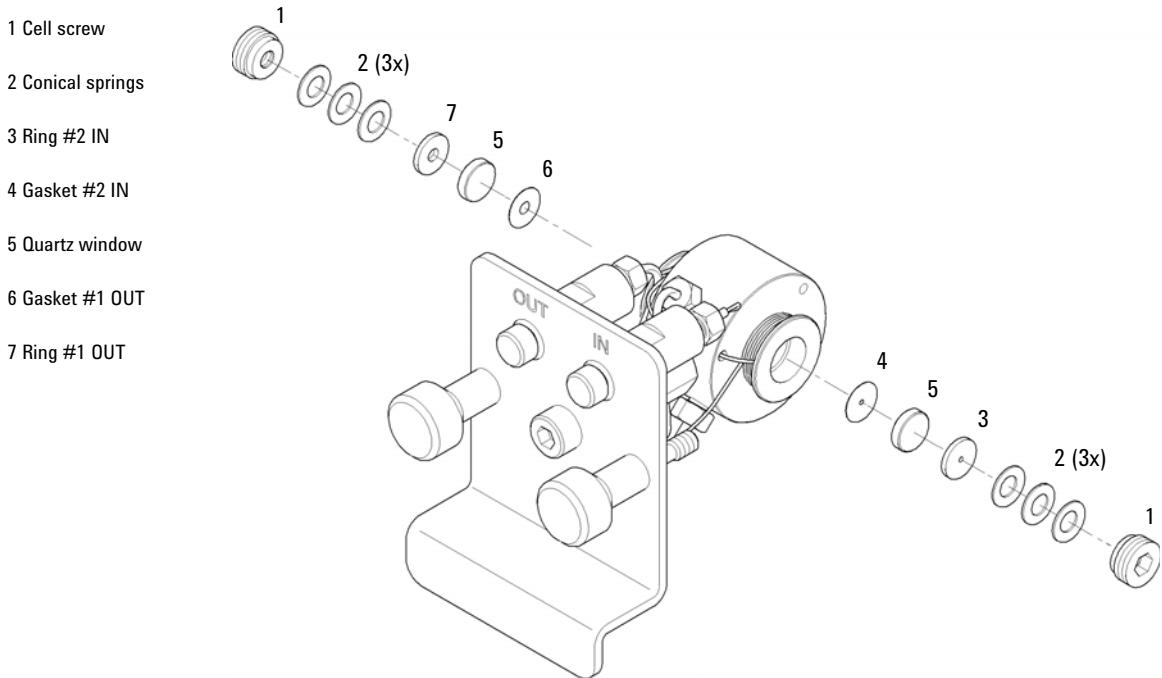


Figure 65 Standard Flow Cell 10 mm / 14 μ L

- 1 Press the release buttons and remove the lower front cover to gain access to the flow cell area
- 2 Disconnect the inlet and outlet capillaries.
- 3 Unscrew the thumb screws in parallel and remove the flow cell.
- 4 Disassembling the flow cell.
 - a Unscrew the cell screw using a 4 mm hexagonal wrench.
 - b Remove the SST rings using a pair of tweezers.
 - c Use adhesive tape to remove the PEEK ring, the window and the gasket.
 - d Repeat step 1 on page 297 through step 3 on page 297 for the other window. Keep the parts separate, otherwise they could get mixed up.

11 Maintenance and Repair

Variable Wavelength Detector (VWD)

5 Cleaning the flow cell parts

- a Pour isopropanol into the cell hole and wipe clean with a piece of lint-free cloth.
- b Clean the windows with ethanol or methanol. Dry with a piece of lint-free cloth.

6 Reassembling the flow cell

- a Hold the flow cell cassette horizontal and place the gasket in position. Ensure that both cell holes can be seen through the holes of gasket.

NOTE

Always use new gaskets.

- b Place the the window on gasket.

- c Place the PEEK ring on the window.

- d Insert the conical springs. Make sure that the conical springs point towards the window, otherwise the window might break when the cell screw is tightened.

- e Screw the cell screw into the flow cell and tighten it.

7 Repeat the procedure for the other side of the cell.

8 Reconnect the inlet and outlet capillaries.

9 Test the flow cell for leaks. If there are no leaks, insert the flow cell into the detector.

10 Perform **Wavelength Calibration** to check the correct positioning of the flow cell (“[Wavelength Verification/Calibration](#)” on page 168).

11 Replace the lower front cover.

Using the Cuvette Holder

This cuvette holder can be placed instead of a flow cell in the variable wavelength detector. Standard cuvettes with standards in it, for example, National Institute of Standards & Technology (NIST) holmium oxide solution standard, can be fixed in it.

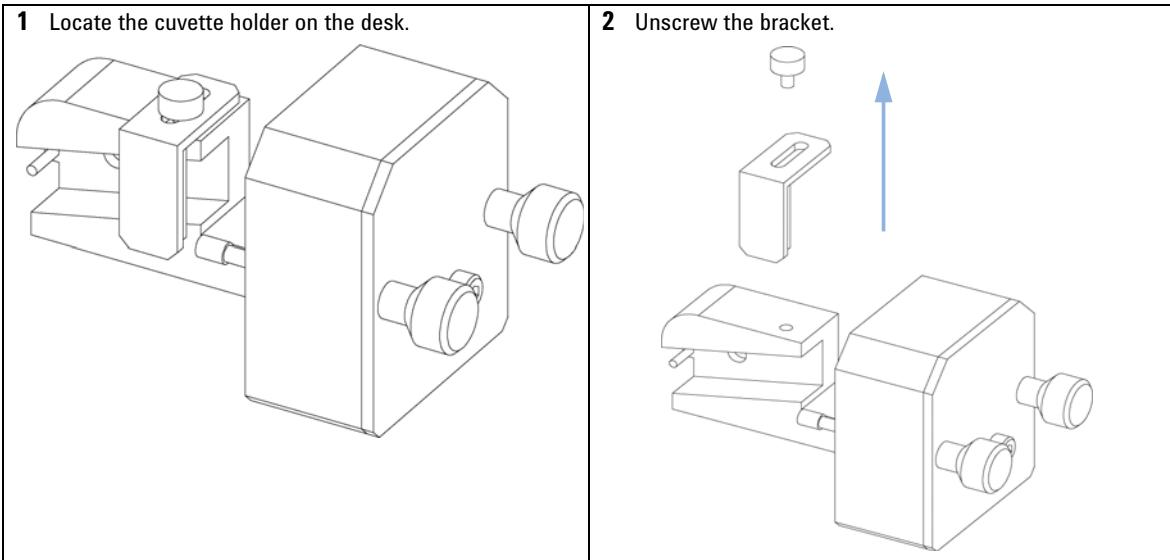
This can be used for wavelength verifications.

When If your own standard should be used to checkout the instrument.

Tools required **Description**
None

Parts required **p/n** **Description**
G1314-60200 Cuvette Holder
Cuvette with the "standard", e.g. NIST certified holmium oxide sample

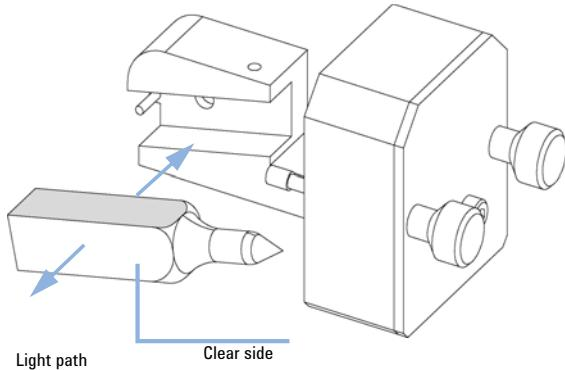
Preparations Remove the normal flow cell.
Have cuvette with standard available.



11 Maintenance and Repair

Variable Wavelength Detector (VWD)

- 3** Insert the cuvette with the sample into the holder. The clear side of the cuvette must be visible.



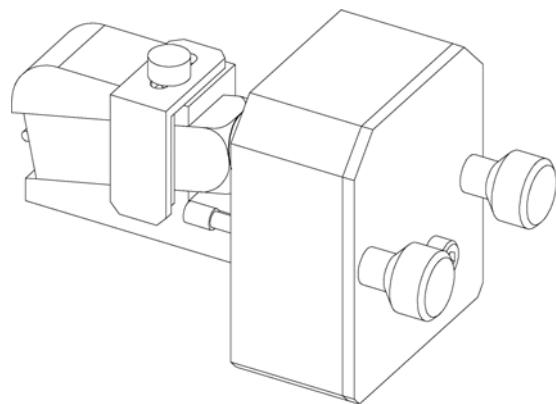
- 4** Reset the lamp counter as described in the user interface documentation.

- 5** Turn the lamp ON.

- 6** Give the lamp more than 10 min to warm-up.

- 7** Perform **Wavelength Verification/Calibration** to check the correct positioning of the lamp.

- 8** Replace the bracket and fix the cuvette.



Next Steps:

- 9** Install the cuvette holder in the instrument.

- 10** Perform your verification.

Correcting Leaks

When If a leakage has occurred in the flow cell area or at the capillary connections

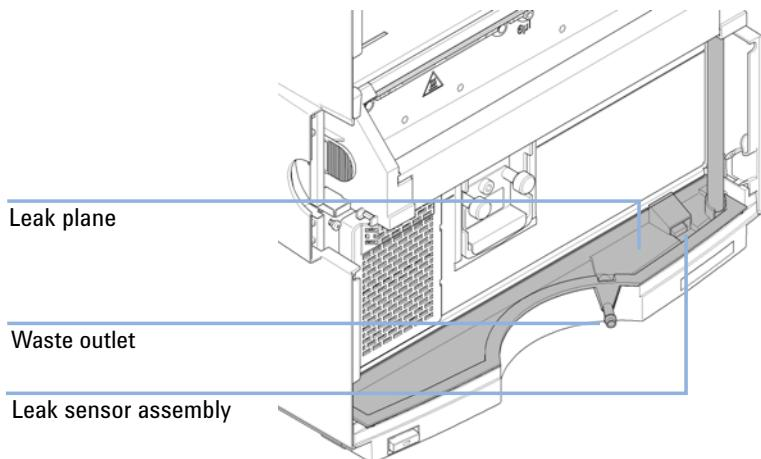
Tools required

Description
Tissue
Wrench, 1/4 inch
for capillary connections

Parts required

Description
None

- 1 Remove the lower front cover.
- 2 Use tissue to dry the leak sensor area.
- 3 Observe the capillary connections and the flow cell area for leaks and correct, if required.
- 4 Replace the front cover.



Diode Array Detector (DAD)

Overview of Maintenance

The following pages describe maintenance (simple repairs) of the detector that can be carried out without opening the main cover.

Table 47 Overview of Maintenance

Procedure	Typical Frequency	Notes
Cleaning of module	If required.	
Deuterium lamp or tungsten lamp exchange	If noise and/or drift exceeds your application limits or lamp does not ignite.	An intensity test should be performed after replacement.
Flow cell exchange	If application requires a different flow cell type.	A holmium or wavelength calibration test should be performed after replacement.
Flow cell parts Cleaning or exchange	If leaking or if intensity drops due to contaminated flow cell windows.	A pressure tightness test should be done after repair.
Holmium oxide filter Cleaning or exchange	If contaminated.	A holmium or wavelength calibration test should be performed after replacement.
Leak sensor drying	If leak has occurred.	Check for leaks.
Leak handling System replacement	If broken or corroded.	Check for leaks.

Cleaning the Module

The module case should be kept clean. Cleaning should be done with a soft cloth slightly dampened with water or a solution of water and mild detergent. Do not use an excessively damp cloth allowing liquid to drip into the module.

WARNING

Liquid dripping into the electronic compartment of your module can cause shock hazard and damage the module

- Do not use an excessively damp cloth during cleaning.
 - Drain all solvent lines before opening any connections in the flow path.
-

Exchanging a Lamp

When If noise or drift exceeds application limits or lamp does not ignite

Tools required **Description**
Screwdriver, Pozidriv #1 PT3

Parts required **#** **p/n** **Description**
1 2140-0820 Longlife Deuterium lamp "C" (with black cover and RFID tag)
OR 1 G1103-60001 Tungsten lamp

WARNING



Eye damage by detector light

Eye damage may result from directly viewing the light produced by the deuterium lamp used in this product.

→ Always turn the deuterium lamp off before removing it.

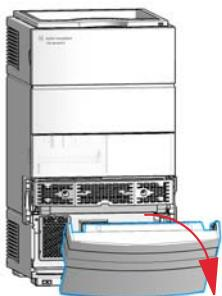
WARNING

Injury by touching hot lamp

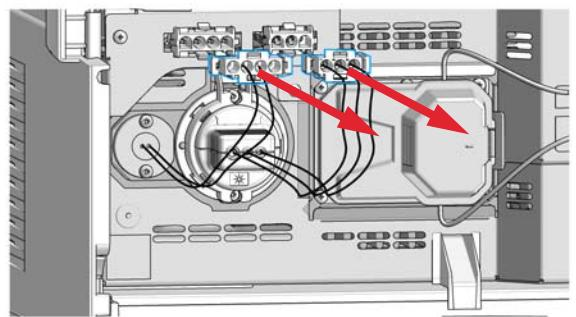
If the detector has been in use, the lamp may be hot.

→ If so, wait for lamp to cool down.

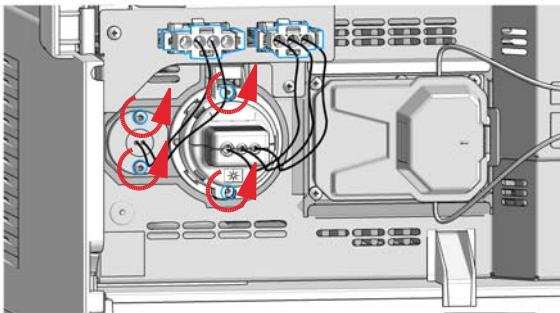
- 1** Remove the lower front cover.



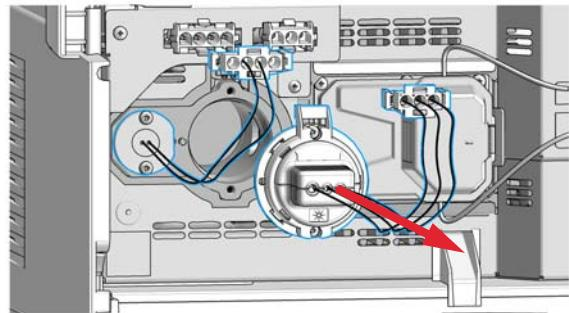
- 2** Unplug the lamp connector as required.



- 3** Unscrew (do not remove) the four lamp screws (Pozidriv) as required.



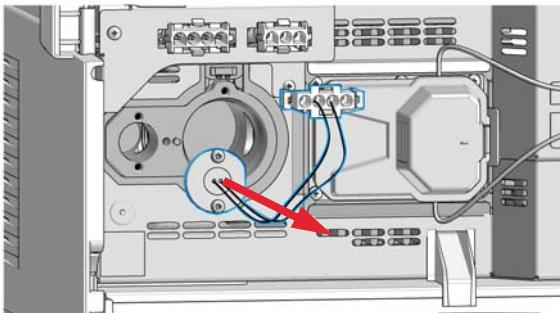
- 4** Remove the deuterium lamp and place it on a clean place.



NOTE

Do not touch the glass bulb with your fingers. It may reduce the light output.

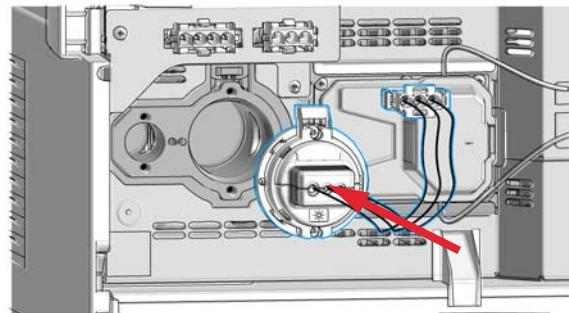
- 5** Remove the Vis-lamp and place it on a clean place.



NOTE

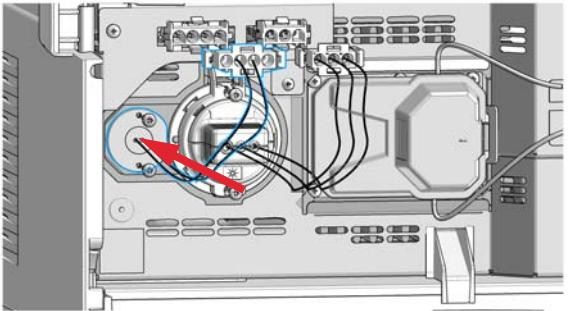
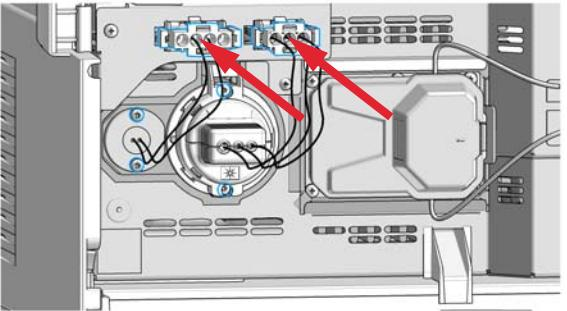
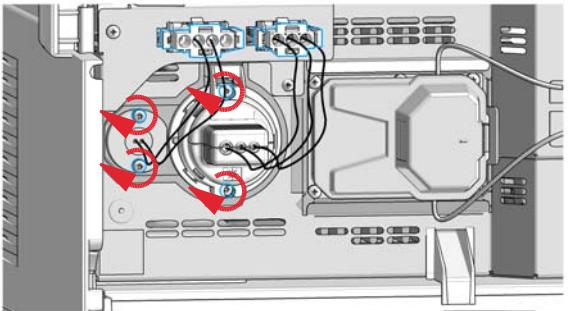
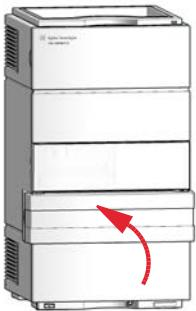
Do not touch the glass bulb with your fingers. It may reduce the light output.

- 6** Insert the deuterium lamp (RFID tag on the top side).



11 Maintenance and Repair

Diode Array Detector (DAD)

<p>7 Insert the Vis-lamp (flat side to the right).</p> 	<p>8 Reconnect the lamp connector as required.</p> 
<p>9 Fix the screws.</p> 	<p>10 Place the lamp cables in the lamp cover.</p>
<p>11 Close the lower front cover.</p> 	<p>Next Steps:</p> <p>12 Reset the lamp counter as described in the user interface documentation (lamps with I.D. tag cannot be reset).</p> <p>13 Turn the lamp on and give the lamp 10 minutes to warm up.</p> <p>14 Perform a “Wavelength Verification and Recalibration” on page 188 or a “Holmium Oxide Test” on page 180 to check the correct positioning of the UV-lamp.</p> <p>15 Perform an “Intensity Test” on page 177.</p>

Remove and Install a Flow Cell

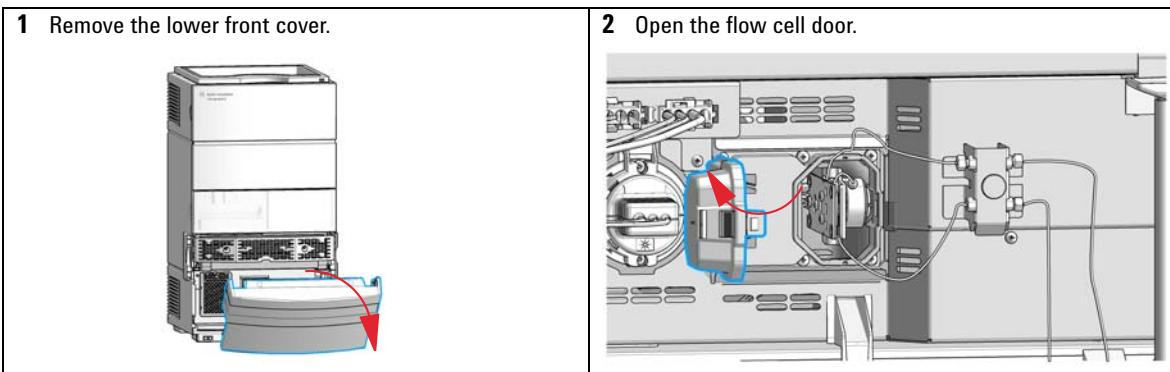
When If an application needs a different type of flow cell or the flow cell needs repair.

Tools required

Description
Wrench, 1/4 inch
for capillary connections

Parts required	p/n	Description
	G1315-60022	Standard flow cell, 10 mm, 13 µL, 120 bar (12 MPa)
	G1315-60025	Semi-micro flow cell, 6 mm, 5 µL, 120 bar (12 MPa)
	G1315-60024	Micro flow cell, 3 mm, 2 µL, 120 bar (12 MPa)
	G1315-60015	High pressure flow cell, 6 mm, 1.7 µL, 400 bar (40 MPa) Nano flow cell

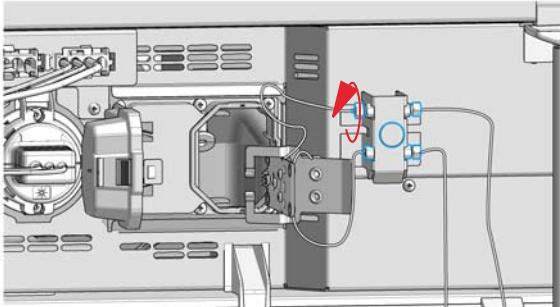
Preparations Turn the lamp(s) off.



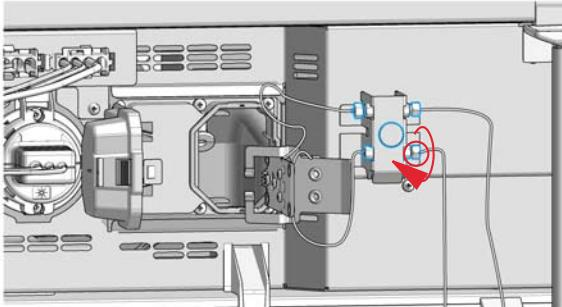
11 Maintenance and Repair

Diode Array Detector (DAD)

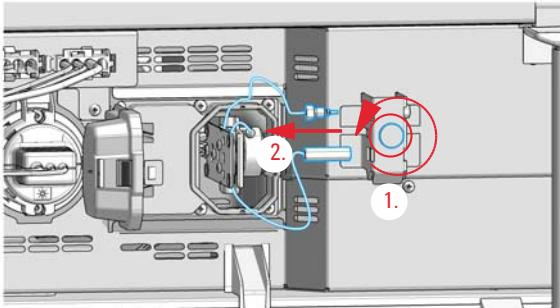
- 3** Disconnect the flow cell inlet capillary (top) from the union.



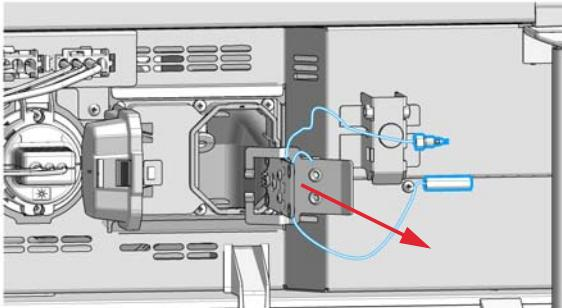
- 4** Disconnect the waste tubing (bottom) from the union.



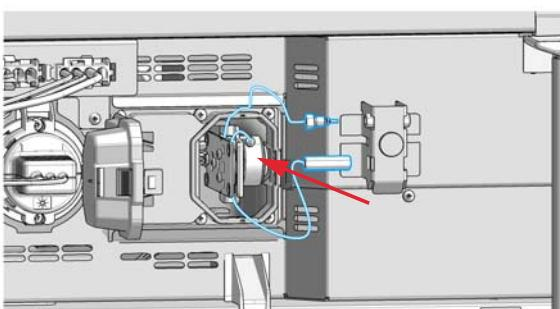
- 5** Loosen the thumb screw and remove the flow cell outlet capillary (bottom) with the union.



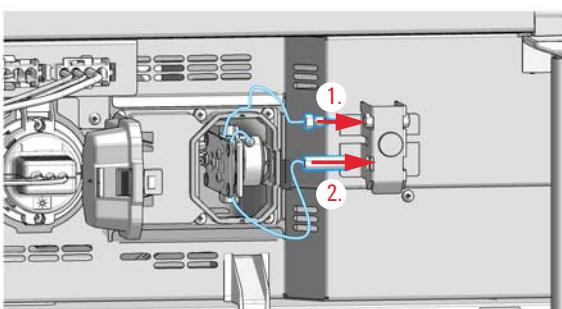
- 6** Remove the flow cell while pressing the flow cell holder.



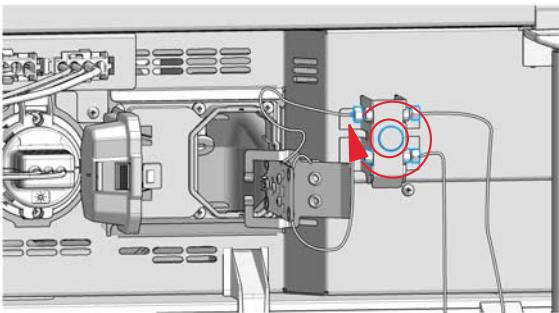
- 7** Insert the flow cell while pressing the flow cell holder.



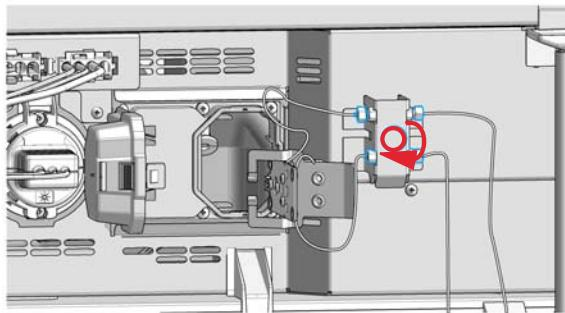
- 8** Insert the flow cell capillaries into the union holder (top is inlet, bottom is outlet).



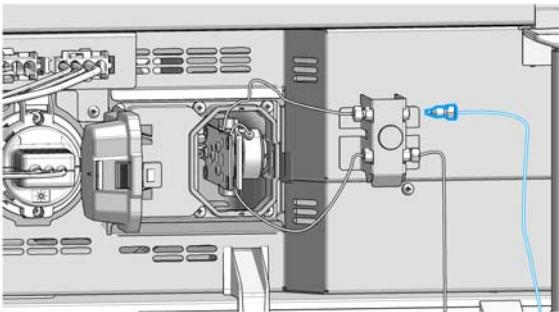
9 Tighten the thumb screw



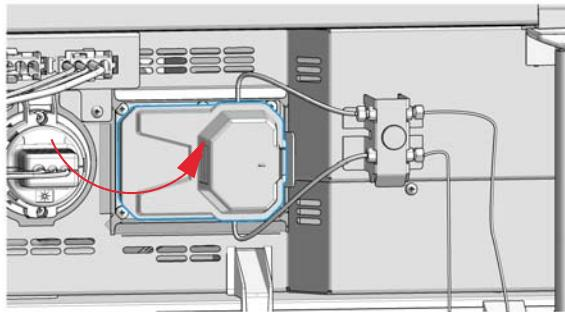
10 Reconnect the waste tubing (bottom) to the union



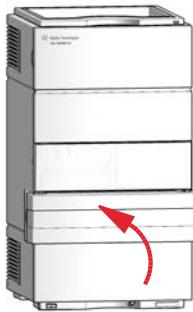
11 Reconnect the flow cell inlet capillary (top) to the union.



12 Close the flow cell door.



13 Close the lower front cover.

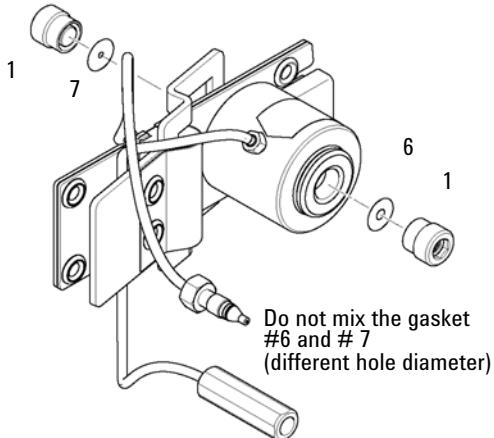


14 Perform a Wavelength Verification and Calibration or a Holmium Oxide Test to check the correct positioning of the flow cell.

Maintenance of the Flow Cell

When	If the flow cell needs repair due to leaks or contaminations (reduced light throughput)
Tools required	Description Wrench, 1/4 inch for capillary connections Hexagonal key, 4 mm Toothpick
Parts required	Description For parts, see “ Standard Flow Cell ” on page 352.
Preparations	Turn the flow off. Remove the front cover. Remove the flow cell, see “ Remove and Install a Flow Cell ” on page 307.

- 1** Use a 4 mm hex key to unscrew the window assembly [1] and remove the gasket [2] from the cell body.



NOTE

Carefully take one of the gaskets (#6 back or # 7 front) and insert it into the cell body.

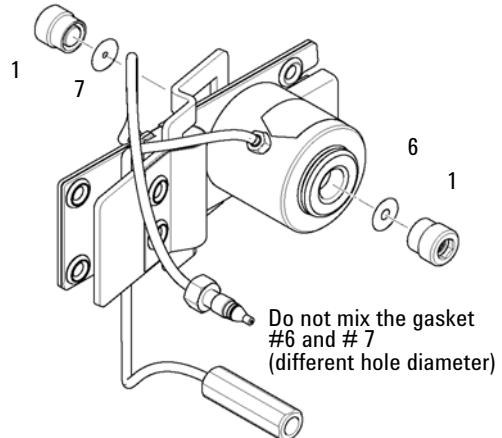
Do not mix the gasket #6 and # 7.

Gasket # 7 has the smaller hole and must be on the light entrance side.

Verify that the gasket is positioned flat on the bottom and the light path is not blocked.

If you removed all individual parts from the window assembly refer to the figures in "[Standard Flow Cell](#)" on page 352 for the correct orientation of the parts.

- 2** Insert the window assembly [1] into the cell body.



11 Maintenance and Repair

Diode Array Detector (DAD)

Next Steps:

- 3** Using a 4-mm hex key, tighten the window screw hand tight plus a quarter turn.
- 4** Reconnect the capillaries, see “[Remove and Install a Flow Cell](#)” on page 307.
- 5** Perform a leak test.
- 6** Insert the flow cell.
- 7** Replace the front cover
- 8** Perform a “[Wavelength Verification and Recalibration](#)” on page 188 or a “[Holmium Oxide Test](#)” on page 180 to check the correct positioning of the flow cell.

Replacing Capillaries on a Standard Flow Cell

When If the capillary is blocked

Tools required

Description
Wrench, 1/4 inch for capillary connections
Wrench, 4 mm (for capillary connections)
Screwdriver, Pozidriv #1 PT3

Parts required

Description
For parts see “ Standard Flow Cell ” on page 352.

Preparations

Turn the lamp(s) off.
Remove the front cover.
Remove the flow cell, see “ Remove and Install a Flow Cell ” on page 307.

NOTE All descriptions in this procedure are based on the default orientation of the cell (as it is manufactured). The heat exchanger/capillary and the cell body can be fixed mirror symmetrically to have both capillaries routed to the bottom or to the top (depending on the routing of the capillaries to the column).

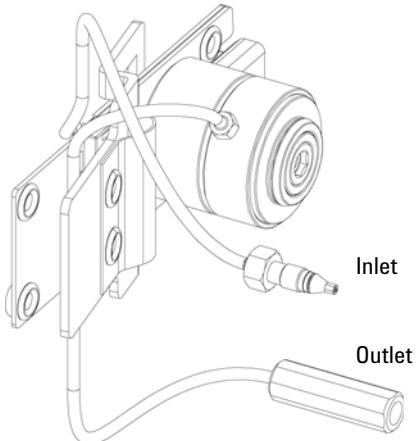
NOTE The fittings at the flow cell body are special types for low dead volumes and not compatible with other fittings.

When retightening the fittings, make sure that they are carefully tightened (handtight plus 1/4 turn with a wrench). Otherwise damage of the flow cell body or blockage may result.

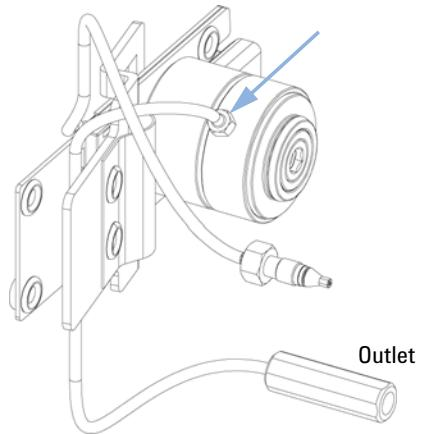
11 Maintenance and Repair

Diode Array Detector (DAD)

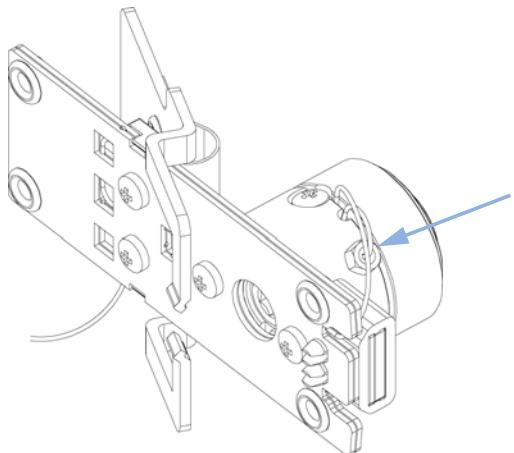
- 1** Identify the inlet and outlet capillaries. To replace the inlet capillary, continue with step "To replace the inlet capillary, use a 4-mm wrench for the fitting."



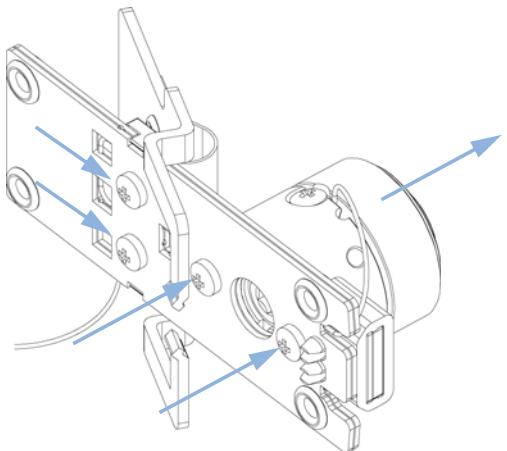
- 2** After replacing the outlet capillary, fix it handtight first. Then do a 1/4 turn with a 4-mm wrench.



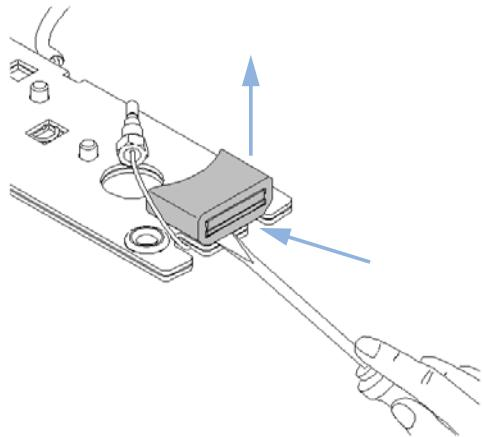
- 3** To replace the inlet capillary, use a 4-mm wrench for the fitting.



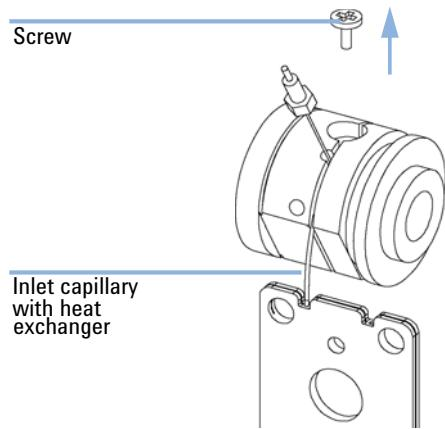
- 4** Unscrew the cell body from the heat exchanger and the heat exchanger from the clamp unit.



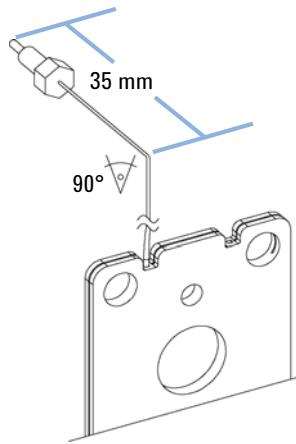
- 5 Use a small flat screw driver to carefully lift off the I.D. tag. Shown is the default orientation. See Note at the beginning of this section.



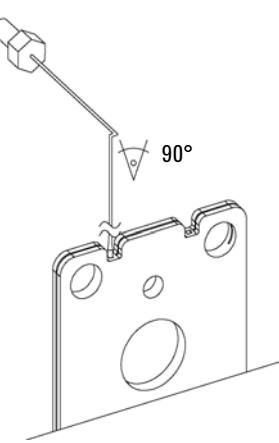
- 6 Unscrew the fixing screw and unwrap the inlet capillary from the groove in the flow cell body.



- 7 Take the new inlet capillary and bend it 90° about 35 mm from its end.



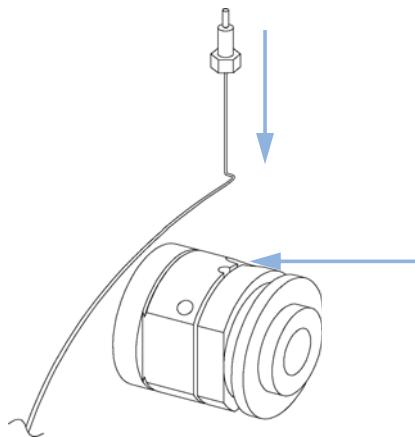
- 8 Bend the capillary again by 90° as shown below.



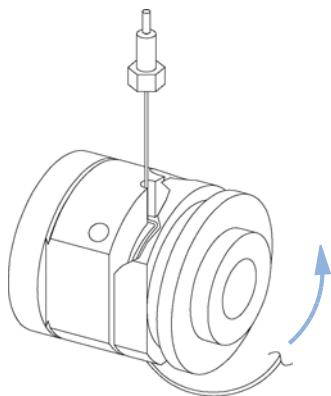
11 Maintenance and Repair

Diode Array Detector (DAD)

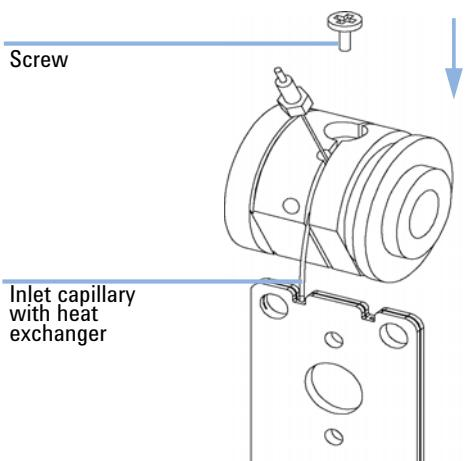
- 9** Insert the capillary into the hole between fixing screw and the inlet fitting.



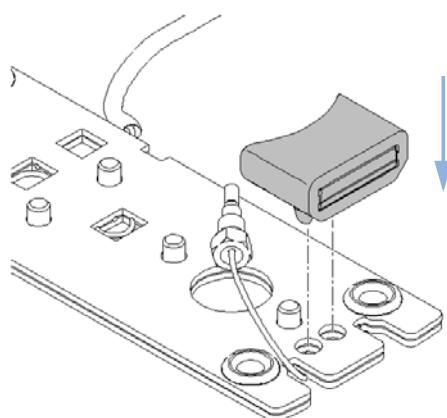
- 10** The capillary lays in the groove and should be tied around the body (in the groove) 5 times.



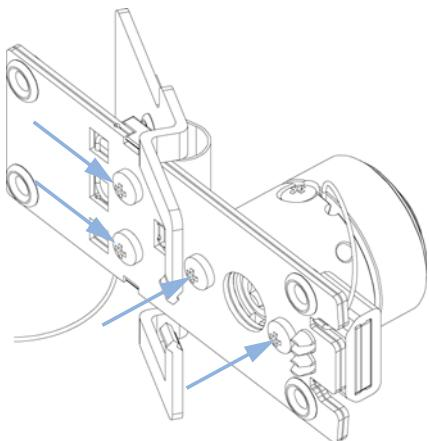
- 11** Insert the fixing screw, so that the capillary cannot leave the groove.



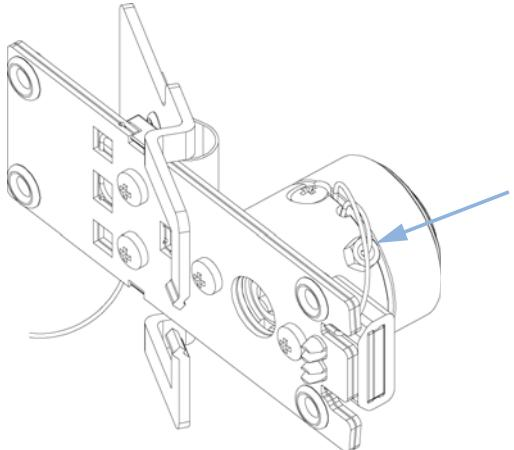
- 12** Carefully insert the I.D. tag into the new heat exchanger. Shown is the default orientation. See Note at the beginning of this section.



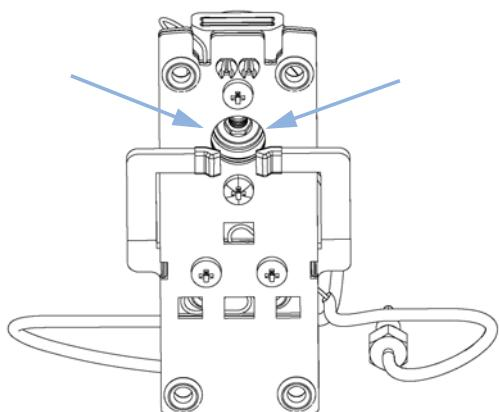
- 13** Fix the heat exchanger to the clamp unit and the flow cell body to the heat exchanger.



- 14** Fix the inlet capillary to the flow cell body handtight first. Then do a 1/4 turn with a 4-mm wrench.



- 15** Check for a centered holder vs. hole. If required adjust with the holder screws.



Next Steps:

16 Reconnect the capillaries, see “[Remove and Install a Flow Cell](#)” on page 307.

17 Perform a leak test.

18 Insert the flow cell.

19 Replace the front cover.

20 Perform a “[Wavelength Verification and Recalibration](#)” on page 188 or a “[Holmium Oxide Test](#)” on page 180 to check the correct positioning of the flow cell.

Cleaning or Exchanging the Holmium Oxide Filter

When If holmium oxide filter is contaminated

Tools required

Screwdriver, Pozidriv #1 PT3
Screwdriver, flat blade
Wrench, 1/4 inch for capillary connections
Pair of tweezers

Parts required

#	p/n	Description
1	79880-22711	Holmium oxide filter

Preparations

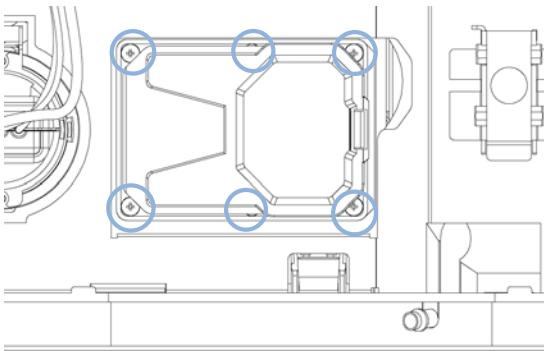
- Turn the lamp(s) off.
- Remove the front cover.
- Remove the flow cell, see “[Remove and Install a Flow Cell](#)” on page 307.

NOTE

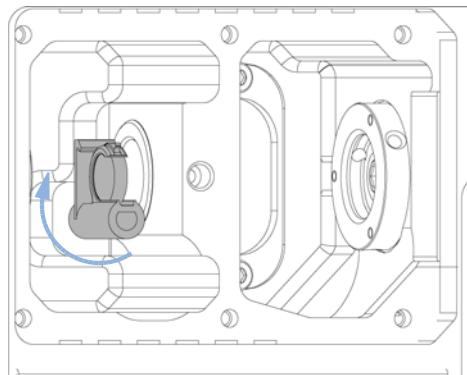
See also “[Declaration of Conformity for HOX2 Filter](#)” on page 389.

The glass tends to build a film on its surface even under normal environmental conditions. This is a phenomenon, which can be found also on the surface of several other glasses and has something to do with the composition of the glass. There is no indication, that the film has an influence on the measurement. Even in the case of a thick film, which scatters the light remarkably, no shift of the peak positions is to be expected. A slight change in the absorbance might be possible. Other components within the light path (lenses, windows,...) are also changing their behavior over the time.

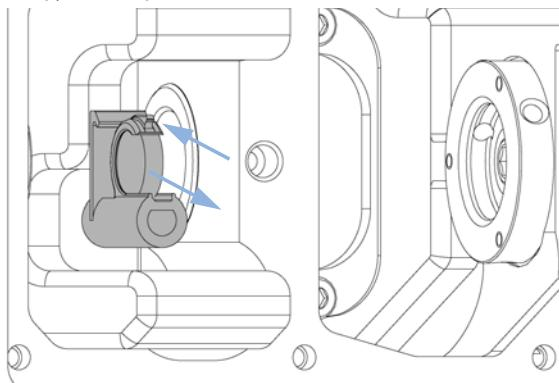
- 1** Unscrew the six screws and remove the flow cell cover.



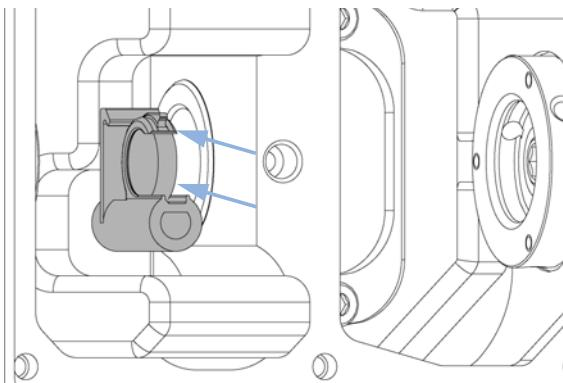
- 2** If not already in this position, move the filter up.



- 3** While releasing the holder with a screw driver (at the top), carefully remove the holmium oxide filter.



- 4** While releasing the holder with a screw driver, carefully insert the holmium oxide filter.



NOTE

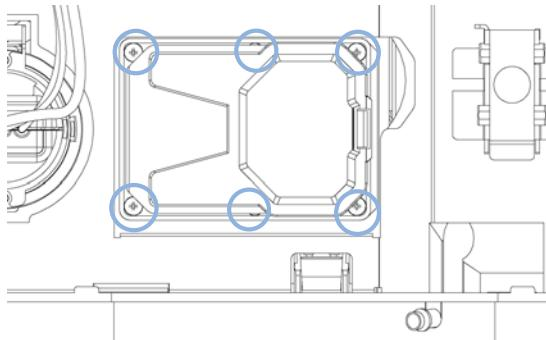
Do not scratch the holmium oxide filter.

The holmium oxide filter can be cleaned with alcohol and a lint-free cloth.

11 Maintenance and Repair

Diode Array Detector (DAD)

- 5** Replace the flow cell cover and fix the six screws.



Next Steps:

- 6** Perform a holmium oxide test, see "[Holmium Oxide Test](#)" on page 180 to check the proper function of the holmium oxide filter.
- 7** Insert the flow cell, see "[Remove and Install a Flow Cell](#)" on page 307.
- 8** Replace the front cover.
- 9** Turn on the flow.

Correcting Leaks

When If a leakage has occurred in the flow cell area or at the heat exchanger or at the capillary connections

Tools required	p/n	Description
		Tissue
		Wrench, 1/4 inch for capillary connections
	5043-0915	Fitting mounting tool for bio-inert capillaries

Preparations Remove the front cover.

- 1 Use tissue to dry the leak sensor area and the leak pan.
- 2 Observe the capillary connections and the flow cell area for leaks and correct, if required.

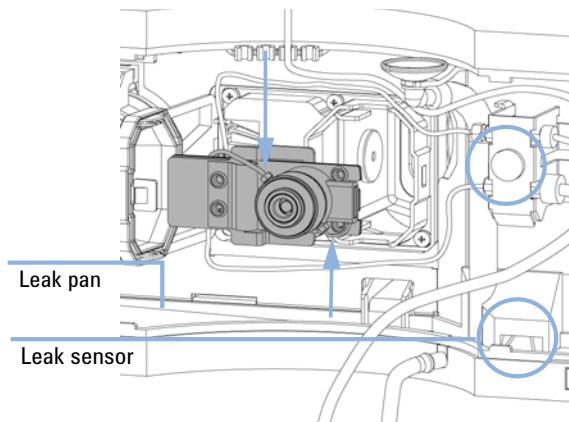


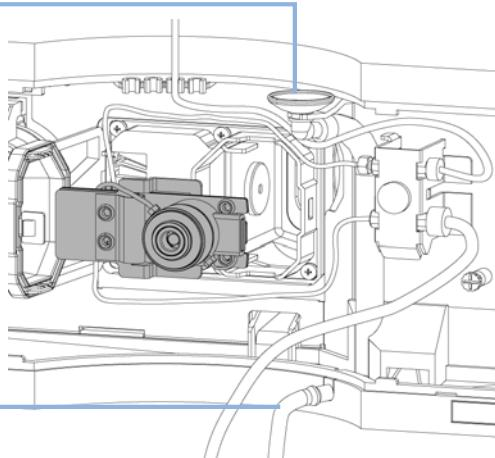
Figure 66 Observing for Leaks

- 3 Replace the front cover.

Replacing Leak Handling System Parts

When	If the parts are corroded or broken		
Tools required	None		
Parts required	#	p/n	Description
	1	5041-8388	Leak funnel
	1	5041-8389	Leak funnel holder
	1	5062-2463	Corrugated tubing, PP, 6.5 mm id, 5 m
Preparations	Remove the front cover.		
	<ol style="list-style-type: none">1 Pull the leak funnel out of the leak funnel holder.2 Pull out the leak funnel with the tubing.3 Insert the leak funnel with the tubing in its position.4 Insert the leak funnel into the leak funnel holder.		

Leak funnel
and holder



Leak tubing

Figure 67 Replacing Leak Handling System Parts

- 5 Replace the front cover.

Algae Growth in HPLC Systems

The presence of algae in HPLC systems can cause a variety of problems that may be incorrectly diagnosed as instrument or application problems. Algae grow in aqueous media, preferably in a pH range of 4 – 8. Their growth is accelerated by buffers, for example phosphate or acetate. Since algae grow through photosynthesis, light also stimulates their growth. Small algae growth is seen even in distilled water after some time.

Instrumental problems associated with algae

Algae deposit and grow everywhere within the HPLC system causing:

- deposits on ball valves, inlet or outlet, resulting in unstable flow or total failure of the pump.
- small-pore solvent inlet filters to plug, resulting in unstable flow or total failure of the pump.
- small-pore high pressure solvent filters (usually placed before the injector) to plug, resulting in high system pressure.
- column filters to plug, resulting in high system pressure.
- flow cell windows of detectors to become dirty, resulting in higher noise levels. Since the detector is the last module in the flow path, this problem is less common.

Symptoms observed with the Agilent 1220 Infinity II LC

In contrast to other HPLC systems, such as HP 1090 and HP 1050 Series, which use helium degassing, algae have a better chance to grow in systems such as the Agilent 1220 Infinity II LC, where helium is not used for degassing (most algae need oxygen and light for growth).

The presence of algae in the Agilent 1220 Infinity II LC can cause the following to occur:

- Blockage of PTFE frits (pack of 5) (01018-22707) (purge valve assembly) and column filter, causing increased system pressure. Algae appear as white or yellowish-white deposits on filters. Typically, black particles from the normal wear of the piston seals do not cause the PTFE frits to block over short-term use.
- Short lifetime of solvent filters (bottle head assembly). A blocked solvent filter in the bottle, especially when only partly blocked, is more difficult to identify and may show up as problems with gradient performance, intermittent pressure fluctuations etc.
- Algae growth may also be the possible source of failures of the ball valves and other components in the flow path.

Preventing and/or reducing the algae problem

- Always use freshly prepared solvents, especially use demineralized water that has been filtered through about 0.2 µm filters.
- Never leave mobile phase in the instrument for several days without flow.
- Always discard old mobile phase.
- Use the Solvent bottle, amber (9301-1450) supplied with the instrument for your aqueous mobile phase.
- If possible, add a few mg/l sodium azide or a few percent organic solvent to the aqueous mobile phase.

Replacing the Module's Firmware

- When**
- The installation of newer firmware might be necessary
- if a newer version solves problems of older versions or
 - to keep all systems on the same (validated) revision.
- The installation of older firmware might be necessary
- to keep all systems on the same (validated) revision or
 - if a new module with newer firmware is added to a system or
 - if third party control software requires a special version.

Tools required	Description
OR	LAN/RS-232 Firmware Update Tool
OR	Agilent Lab Advisor software
	Instant Pilot G4208A (only if supported by module)

Parts required	#	Description
	1	Firmware, tools and documentation from Agilent web site

Preparations Read update documentation provided with the Firmware Update Tool.

To upgrade/downgrade the module's firmware carry out the following steps:

- 1 Download the required module firmware, the latest LAN/RS-232 FW Update Tool and the documentation from the Agilent web.
 - <http://www.agilent.com/en-us/firmwareDownload?whid=69761>
- 2 For loading the firmware into the module follow the instructions in the documentation.

Module Specific Information

There is no specific information for this module.

11 Maintenance and Repair

Replacing the Module's Firmware

12

Parts for Maintenance and Repair

1220 Infinity II LC System	328
System Parts	328
Mainboard overview	329
Fuses	330
Solvent Delivery System	331
Injection System	340
Manual Injector	340
Autosampler	342
Column Oven	347
Detector	348
Variable Wavelength Detector (VWD)	348
Diode Array Detector (DAD)	352

This chapter provides information on parts for maintenance and repair.



Agilent Technologies

12 Parts for Maintenance and Repair

1220 Infinity II LC System

1220 Infinity II LC System

System Parts

System part numbers

p/n	Description
0950-4997	Power supply
G4280-65850	Agilent 1220 Infinity II LC main board
G4280-65800	Agilent 1220 Infinity II LC VL main board
G4280-68723	Cabinet kit
G4280-60056	Side cover kit, walk-up
G4280-60102	Front door, top
G4280-60001	Front door, bottom
G4280-60014	Main Power Switch and Leak Shutdown Board
G4280-65802	FSL board (Status LED board)
5067-5378	Connecting tube, DCGV to PIV
G4280-80004	Fan
8121-1833	Power Switch Cable
G4900-60800	Primary Inlet Filter Assy
G4280-81602	Cable, status LED
G4280-81620	Temp Sensor Cable
G4280-40007	Light pipe, status
G4280-40016	Power Switch Coupler ZL
G4280-44013	Leak plane, man. inj.
G4280-44500	Leak panel, bottom
G4280-44501	Leak plane, pump
G4280-44502	Leak plane, ALS
G4280-44016	Holder, temp. sensor
5061-3356	Leak Sensor Assembly

Mainboard overview

Pump module connectors:

1. Instrument Status LED
2. Pressure Sensor
3. Dual Channes Gradient Valve
4. Pump Drive Encoder
5. Pump Fan
6. Solvent Selection Valve (SSV)
7. Active Inlet Valve (not used)
8. Degasser Assembly Connectors
9. Pump Drive

Sampler module connectors:

10. Connector for Manual Injector
- 11 Connectors for autosampler transport assembly

12. Heater assembly connector

VWD connectors:

13. Grating motor
14. Filter motor
15. Grating pos. sensor
16. Filter pos. sensor
17. Grounding connector lamp
18. Sampler diode
19. Reference diode
20. D2 lamp
21. VWD fan
22. VWD heater sensor
23. VWD heater

Power supply sensors

24. Power fail connector
25. Power switch control
26. Power and ground

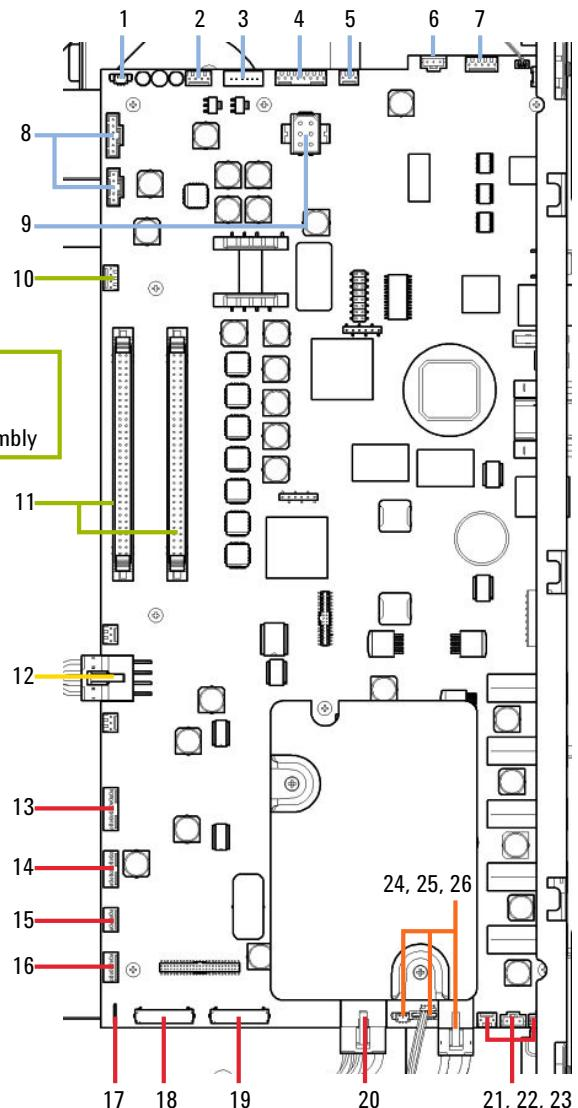


Figure 68 Connectors of the Agilent 1220 Infinity II Instrument Mainboard

12 Parts for Maintenance and Repair

1220 Infinity II LC System

Fuses

5 mainboard fuses

p/n	Description
2110-1417	Fuse 3.15 A 250 V

NOTE

Next to each fuse is an LED. Red LED indicates the fuse is blown. If one of the fuses is blown, the green LED of the power switch flashes.

- Fuse F1 (degasser, pump, injector motors)
- Fuse F2 (injector sensors, column oven, ext. 24 V connector)
- Fuse F3 (processor core, +5 V, +15 V, -15 V supply on mainboard)
- Fuse F4 (VWD incl. D2-lamp)
- Fuse F5 (VWD heater, fan)

1 fuse netfilter

p/n	Description
2110-1004	Fuse 10 A t

Solvent Delivery System

Pump Main Repair Parts

p/n	Description
G1311-60001	Pump drive
G1311-60065	Damper Isocratic/Quaternary Pump 600 bar
79835-60005	Damper Quaternary Pump 400 bar
G4280-60004	Dual-channel gradient valve
3160-1017	Fan 92 mm x 92 mm
G4280-68708	SSV

12 Parts for Maintenance and Repair

Solvent Delivery System

Pump Head Assembly Without Seal Wash

Item	p/n	Description
	G1312-60056	Pump Head 1200 SL without Seal Wash
1	5063-6586	Sapphire piston
2	G1311-60002	Piston housing
3	5067-1560	Support Ring SL, no seal wash
4	5062-2484	Gasket, seal wash (pack of 6)
5	5042-8952	Seal holder
6	5063-6589	Piston seal PTFE, carbon filled, black (pack of 2), default
OR	0905-1420	PE seals (pack of 2)
7	G1311-25200	Pump chamber housing
8	G1312-60066	Passive inlet valve 1220/1260
9	G1312-60067	Outlet valve 1220/1260
10	5042-1303	Lock screw
11	G4280-60061	Purge valve
12	0515-2118	Pump head screw (M5, 60 mm)

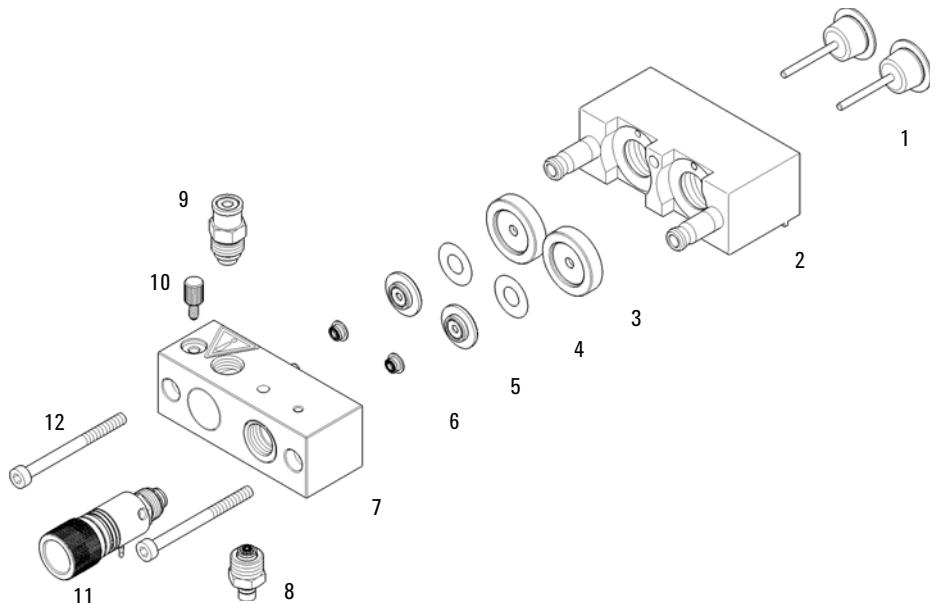


Figure 69 Pump head assembly without seal wash option

The Pump Head 1200 SL without Seal Wash (G1312-60056) includes items 1-7, 10 and 12.

12 Parts for Maintenance and Repair

Solvent Delivery System

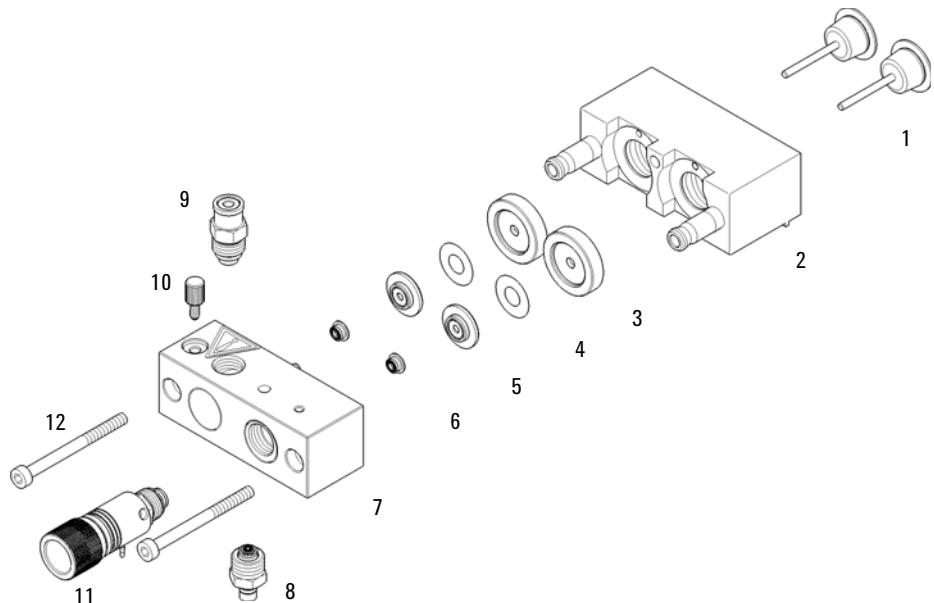
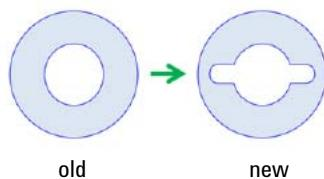


Figure 70 Pump head assembly without seal wash option

NOTE

The design of the seal wash gasket has changed, see below.



Pump Head Assembly With Seal Wash Option

Item	p/n	Description
	G1312-60045	Pump head assembly with seal wash
1	5063-6586	Sapphire piston
2	G1311-60002	Piston housing
3	01018-60027	Support ring seal wash
4	0905-1175	Wash seal (PTFE)
OR	0905-1718	Wash Seal PE
	5065-9978	Tubing, 1 mm i.d., 3 mm o.d., silicone, 5 m
5	5062-2484	Gasket, seal wash (pack of 6)
6	5042-8952	Seal holder
7	5063-6589	Piston seal PTFE, carbon filled, black (pack of 2), default
	0905-1420	PE seals (pack of 2)
8	G1311-25200	Pump chamber housing
9	G1312-60066	Passive inlet valve 1220/1260
10	G1312-60067	Outlet valve 1220/1260
11	5042-1303	Lock screw
12	G4280-60061	Purge valve
13	0515-2118	Pump head screw (M5, 60 mm)
	01018-23702	Insert tool

The Pump head assembly with seal wash (G1312-60045) includes items 1-8, 11 and 13.

NOTE

Active seal wash is not supported for the 1220 Infinity II LC, only continuous seal wash.

12 Parts for Maintenance and Repair

Solvent Delivery System

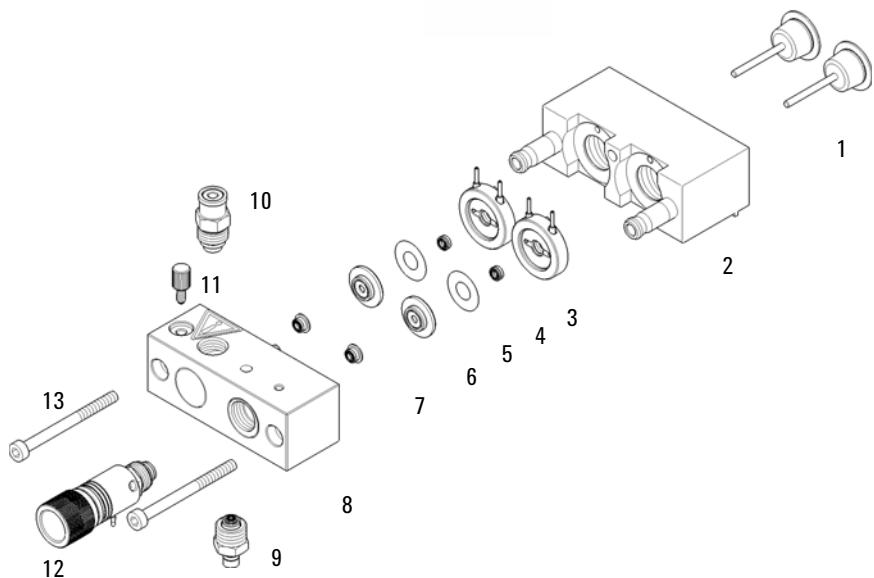
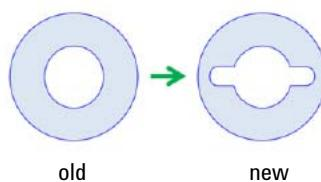


Figure 71 Pump Head with Seal Wash Option

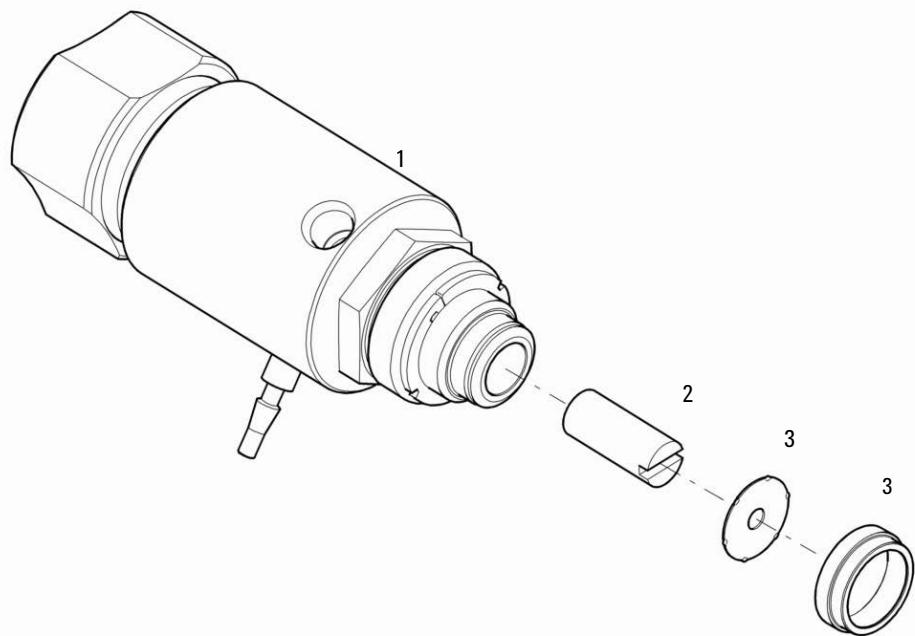
NOTE

The design of the seal wash gasket has changed, see below.



Purge Valve Assembly

Item	p/n	Description
	G4280-60061	Purge valve
1		Valve body
2	01018-22707	PTFE frits (pack of 5)
3	5067-4728	Seal Cap Assembly



Degasser Unit

p/n	Description
G4280-60070	Degassing unit
G1311-60007	Vacuum Pump for Quaternary Pump Degasser

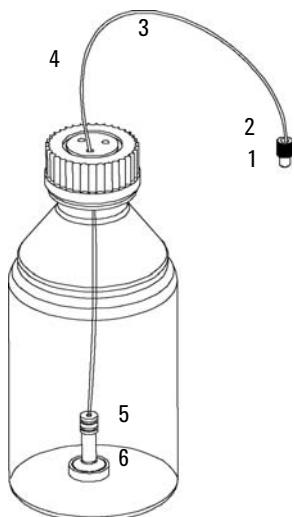
12 Parts for Maintenance and Repair

Solvent Delivery System

p/n	Description
G4280-67300	Vacuum Tubing Kit
5067-4734	Degasser Chamber (1 Channel) for G1311-60070

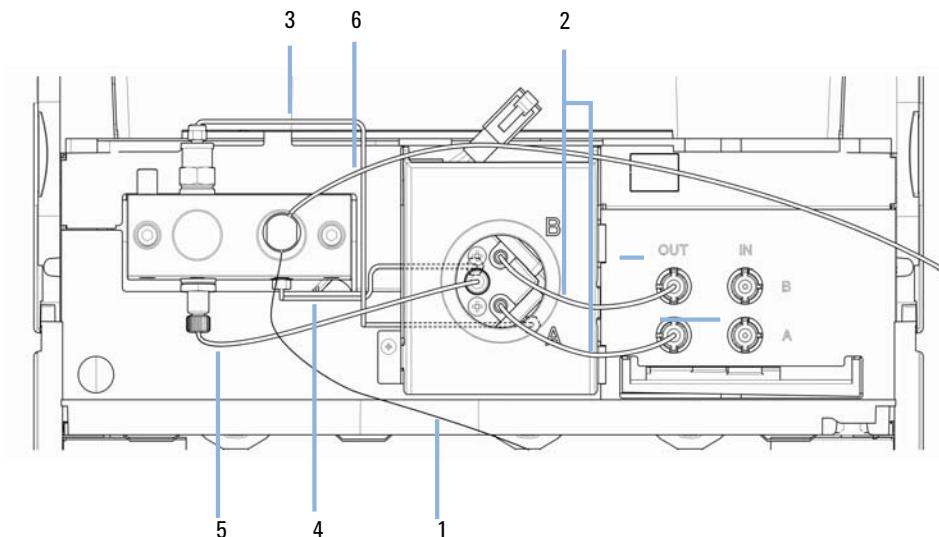
Bottle Head Assembly

Item	p/n	Description
	G1311-60003	Bottle-head assembly
1	5063-6598	Ferrules with lock ring (10/Pk)
2	5063-6599	Tube screw (10/Pk)
3		Wire marker
4	5062-2483	Tube PTFE 1.5 mm x 5 m, 3 mm od
5	5062-8517	Frit adapter (pack of 4)
6	5041-2168	Solvent inlet filter, 20 µm pore size



Hydraulic Path

Item	p/n	Description
1	G1312-67305	Outlet capillary, pump to injector
	G1311-60003	Bottle head assembly, bottle to passive inlet valve or vacuum degasser
2	G4280-60034	Solvent tube, vacuum degasser to DCGV
3	G4280-81300	Capillary, plunger 1 to damper
4	G4280-81301	Capillary, damper to plunger 2
	5067-5378	Connecting tube, DCGV to PIV
6	5062-2461	Waste tube, 5 m (reorder pack)

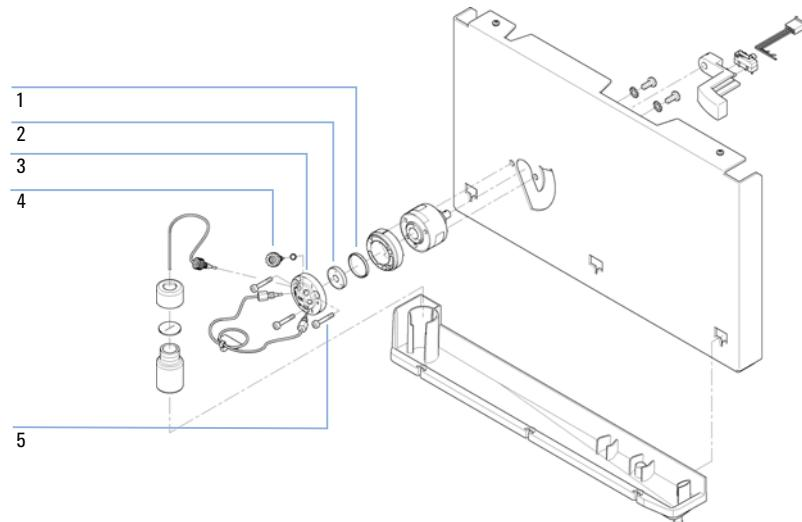


Injection System

Manual Injector

Injection Valve Assembly

Item	p/n	Description
	5067-4202	Manual injection valve 600 bar, complete including loop and needle port
	1535-4045	Isolation seal
2	5068-0082	Rotor seal, PEEK
3	0101-1417	Stator head
4	5067-1581	Needle port
5	5068-0018	Stator screws
	8710-0060	Hex-key wrench, 9/64 inch



Sample Loops

Sample loops stainless steel

p/n	Description
0101-1248	Sample loop 5 µL
0100-1923	Sample loop 10 µL
0100-1922	Sample loop 20 µL
0100-1924	Sample loop 50 µL
0100-1921	Sample loop 100 µL
0101-1247	Sample loop 200 µL
0101-1246	Sample loop 500 µL
0101-1245	Sample loop 1 mL
0101-1244	Sample loop 2 mL

Sample loops PEEK

p/n	Description
0101-1241	Sample loop 5 µL
0101-1240	Sample loop 10 µL
0101-1239	Sample loop 20 µL
0101-1238	Sample loop 50 µL
0101-1242	Sample loop 100 µL
0101-1227	Sample loop 200 µL
0101-1236	Sample loop 500 µL
0101-1235	Sample loop 1 mL
0101-1234	Sample loop 2 mL

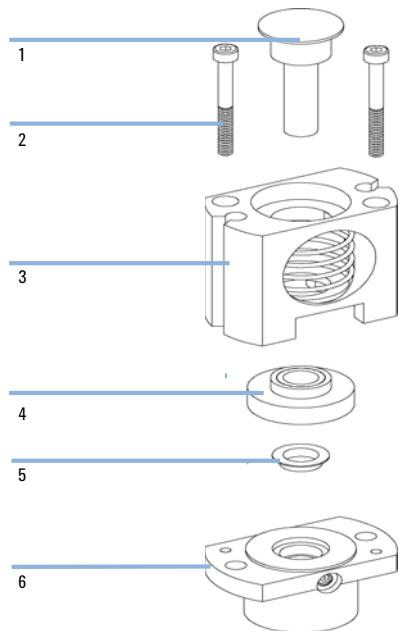
Autosampler

Autosampler Main Assemblies

p/n	Description
G1329-60009	Transport assembly
G4280-60027	Sampling unit assembly (excluding injection valve and analytical head)
01078-60003	Analytical head assembly, 100 µL
0101-1422	Injection valve
G1313-44510	Vial tray
G1313-60010	Gripper assembly
G4280-87304	Waste capillary
G4280-81615	Cable, sampling unit
G4280-81616	Cable, sample transport
5067-1581	Needle port

Analytical Head Assembly

Item	p/n	Description
	01078-60003	Analytical head assembly, 100 µL
1	5063-6586	Sapphire piston
2	0515-0850	Screw M4, 40 mm long
3	01078-23202	Adapter
4	5001-3739	Support Seal assembly
5	5063-6589	Metering seal (pack of 2) for 100 µL analytical head
6	01078-27710	Head body
	0515-2118	Screw M5, 60 mm long, for mounting assembly

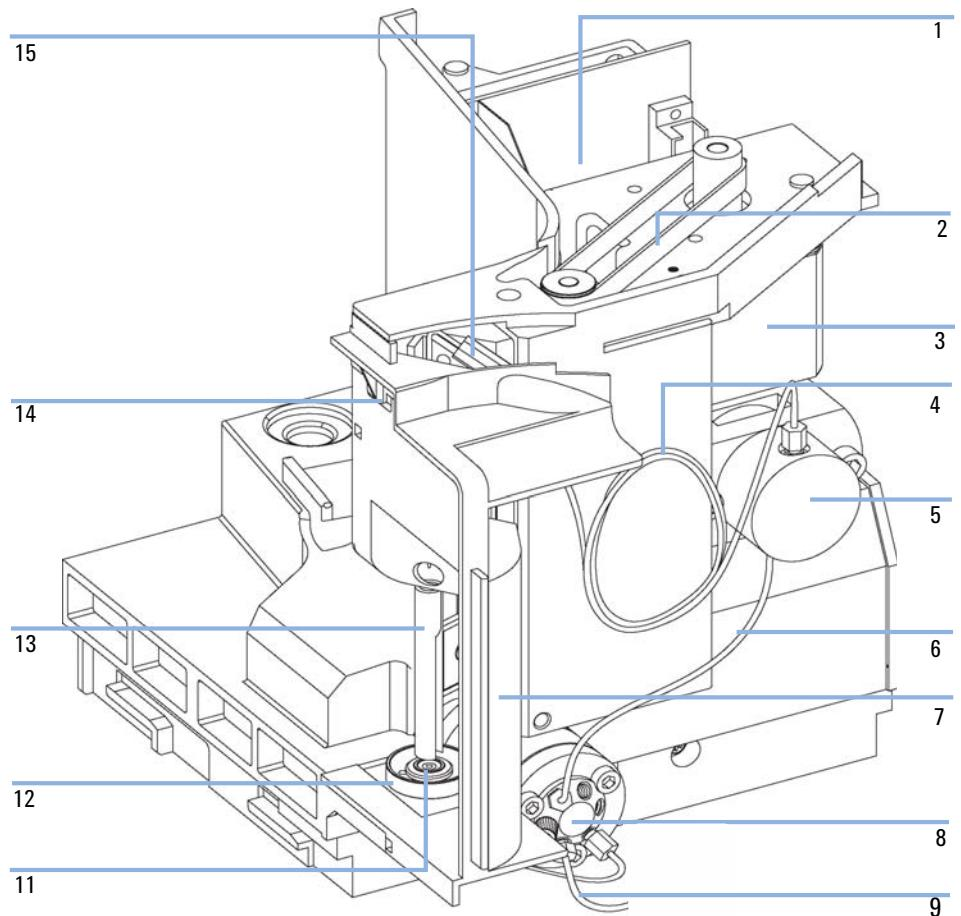


12 Parts for Maintenance and Repair

Injection System

Sampling Unit Assembly

Item	p/n	Description
	G4280-60027	Sampling unit assembly (excluding injection valve and analytical head)
1	G1313-65803	SUD board
2	1500-0697	Belt gear for metering unit and needle arm
3	5062-8590	Stepper motor for metering unit and needle arm
4	01078-87302	Loop capillary (100 µL)
5	01078-60003	Analytical head assembly, 100 µL
6	G1313-87301	Capillary, injection valve to analytical head (160 mm × 0.25 mm)
7	G1329-44115	Safety cover
8	0101-1422	Injection valve
9	G1313-87300	Waste tube injection valve assy (120 mm)
11	G1313-87101	Needle-seat assy (0.17 mm i.d. 2.3 µL)
12	G1313-43204	Seat adapter
13	G1313-44106	Safety flap
14	G1313-68715	Flex board
15	G1313-87201	Needle assembly
	G1313-68713	Clamp Kit (includes needle clamp and 2 x clamp screw)

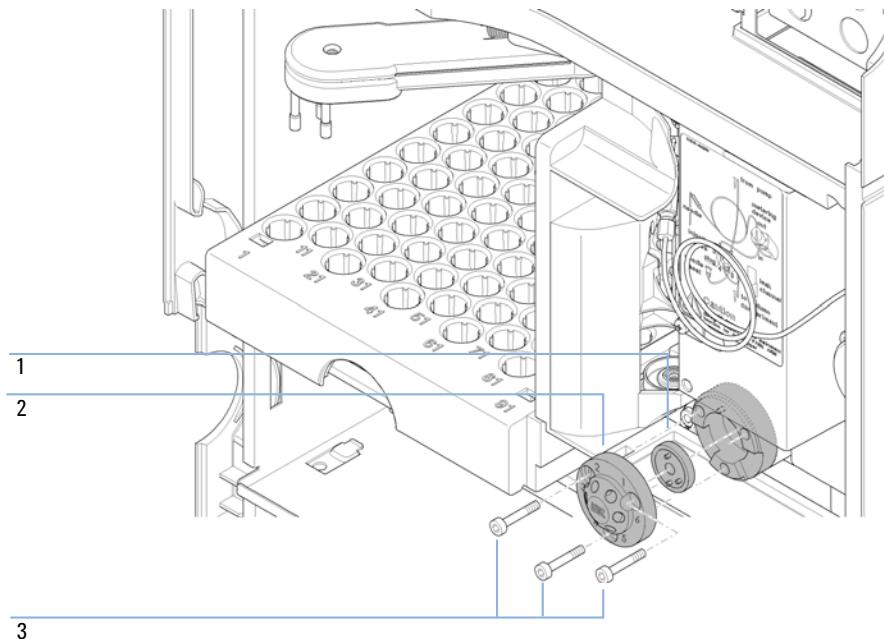


12 Parts for Maintenance and Repair

Injection System

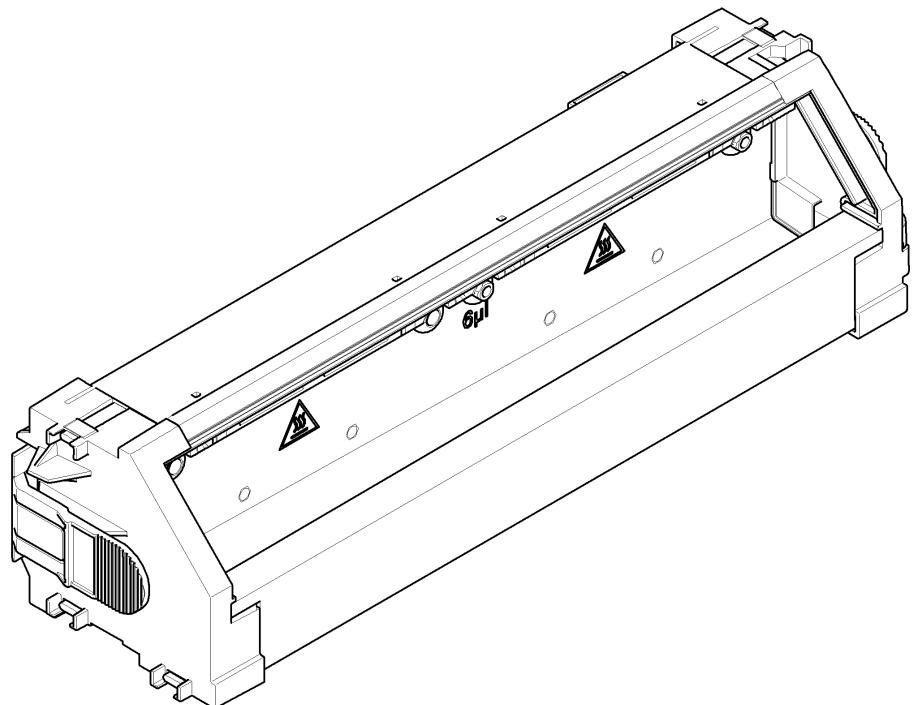
Injection Valve Assembly

Item	p/n	Description
	0101-1422	Injection valve
	0100-1852	Isolation seal
1	0101-1416	Rotor seal (PEEK)
2	0101-1417	Stator head
3	1535-4857	Stator screws



Column Oven

p/n	Description
G4280-60040	Complete column oven assembly
G4280-60017	Heater door assembly



12 Parts for Maintenance and Repair

Detector

Detector

Variable Wavelength Detector (VWD)

VWD Main Repair Parts

p/n	Description
G1314-60061	Complete optical unit assembly
G1314-69061	Optical Unit (exchange part)
3160-1016	Fan 119 mm x 119 mm
G4280-81607	Lamp cable
G4280-00013	Z-PLANE TOP VWD
G1314-60101	Deuterium lamp (with RFID tag)
G1314-60100	Deuterium lamp
G1314-60186	Standard flow cell 10 mm, 14 µL, 40 bar (with RFID tag)
G1314-60187	Micro flow cell 3 mm, 2 µL, 120 bar (with RFID tag)
G1314-60183	Semi-micro flow cell 6 mm, 5 µL (with RFID tag)
G1314-60182	High pressure flow cell 10 mm, 14 µL, 400 bar (with RFID tag)

Standard Flow Cell 10 mm / 14 µL

Item	p/n	Description
	G1314-60086	Standard flow cell, 10 mm, 14 µL, 40 bar
	5062-8522	Capillary column - detector PEEK 600 mm lg, 0.17 mm i.d., 1/16 inch o.d.

Item	p/n	Description
	G1314-65061	Cell Repair Kit, includes 2x Gasket #1, 2x Gasket #2, 2x Window Quartz
1	G1314-65062	Cell screw kit
2	79853-29100	Conical spring kit, 10/pk
3	G1314-65066	Ring #2 kit (IN small hole, i.d. 1 mm) PEEK, 2/pk
4	G1314-65064	Gaskets #2 IN (small hole i.d. 1 mm), KAPTON 10/pk
5	79853-68742	Window quartz kit, 2/pk
6	G1314-65063	Gasket #1 kit (OUT large hole, i.d. 2.4 mm) KAPTON, 2/pk
7	G1314-65065	Ring #1 kit (OUT large hole, i.d. 2.4 mm) PEEK, 2/pk

- 1 Cell screw
- 2 Conical springs
- 3 Ring #2 IN
- 4 Gasket #2 IN
- 5 Quartz window
- 6 Gasket #1 OUT
- 7 Ring #1 OUT

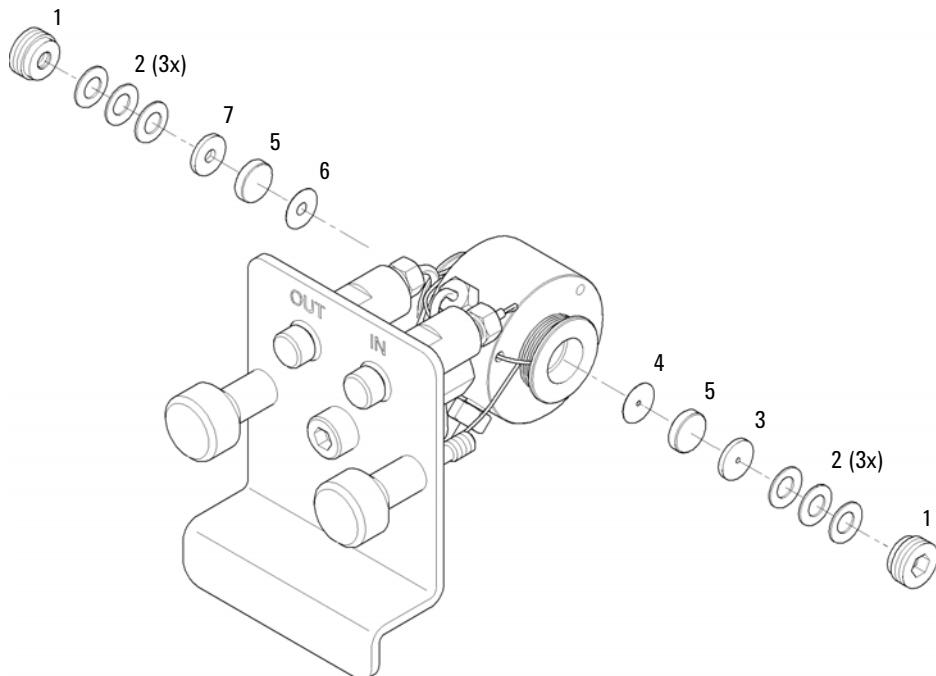


Figure 72 Standard Flow Cell 10 mm / 14 μL

12 Parts for Maintenance and Repair

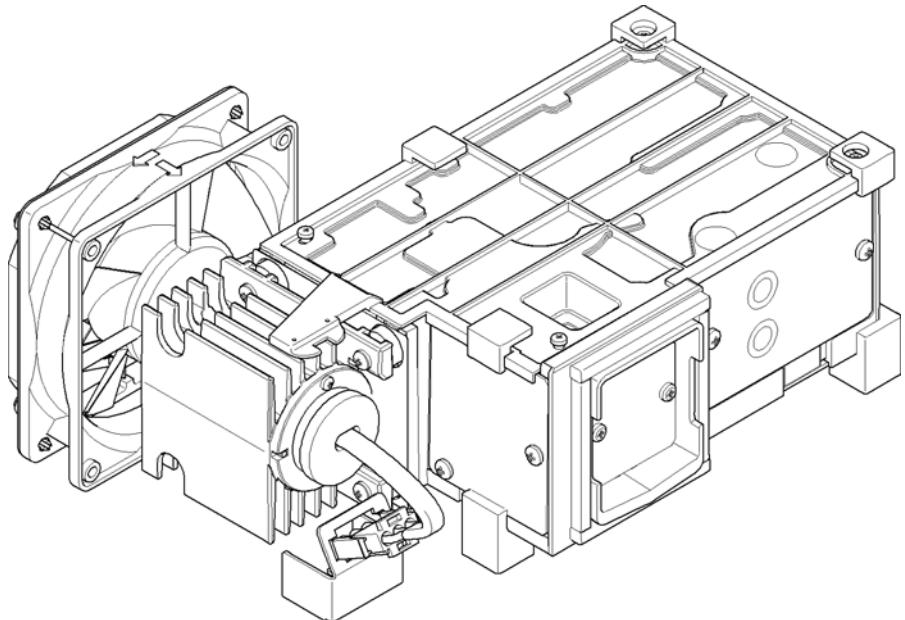
Detector

Detector Lamp

p/n	Description
G1314-60100	Deuterium lamp

Optical Unit and Fan Assembly

p/n	Description
G1314-60061	Complete optical unit assembly
G4280-80004	Fan
G1314-60113	Heater Assembly
G1314-81010	Heater Cable
G1314-67000	Heater Interface Board Kit (includes 4 rivets)
G1314-65802	VWD temp. sensor board
G4280-81623	Cable Front end VWD Reference
G4280-81619	Cable Front end VWD Sample



NOTE

Repairs to the optical unit require specialist knowledge.

12 Parts for Maintenance and Repair

Detector

Diode Array Detector (DAD)

Standard Flow Cell

Item	p/n	Description
	G1315-60022	Standard flow cell, 10 mm, 13 µL, 120 bar (12 MPa)
1	79883-22402	Window screw
2	5062-8553	Washer kit (10/pk)
3	79883-28801	Compression washer
4	79883-22301	Window holder
5	1000-0488	Quartz window
6	G1315-68711	Gasket BACK (PTFE), 2.3 mm hole, outlet side (12/pk)
7	G1315-68710	Gasket FRONT (PTFE), 1.3 mm hole, inlet side (12/pk)
8		Window assembly (comprises window screw, spring washers, compression washer, window holder and quartz window)
	G1315-87321	Capillary IN (0.17 mm, 590 mm lg) including heat exchanger
10	G1315-87302	Capillary OUT (0.17 mm, 200 mm lg)
11	G1315-84910	Clamp unit
	0515-1056	Screw M 2.5, 4 mm lg for cell body/clamp
	5022-2184	Union ZDV
	G1315-68712	Cell repair kit STD includes window screw kit, 4 mm hexagonal wrench and seal kit
	79883-68703	Window screw kit, includes 2 quartz windows, 2 compression washers, 2 window holders, 2 window screws and 10 washers

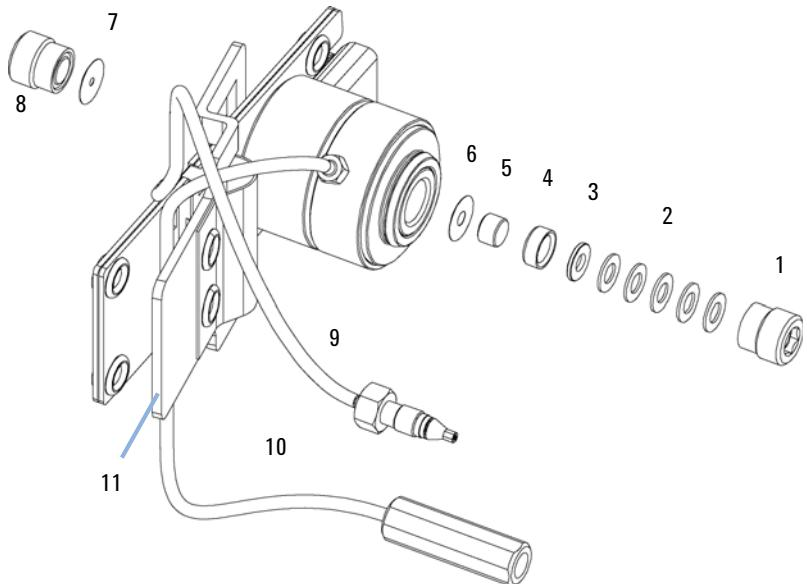


Figure 73 Standard Flow Cell Parts

NOTE

Gaskets # 6 and #7 have different hole diameters.

- 1 - window screw
- 2 - spring washers
- 3 - compression washer
- 4 - window holder
- 5 - quartz window
- 6 - Gasket

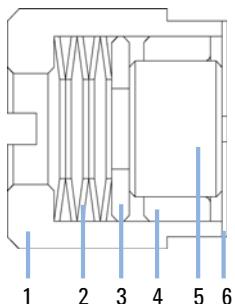


Figure 74 Orientation of Spring Washers

12 Parts for Maintenance and Repair

Detector

Detector Lamps

p/n	Description
2140-0820	Longlife Deuterium lamp "C" (with black cover and RFID tag)
G1103-60001	Tungsten lamp

13

Upgrading the 1220 Infinity II LC

Oven Upgrade 356

Install the (optional) External Tray 359

Install the Active Inlet Valve 363

This chapter provides information for upgrading the LC system.



Agilent Technologies

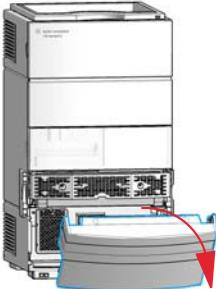
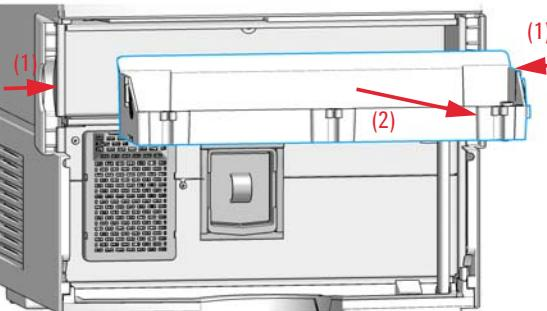
355

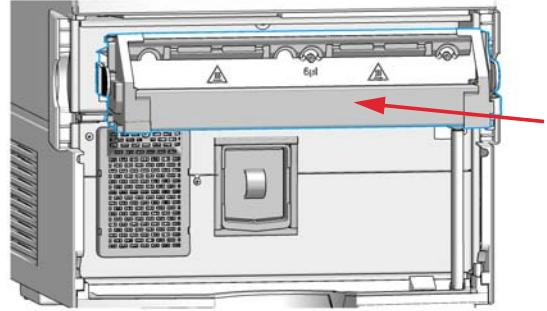
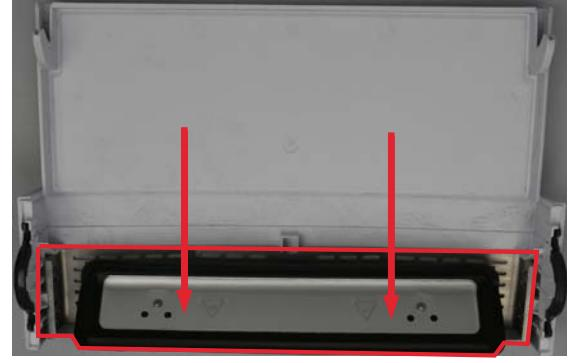
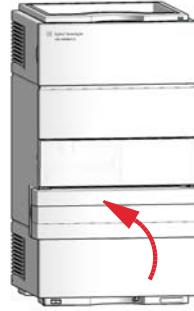
Oven Upgrade

Parts required	p/n	Description
	G4297A	1220 Infinity Oven Kit

Software required Lab Advisor Software

Preparations Switch off the instrument.

1 Remove the lower front cover. 	2 Disconnect the column and remove it ("Remove a Column" on page 55).
3 Remove the leak drain tube.  Leak drain tube	4 Press the knurled portion at either side of the column tray inwards (1) and remove the column tray (2). 

<p>5 Unpack the oven upgrade kit and separate the two parts.</p>	<p>6 Click the oven into position in place of the column tray.</p>  <p>The electrical connection to the oven is made automatically.</p>
<p>7 Replace the leak drain tube.</p>	<p>8 Click the oven insulation into place in the lower front cover, with the cutout in the oven insulation support at the bottom.</p> 
<p>9 Install the column and reconnect the capillaries ("Install a Column" on page 53).</p>	<p>10 Close the lower front cover.</p> 

13 Upgrading the 1220 Infinity II LC

Oven Upgrade

Next Steps:

11 Configure the oven in the Lab Advisor software.

12 Power cycle the instrument and start your Chromatographic Data System.

13 Use Auto configure and the instrument now should be registered under its new main assembly number.

Install the (optional) External Tray

Preparations

Remove all solvent bottles from the solvent bottle cabinet.

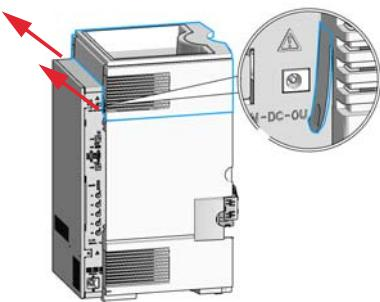
WARNING

Sharp metal edges

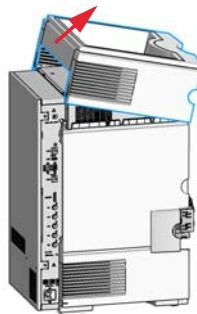
Sharp-edged parts of the equipment may cause injuries.

- To prevent personal injury, be careful when getting in contact with sharp metal areas.

- 1 Lift the two clips at the rear of the instrument to remove the top cover.



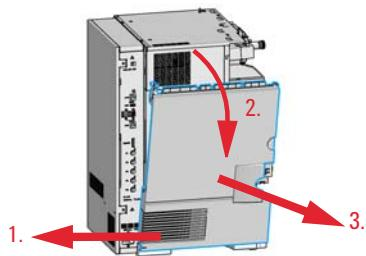
- 2 Remove the top cover.



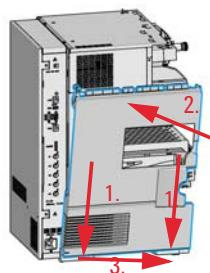
13 Upgrading the 1220 Infinity II LC

Install the (optional) External Tray

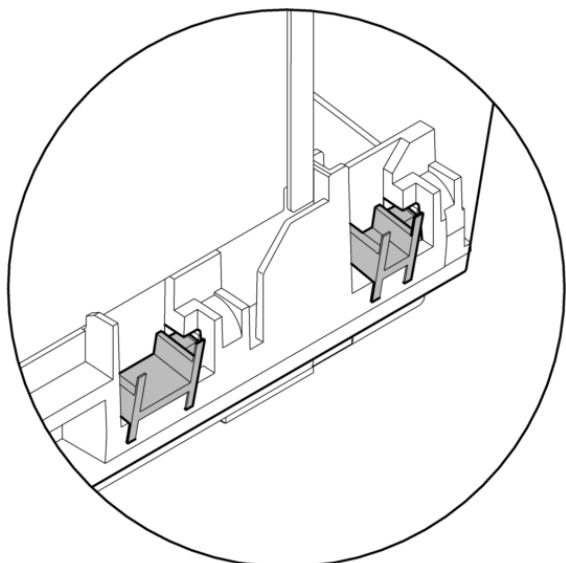
- 3** Remove the left side cover by pushing it carefully towards the rear (1) to get it out of the holders at the bottom plate and the metal frame. Once it is free of the holders, you can swing it towards the side (2) and pull it finally out (3).



- 4** Take the new side cover with the open area, fit it into the holders (see detail view below) at the bottom plate (1) and swing it to the metal cover (2). Push the side cover to the front (3) and verify that the holders on the metal plate catch the side cover correctly.



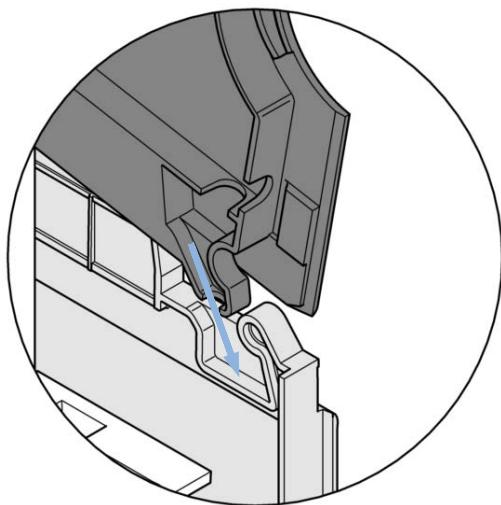
Detail View of the holder mechanism between side cover and bottom plate:



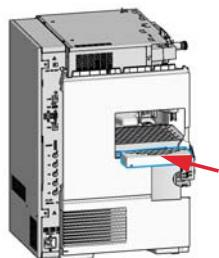
- 5 Hook in the top cover at the front first.

NOTE

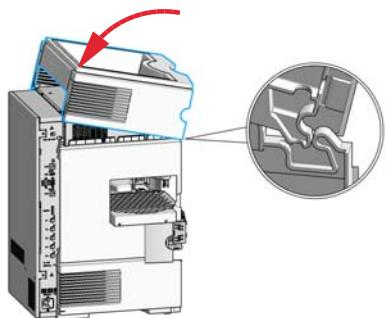
Pay attention to the capillary connected at the pump's outlet ball valve.



- 7 Install the external tray to the module by attaching the fixing rods (attached to it) into the two holes on the left side of the autosampler until you hear a click.



- 6 Carefully swing down the top cover to lock into the rear clips.



- 8 Remove the two rubber plugs that are located at the upper and the lower part of the tray base.

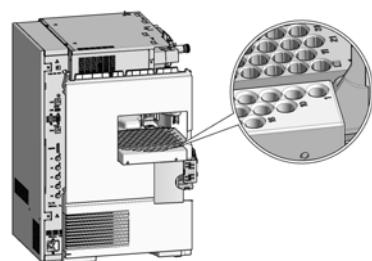


13 Upgrading the 1220 Infinity II LC

Install the (optional) External Tray

9 Install the external tray into the autosampler tray base.

10 Verify the position of the tray and correct it if necessary.
It has to be flush with the inner side of the side cover.



Install the Active Inlet Valve

Tools required**p/n****Description**

Wrench, 14 mm

5067-5688

Torque wrench 1 – 25 Nm with 14 mm wrench

Parts required**p/n****Description**

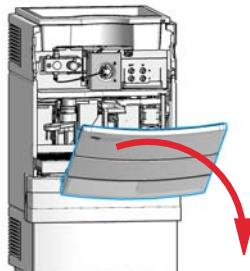
G5699B

1220 Infinity II AIV Upgrade Kit

Preparations

- Switch off the instrument at the main power switch.
- Use an optional solvent shutoff valve or lift up solvent filters in solvent reservoirs to avoid leakages.

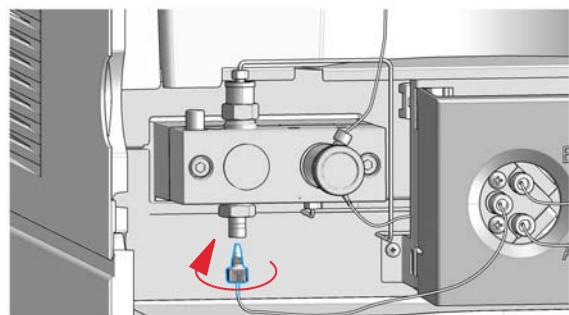
- 1**
- Remove the upper front cover.



- 2**
- Disconnect the solvent inlet tube from the inlet valve.

NOTE

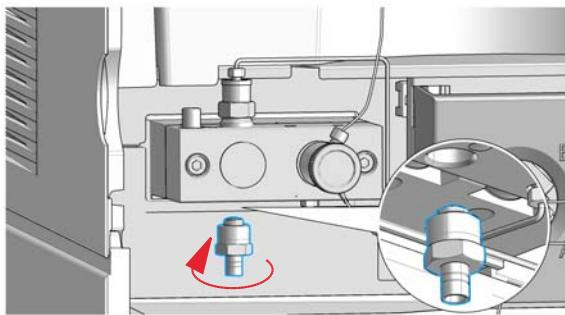
Beware of leaking solvents due to hydrostatic pressure.



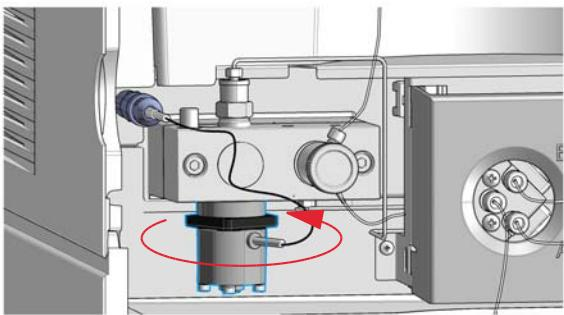
13 Upgrading the 1220 Infinity II LC

Install the Active Inlet Valve

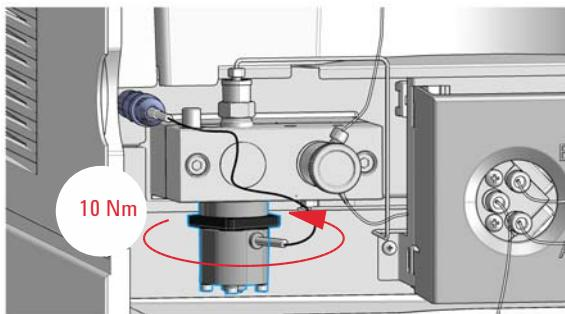
- 3** Using a 14 mm wrench, loosen the passive inlet valve and remove the valve from the pump head.



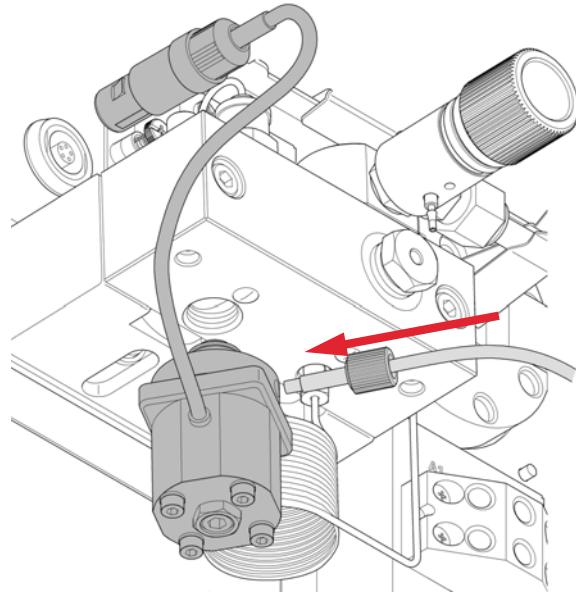
- 4** Insert the active inlet valve into the pump head. Using the 14 mm, turn the nut until it is hand tight. Position the valve such that the solvent inlet tube connection points towards the front.



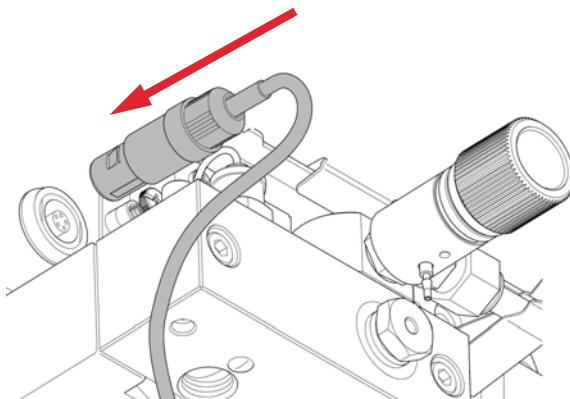
- 5** Tighten the active inlet valve using a 14 mm wrench (approx. 10 Nm).



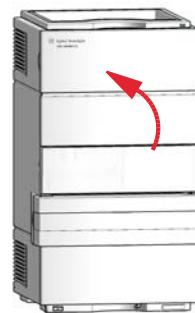
- 6** Use the new solvent tubing provided in update kit to connect the flow path between AIV and DCGV.



- 7** Connect the active inlet valve cable to the connector in the Z-panel.



- 8** Close the upper front cover.



13 Upgrading the 1220 Infinity II LC

Install the Active Inlet Valve

14 Identifying Cables

- Cable Overview [368](#)
- Analog Cables [370](#)
- Remote Cables [372](#)
- BCD Cables [375](#)
- CAN/LAN Cables [377](#)
- Agilent 1200 module to PC [378](#)

This chapter provides information on cables used with the Agilent 1200 Infinity Series modules.



14 Identifying Cables

Cable Overview

Cable Overview

NOTE

Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.

Analog cables

p/n	Description
35900-60750	Agilent module to 3394/6 integrators
35900-60750	Agilent 35900A A/D converter
01046-60105	Analog cable (BNC to general purpose, spade lugs)

Remote cables

p/n	Description
03394-60600	Agilent module to 3396A Series I integrators
	3396 Series II / 3395A integrator, see details in section “ Remote Cables ” on page 372
03396-61010	Agilent module to 3396 Series III / 3395B integrators
5061-3378	Remote Cable
01046-60201	Agilent module to general purpose

BCD cables

p/n	Description
03396-60560	Agilent module to 3396 integrators
G1351-81600	Agilent module to general purpose

CAN cables

p/n	Description
5181-1516	CAN cable, Agilent module to module, 0.5 m
5181-1519	CAN cable, Agilent module to module, 1 m

LAN cables

p/n	Description
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5023-0202	Twisted pair network cable, shielded, 7 m (for point to point connection)

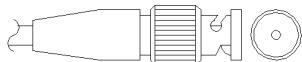
RS-232 cables

p/n	Description
G1530-60600	RS-232 cable, 2 m
RS232-61601	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It's also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3, 3-2, 4-6, 5-5, 6-4, 7-8, 8-7, 9-9.
5181-1561	RS-232 cable, 8 m

14 Identifying Cables

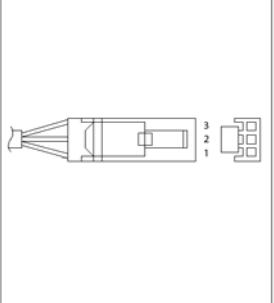
Analog Cables

Analog Cables



One end of these cables provides a BNC connector to be connected to Agilent modules. The other end depends on the instrument to which connection is being made.

Agilent Module to 3394/6 Integrators

p/n 35900-60750	Pin 3394/6	Pin Agilent module	Signal Name
	1		Not connected
	2	Shield	Analog -
	3	Center	Analog +

Agilent Module to BNC Connector

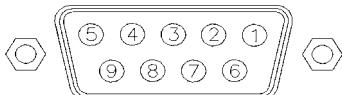
p/n 8120-1840	Pin BNC	Pin Agilent module	Signal Name
	Shield	Shield	Analog -
	Center	Center	Analog +

Agilent Module to General Purpose

14 Identifying Cables

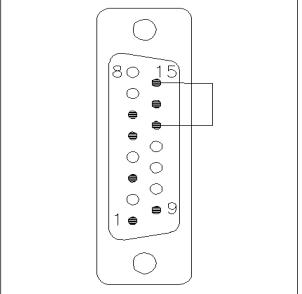
Remote Cables

Remote Cables



One end of these cables provides a Agilent Technologies APG (Analytical Products Group) remote connector to be connected to Agilent modules. The other end depends on the instrument to be connected to.

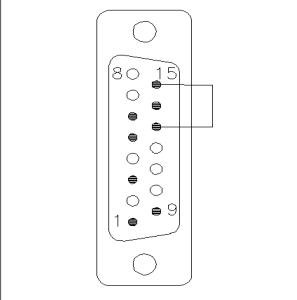
Agilent Module to 3396A Integrators

p/n 03394-60600	Pin 3396A	Pin Agilent module	Signal Name	Active (TTL)
	9	1 - White	Digital ground	
	NC	2 - Brown	Prepare run	Low
	3	3 - Gray	Start	Low
	NC	4 - Blue	Shut down	Low
	NC	5 - Pink	Not connected	
	NC	6 - Yellow	Power on	High
	5,14	7 - Red	Ready	High
	1	8 - Green	Stop	Low
	NC	9 - Black	Start request	Low
	13, 15		Not connected	

Agilent Module to 3396 Series II / 3395A Integrators

Use the cable Agilent module to 3396A Series I integrators (03394-60600) and cut pin #5 on the integrator side. Otherwise the integrator prints START; not ready.

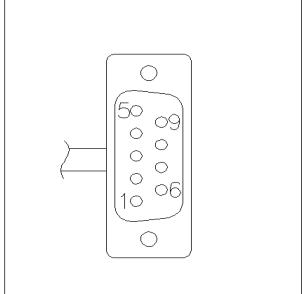
Agilent Module to 3396 Series III / 3395B Integrators

p/n 03396-61010	Pin 33XX	Pin Agilent module	Signal Name	Active (TTL)
	9	1 - White	Digital ground	
	NC	2 - Brown	Prepare run	Low
	3	3 - Gray	Start	Low
	NC	4 - Blue	Shut down	Low
	NC	5 - Pink	Not connected	
	NC	6 - Yellow	Power on	High
	14	7 - Red	Ready	High
	4	8 - Green	Stop	Low
	NC	9 - Black	Start request	Low
	13, 15		Not connected	

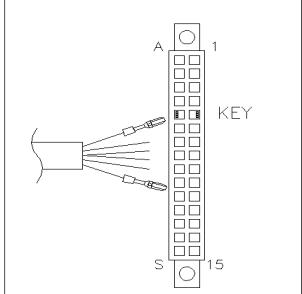
14 Identifying Cables

Remote Cables

Agilent Module to Agilent 35900 A/D Converters

p/n 5061-3378	Pin 35900 A/D	Pin Agilent module	Signal Name	Active (TTL)
	1 - White	1 - White	Digital ground	
	2 - Brown	2 - Brown	Prepare run	Low
	3 - Gray	3 - Gray	Start	Low
	4 - Blue	4 - Blue	Shut down	Low
	5 - Pink	5 - Pink	Not connected	
	6 - Yellow	6 - Yellow	Power on	High
	7 - Red	7 - Red	Ready	High
	8 - Green	8 - Green	Stop	Low
	9 - Black	9 - Black	Start request	Low

Agilent Module to General Purpose

p/n 01046-60201	Wire Color	Pin Agilent module	Signal Name	Active (TTL)
	White	1	Digital ground	
	Brown	2	Prepare run	Low
	Gray	3	Start	Low
	Blue	4	Shut down	Low
	Pink	5	Not connected	
	Yellow	6	Power on	High
	Red	7	Ready	High
	Green	8	Stop	Low
	Black	9	Start request	Low

BCD Cables



One end of these cables provides a 15-pin BCD connector to be connected to the Agilent modules. The other end depends on the instrument to be connected to

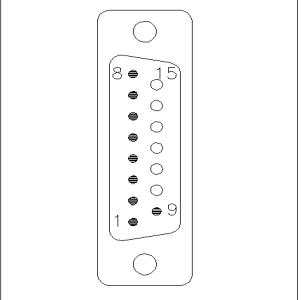
Agilent Module to General Purpose

p/n G1351-81600	Wire Color	Pin Agilent module	Signal Name	BCD Digit
	Green	1	BCD 5	20
	Violet	2	BCD 7	80
	Blue	3	BCD 6	40
	Yellow	4	BCD 4	10
	Black	5	BCD 0	1
	Orange	6	BCD 3	8
	Red	7	BCD 2	4
	Brown	8	BCD 1	2
	Gray	9	Digital ground	Gray
	Gray/pink	10	BCD 11	800
	Red/blue	11	BCD 10	400
	White/green	12	BCD 9	200
	Brown/green	13	BCD 8	100
	not connected	14		
	not connected	15	+ 5 V	Low

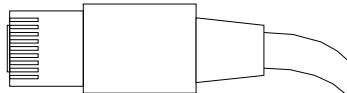
14 Identifying Cables

BCD Cables

Agilent Module to 3396 Integrators

p/n 03396-60560	Pin 3396	Pin Agilent module	Signal Name	BCD Digit
	1	1	BCD 5	20
	2	2	BCD 7	80
	3	3	BCD 6	40
	4	4	BCD 4	10
	5	5	BCD0	1
	6	6	BCD 3	8
	7	7	BCD 2	4
	8	8	BCD 1	2
	9	9	Digital ground	
	NC	15	+ 5 V	Low

CAN/LAN Cables



Both ends of this cable provide a modular plug to be connected to Agilent modules CAN or LAN connectors.

CAN Cables

p/n	Description
5181-1516	CAN cable, Agilent module to module, 0.5 m
5181-1519	CAN cable, Agilent module to module, 1 m

LAN Cables

p/n	Description
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5023-0202	Twisted pair network cable, shielded, 7 m (for point to point connection)

14 Identifying Cables

Agilent 1200 module to PC

Agilent 1200 module to PC

p/n	Description
RS232-61601	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It's also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3, 3-2, 4-6, 5-5, 6-4, 7-8, 8-7, 9-9.
5181-1561	RS-232 cable, 8 m

15 Appendix

- General Safety Information 380
- Solvent Information 383
- Radio Interference 385
- UV Radiation 386
- Sound Emission 387
- Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC) 388
- Declaration of Conformity for HOX2 Filter 389
- Agilent Technologies on Internet 390

This chapter provides addition information on safety, legal and web.



Agilent Technologies

General Safety Information

General Safety Information

The following general safety precautions must be observed during all phases of operation, service, and repair of this instrument. Failure to comply with these precautions or with specific warnings elsewhere in this manual violates safety standards of design, manufacture, and intended use of the instrument. Agilent Technologies assumes no liability for the customer's failure to comply with these requirements.

WARNING

Ensure the proper usage of the equipment.

The protection provided by the equipment may be impaired.

- The operator of this instrument is advised to use the equipment in a manner as specified in this manual.
-

Safety Standards

This is a Safety Class I instrument (provided with terminal for protective earthing) and has been manufactured and tested according to international safety standards.

Operation

Before applying power, comply with the installation section. Additionally the following must be observed.

Do not remove instrument covers when operating. Before the instrument is switched on, all protective earth terminals, extension cords, auto-transformers, and devices connected to it must be connected to a protective earth via a ground socket. Any interruption of the protective earth grounding will cause a potential shock hazard that could result in serious personal injury. Whenever it is likely that the protection has been impaired,

the instrument must be made inoperative and be secured against any intended operation.

Make sure that only fuses with the required rated current and of the specified type (normal blow, time delay, and so on) are used for replacement. The use of repaired fuses and the short-circuiting of fuse holders must be avoided.

Some adjustments described in the manual, are made with power supplied to the instrument, and protective covers removed. Energy available at many points may, if contacted, result in personal injury.

Any adjustment, maintenance, and repair of the opened instrument under voltage should be avoided whenever possible. When inevitable, this has to be carried out by a skilled person who is aware of the hazard involved. Do not attempt internal service or adjustment unless another person, capable of rendering first aid and resuscitation, is present. Do not replace components with power cable connected.

Do not operate the instrument in the presence of flammable gases or fumes. Operation of any electrical instrument in such an environment constitutes a definite safety hazard.

Do not install substitute parts or make any unauthorized modification to the instrument.

Capacitors inside the instrument may still be charged, even though the instrument has been disconnected from its source of supply. Dangerous voltages, capable of causing serious personal injury, are present in this instrument. Use extreme caution when handling, testing and adjusting.

When working with solvents, observe appropriate safety procedures (for example, goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet by the solvent vendor, especially when toxic or hazardous solvents are used.

15 Appendix

General Safety Information

Safety Symbols

Table 48 Safety Symbols

Symbol	Description
	The apparatus is marked with this symbol when the user should refer to the instruction manual in order to protect risk of harm to the operator and to protect the apparatus against damage.
	Indicates dangerous voltages.
	Indicates a protected ground terminal.
	Indicates eye damage may result from directly viewing the light produced by the deuterium lamp used in this product.
	The apparatus is marked with this symbol when hot surfaces are available and the user should not touch it when heated up.

WARNING

A WARNING

alerts you to situations that could cause physical injury or death.

- Do not proceed beyond a warning until you have fully understood and met the indicated conditions.

CAUTION

A CAUTION

alerts you to situations that could cause loss of data, or damage of equipment.

- Do not proceed beyond a caution until you have fully understood and met the indicated conditions.

Solvent Information

Observe the following recommendations on the use of solvents.

Flow Cell

Avoid the use of alkaline solutions ($\text{pH} > 9.5$) which can attack quartz and thus impair the optical properties of the flow cell.

Prevent any crystallization of buffer solutions. This will lead into a blockage/damage of the flow cell.

If the flow cell is transported while temperatures are below $5\text{ }^{\circ}\text{C}$, it must be assured that the cell is filled with alcohol.

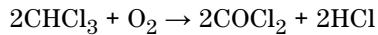
Aqueous solvents in the flow cell can built up algae. Therefore do not leave aqueous solvents sitting in the flow cell. Add small % of organic solvents (for example, acetonitrile or methanol $\sim 5\text{ \%}$).

Solvents

Brown glass ware can avoid growth of algae.

Always filter solvents, small particles can permanently block the capillaries.
Avoid the use of the following steel-corrosive solvents:

- Solutions of alkali halides and their respective acids (for example, lithium iodide, potassium chloride, and so on).
- High concentrations of inorganic acids like nitric acid, sulfuric acid especially at higher temperatures (replace, if your chromatography method allows, by phosphoric acid or phosphate buffer which are less corrosive against stainless steel).
- Halogenated solvents or mixtures which form radicals and/or acids, for example:



This reaction, in which stainless steel probably acts as a catalyst, occurs quickly with dried chloroform if the drying process removes the stabilizing alcohol.

15 Appendix

Solvent Information

- Chromatographic grade ethers, which can contain peroxides (for example, THF, dioxane, di-isopropylether) such ethers should be filtered through dry aluminium oxide which adsorbs the peroxides.
- Solutions of organic acids (acetic acid, formic acid, and so on) in organic solvents. For example, a 1 % solution of acetic acid in methanol will attack steel.
- Solutions containing strong complexing agents (for example, EDTA, ethylene diamine tetra-acetic acid).
- Mixtures of carbon tetrachloride with 2-propanol or THF.

Radio Interference

Cables supplied by Agilent Technologies are screened to provide optimized protection against radio interference. All cables are in compliance with safety or EMC regulations.

Test and Measurement

If test and measurement equipment is operated with unscreened cables, or used for measurements on open set-ups, the user has to assure that under operating conditions the radio interference limits are still met within the premises.

UV Radiation

Emissions of ultraviolet radiation (200 – 315 nm) from this product is limited such that radiant exposure incident upon the unprotected skin or eye of operator or service personnel is limited to the following TLVs (Threshold Limit Values) according to the American Conference of Governmental Industrial Hygienists:

Table 49 UV radiation limits

Exposure/day	Effective irradiance
8 h	0.1 $\mu\text{W}/\text{cm}^2$
10 min	5.0 $\mu\text{W}/\text{cm}^2$

Typically the radiation values are much smaller than these limits:

Table 50 UV radiation typical values

Position	Effective irradiance
Lamp installed, 50 cm distance	average $0.016 \mu\text{W}/\text{cm}^2$
Lamp installed, 50 cm distance	maximum $0.14 \mu\text{W}/\text{cm}^2$

Sound Emission

Manufacturer's Declaration

This statement is provided to comply with the requirements of the German Sound Emission Directive of 18 January 1991.

This product has a sound pressure emission (at the operator position) < 70 dB.

- Sound Pressure $L_p < 70$ dB (A)
- At Operator Position
- Normal Operation
- According to ISO 7779:1988/EN 27779/1991 (Type Test)

15 Appendix

Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC)

Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC)

Abstract

The Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC), adopted by EU Commission on 13 February 2003, is introducing producer responsibility on all electric and electronic appliances starting with 13 August 2005.

NOTE

This product complies with the WEEE Directive (2002/96/EC) marking requirements. The affixed label indicates that you must not discard this electrical/electronic product in domestic household waste.

Product Category:

With reference to the equipment types in the WEEE Directive Annex I, this product is classed as a Monitoring and Control Instrumentation product.



NOTE

Do not dispose of in domestic household waste

To return unwanted products, contact your local Agilent office, or see <http://www.agilent.com> for more information.

Declaration of Conformity for HOX2 Filter

Declaration of Conformity					
<p>We herewith inform you that the</p> <p style="text-align: center;">Holmium Oxide Glass Filter</p> <p>used in Agilent's absorbance detectors listed in the table below meets the requirements of National Institute of Standards and Technology (NIST) to be applied as certified wavelength standard.</p> <p>According to the publication of NIST in J. Res. Natl. Inst. Stand. Technol. 112, 303-306 (2007) the holmium oxide glass filters are inherently stable with respect to the wavelength scale and need no recertification. The expanded uncertainty of the certified wavelength values is 0.2 nm.</p> <p>Agilent Technologies guarantees, as required by NIST, that the material of the filters is holmium oxide glass representing the inherently existent holmium oxide absorption bands.</p> <p>Test wavelengths:</p> <p>Where "x" can be any alphanumeric character</p>					
Product Number	Series	Measured Wavelength *	Wavelength Accuracy	Optical Bandwidth	
G1315x, G1365x	1100, 1200, 1260	361.0 nm 418.9 nm 453.7 nm 536.7 nm	+/- 1 nm	2 nm	
G7115x, G7165x	1260				
G1600x, G7100x	CE				
G1314x	1100, 1200, 1260, 1290	360.8nm 418.5nm 536.4nm	+/- 1 nm	6 nm	
G7114x	1260, 1290				
G4286x,..., 94x	1120, 1220				
<small>*) The variation in Measured Wavelength depends on the different Optical Bandwidth.</small>					
<small>28-Oct-2014</small> <small>(Date)</small>					
 <small>(R&D Manager)</small>		 <small>(Quality Manager)</small>			
P/N 89550-90501 	Revision: G Effective by: 28-Oct-2014		<input checked="" type="checkbox"/>		

15 Appendix

Agilent Technologies on Internet

Agilent Technologies on Internet

For the latest information on products and services visit our worldwide web site on the Internet at:

<http://www.agilent.com>

Select Products/Chemical Analysis

Index

GLP features 20

A

absorbance
Beer-Lambert 137

achromat
source lens 121

active inlet valve
installing 363

adapter 251

Agilent Technologies 390

algae growth 97

algea information 383

alignment teaching 156

alignment
gripper 159

alpha and beta line 188

ambient non-operating temperature 19

ambient operating temperature 19

analog

cable 370

analytical head 108

arm down 157

arm up 157

array 122

ASTM

environmental conditions 18
noise test (ChemStation only) 184

AUTO mode 95

automatic configuration with Bootp 70

autosampler

EMF counters 12
introduction 103

simple repairs 271
transprot assembly parts 272

B

ball-screw drive 88

band width 6.5 nm 25

BCD

cable 375

Beer-Lambert (law) 137

bench space 17

beta and alpha line 188

blank scan 170

blockage 151, 206

BootP service

installation 71

restart 78

settings 77

stop 77

Bootp

automatic configuration 70

storing the settings permanently 79

buffer application 96

buffer solution 264

bypass 105

C

cable

analog 370

BCD 375

CAN 377

LAN 377

remote 372

RS-232 378

cables

analog 368

BCD 368

CAN 369

LAN 369

overview 368

remote 368

RS-232 369

calibration

oven 161

CAN

cable 377

caps 111

cell 162

change

gripper 154

needle 154

piston 154

cleaning the autosampler 272

cleaning the module 303

close gripper 157

column oven 116

column

installing 53

removing 55

communications 20

compensation sensor open 197

compensation sensor short 197

composition precision 21, 22

composition range 21, 22

compressibility compensation 21, 22, 93

condensation 18

configurations 10

control

system 20

Index

correction factors for flow cells 138

counter

 autosampler 12

 detector 13

 injection valve 13

 liter 12

 needle movement 13

 seal wear 12

cuvette holder 299

D

D/A converter 191

DAC

 Agilent LabAdvisor 191

 Instant Pilot 192

dark current 163

 test 175

data evaluation 20

declaration of conformity 389

degasser exchange 153

delay volume 21, 91, 91

delivery checklist 30

description 100

detection type 25

detection

 compound classes 132

detector

 EMF counters 13

 features 119

DHCP

 general information 66

 setup 67

diagnostic buffers 144

diagnostic tables 144

dimensions 19

diode

 array 122, 122

 width 26

draw 156

drift (ASTM) and noise 26

drift 25

DSP not running 239

dual plunger in-series design 86

dual-channel gradient valve 264

E

Early maintenance feedback 20

EE 2060 209

electronic waste 388

electrostatic discharge (ESD) 245, 247

EMF

 pump head 12

EMF

 counters, pump 12

encoder missing 203

entrance slit 122

environment 18

error message

 ADC hardware error 229

error messages

 arm movement failed 216

 calibration failed 229

 compensation sensor open 197

 compensation sensor short 197

 diode current leakage 236

 DSP not running 239

 encoder missing 203

 fan failed 198

 filter check failed 230

 filter missing 230

 grating missing 232

 grating/filter motor defective 231

 heater current missing 232

 heater failed 226

 heater power at limit 226

 holmium oxide test failed 233, 238

 illegal temperature value from sensor at air inlet 227

illegal value from sensor on main board 227

index adjustment 203

index limit 204

index missing 204

initialization failed 205, 217

initialization with vial 217

invalid vial position 218

lamp ignition failed 233

leak sensor open 199

leak sensor short 200

leak 199

lost CAN partner 201

MCGV fuse 213

metering home failed 219

missing pressure reading 205

missing vial 219

missing wash vial 220

motor failed 220

motor-drive power 206

needle down failed 221

needle up failed 222

no run data available in device 239

pressure above upper limit 207

pressure below lower limit 207

pressure signal missing 208

pump configuration 209

pump head missing 209

remote timeout 200

safety flap missing 223

servo restart failed 210

shutdown 201

solvent zero counter 214

stroke length 211

temperature limit exceeded 211

temperature out of range 212

timeout 202

uv heater current 237

uv ignition failed 236

uv lamp current 227

uv lamp voltage 228

- valve failed 212
 valve to bypass failed 223
 valve to mainpass failed 224
 vial in gripper 224
 visible lamp current 235
 visible lamp voltage 235
 wait timeout 213
 wavelength calibration failed 238
 wavelength check failed 234
 wavelength recalibration lost 239
 exchange degasser 153
 exchanging
 dual channel gradient valve (DCGV) 264
 injection seal 268
 passive inlet valve 250
 purge valve frit 254
 purge valve 254
 export data 170
- F**
- fan failed 198
 features
 GLP 25
 safety and maintenance 25
 filter motor 167
 filter test 172
 filter/grating test 167
 firmware 144
 updates 325, 325
 upgrade/downgrade 325
 upgrade/downgrade 325
 flow cell
 correction factors 138
 specifications 26
 standard (parts) 348
 support windows 121
 test 184
 types and data 25
 flow path blockage 151
- flow precision 21, 21, 22, 22
 flow range 21, 22
 flow
 unstable 323
 frequency range 19
- G**
- general error messages 197
 gradient formation 21, 22
 gradient valve (DCGV) 264
 grating motor 167
 grating 122, 122
 gripper arm
 repair 288
 gripper fingers 109
 gripper verification 159
 gripper
 change 154
- H**
- half trays 110
 holmium oxide
 declaration of conformity 389
 filter 121
 test 180
 humidity 19
 hydraulic path 86
 hydraulic system 21, 22
- I**
- increased system pressure 324
 index adjustment 203
 index limit 204
 index missing 204
 information 144
 on cuvette holder 299
 on solvents 383
 on UV radiation 386
- initialization failed 205
 initialization mode selection 64
 initialization
 pump 89
 injecting sample 100
 INJECT 100, 102
 injection seal 268
 tefzel 100
 vespel 100
 injection sequence 105
 injection valve 103, 107, 108
 inlet valve 250
 installation check 143
 installation
 delivery checklist 30
 site requirements 16
 installing the autosampler
 sample trays 110
 installing
 active inlet valve 363
 column 53
 oven upgrade kit 356
 restriction capillary 57
 intensity test 165, 177
 Internet 390
 introduction
 optical unit parts 121
- L**
- Lab Advisor 243
 lamp intensity 165
 lamp
 type 25
 lamps 121
 LAN
 automatic configuration with Bootp 70
 cable 377
 initialization mode selection 64

Index

- link configuration selection 69
 - manual configuration with telnet 81
 - manual configuration 80
 - storing the settings permanently 79
 - TCP/IP parameter configuration 62
 - using default 65
 - using stored 64
 - leak sensor open 199
 - leak sensor short 200
 - leak test 145
 - leakage current 163
 - leak
 - correcting 321
 - line frequency 19
 - line voltage 19
 - linear range 26
 - linearity 25
 - link configuration selection 69
 - liquimeter 12
 - LOAD 100, 101
 - lost CAN partner 201
- ## M
- MAC address
 - determine 74
 - mainpass 105
 - maintenance functions 273
 - step commands 155
 - maintenance
 - exchanging lamps 291
 - overview 302
 - replacing firmware 325, 325
 - using the cuvette holder 299
 - make-before-break 100
 - manual configuration
 - of LAN 80
 - manual control 155
 - materials in contact with mobile phase 91, 91
- MCGV fuse 213
 - message
 - ADC hardware error 229
 - calibration failed 229
 - calibration lost 234
 - diode current leakage 236
 - filter check failed 230
 - filter missing 230
 - grating missing 232
 - grating/filter motor defective 231
 - heater current missing 232
 - heater failed 226
 - heater power at limit 226
 - holmium oxide test failed 233, 238
 - illegal temperature value from sensor at air inlet 227
 - illegal value from sensor on main board 227
 - lamp ignition failed 233
 - no run data available in device 239
 - remote timeout 200
 - uv heater current 237
 - uv ignition failed 236
 - uv lamp current 227
 - uv lamp voltage 228
 - visible lamp current 235
 - visible lamp voltage 235
 - wavelength calibration failed 238
 - wavelength check failed 234
 - metering device 107
 - missing pressure reading 205
 - module info 144
 - module options 144
 - motor-drive power 206
 - move arm home 155
 - multi-draw option 103
- ## N
- needle drive 107, 107
 - needle into sample 156
- needle into seat 156
 - needle type 102
 - needle up 155, 156
 - needle
 - change 154
 - needles 102
 - negative absorbance 130
 - noise and drift (ASTM) 26
 - noise, short term 25
 - non-operating altitude 19
 - non-operating temperature 19
 - numbering of vials 110
- ## O
- open gripper 157
 - operating Altitude 19
 - operating temperature 19
 - operational pressure range 21, 22
 - optimization
 - margins for negative absorbance 130
 - of selectivity 130
 - peak width 123
 - sample and reference wavelength 124
 - slit width 127
 - spectra acquisition 129
 - oven calibration 161
 - oven test 160
 - oven upgrade kit
 - installing 356
- ## P
- park arm 155
 - parts for maintenance
 - standard flow cell 348
 - parts identification
 - cables 367
 - parts

- repair 348
peak width (response time) 123
pH range 21, 22
photometric accuracy 138
physical specifications 19, 19
piston chamber 86
piston
 change 154
plateaus, leak test 149
plunger home 155
PM 243
power consumption 19
power
 considerations 16
 cords 17
precision 101
pressure above upper limit 207
pressure below lower limit 207
pressure profile 145
pressure pulsation 21, 22, 93, 95
pressure signal missing 208
pressure too high check 151
pressure, operating range 21, 22
pressure 86
preventive maintenance 243
programmable slit width 26
proportioning valve
 high-speed 86
PTFE frit 254
pump configuration 209
pump failure 323
pump head missing 209
pump head
 reinstalling 262
pump piston 96
pump
 overview 86
pump
 functional principle 88
hints for successful use 96
purge pump 152
purge valve frit 96
purge valve 254
cable 378
run recovery
 no run data available in 239
running the leak test 147
- ## R
- radio interference 385
reassembling the pump head 262
recommended pH range 21, 22
remote
 cable 372
removing
 column 55
 restriction capillary 51
repair procedures
 injection seal 268
repair
 parts 348
repairs
 cleaning the instrument 303
 correction leaks 321
 metering plunger 284
 metering seal 284
 needle assembly 274
 needle-seat assembly 278
 overview of simple repairs 290
 replacing firmware 325, 325
 replacing leak handling system 322
 rotor seal 280
reset 156
response time (peak width) 123
restart without cover 209
restriction capillary
 installing 57
 removing 51
RFID tag 120
rotor seal
 exchange 280
RS-232C
- ## S
- safety class I 380
safety features
 system 20
safety
 general information 380
 standards 19
 symbols 382
sample and reference wavelength 124
sample loops 100
sample scan 170
sample trays 110
 numbering of vial positions 110
sample volume 101
sampling sequence 104
sampling unit 107
sapphire piston 88, 88
scan 170
seal wear counters 12
selectivity optimization 130
servo restart failed 210
setable flow range 21, 22
shutdown 201
signal plots 144
simple repairs
 autosampler 271
site requirements
 bench space 17
 environment 18
 power considerations 16
 power cords 17
slit test 174
slit width 26, 127
solvent delivery system 86

Index

- solvent filters
 checking 248
 cleaning 249
 prevent blockage 97
- solvent information 383
- solvent inlet filters 96
- Solvent Selection Valve 10, 11
- solvent zero counter 214
- specification
 physical 19
- specifications
 diode width 26
 flow cell 26
 linear range 26
 noise and drift (ASTM) 26
 physical 19
 programmable slit width 26
 wavelength accuracy 26
 wavelength range 26
- spectral flatness test 182
- spectral scan 170
- spectra
 acquisition 129
- spectrograph
 diodes per nm 122
- SSV 10, 11
- state info 144
- stator 108
- step commands 155
- stepper motor 107
- storing the settings permanently 79
- stroke length 211
- stroke volume 88, 95
- suppression
 quantifying 130
- system configurations 10
- T**
- TCP/IP parameter configuration 62
- telnet
 configuration 81
- temperature limit exceeded 211
- temperature out of range 212
- temperature sensor 199
- test chromatogram 169, 186
- test function
 D/A converter 191
 DAC 191
- tests
 ASTM noise (ChemStation only) 184
 dark current 175
 filter 172
 flow cell (ChemStation only) 184
 holmium oxide 180
 intensity 177
 slit 174
 spectral flatness (ChemStation only) 182
 test chromatogram 186
 wavelength calibration 168
- theta-axis 109
- timeout 202
- transport assembly 109
- transport mechanism 103
- troubleshooting
 error messages 196
- U**
- URL 390
- using
 the cuvette holder 299
- UV lamp on 170
- UV radiation 386
- V**
- vacuum degasser 96
- valve bypass 155
- valve failed 212
- valve frit 254
- valve mainpass 156
- valve
 proportioning 86
- variable entrance slit 122
- variable reluctance motor 88
- variable stroke volume 95
- verification and recalibration of wavelength 188
- vial numbering 110
- vial racks 103
- vial to seat 155
- vial to tray 156
- vials 103, 111
- voltage range 19
- VWD
 EMF counters 13
- W**
- wait timeout 213
- waste
 electrical and electronic equipment 388
- wavelength recalibration lost 239
- wavelength
 accuracy 26, 26
 calibration 168
 range 190-600 nm 25
 range 26
 verification and recalibration 188
- WEEE directive 388
- weight 19
- X**
- X-axis 109
- Z**
- Z-axis 109

In This Book

This manual contains information on how to use, maintain, and upgrade the Agilent 1220 Infinity II LC System.

The manual contains the following chapters:

- Introduction
- Installation
- Agilent 1220 Infinity II LC Description
- Test Functions and Calibration
- Error Information
- Maintenance
- Parts for Maintenance
- Upgrading the Agilent 1220 Infinity II LC
- Appendix

© Agilent Technologies 2016

Printed in Germany
09/2016



G4280-90020 Rev. B



Agilent Technologies