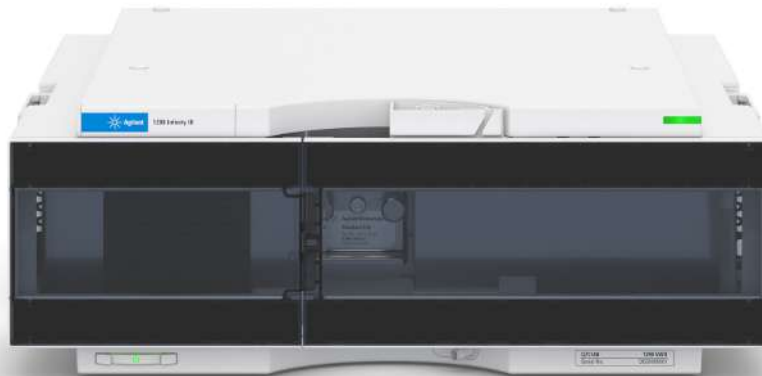




Agilent InfinityLab LC Series
Variable Wavelength Detectors
User Manual



Notices

Document Information

The information in this document also applies to 1260 Infinity II and 1290 Infinity II modules.

Document No: SD-29000240 Rev. E
Edition: 10/2024

Copyright

© Agilent Technologies, Inc.
2014-2024

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.

Agilent Technologies
Hewlett-Packard-Strasse 8
76337 Waldbronn, Germany

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

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

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.

Contents

1 Introduction 8

Overview of the Module 9

Product Description of the 1260 Infinity III Variable Wavelength Detector (G7114A) 10

Features of the 1260 Infinity III Variable Wavelength Detector (G7114A) 11

Product Description of the 1290 Infinity III Variable Wavelength Detector (G7114B) 12

Features of the 1290 Infinity III Variable Wavelength Detector (G7114B) 13

Operating Principle 14

Dual-Wavelength Mode 20

2 Site Requirements and Specifications 21

Site Requirements 22

Specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A) 25

Specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B) 28

Specification Conditions 31

3 Installation 32

Installing Capillaries 33

Handling Leak and Waste 38

Connecting Modules and Control Software 48

Instrument Configuration 49

4 Using the Module 51

General Information 52

Preparation of the System 56

Preparing the Module 66

5	Optimizing the Performance of the Module	75
	Introduction	76
	Match the Flow Cell to the Column	77
	Set the Detector Parameters	81
	Warm up of the Detector	82
6	Diagnostics and Troubleshooting	84
	Diagnostic Features	85
	Overview of Available Tests and Tools	86
	Maintenance and Troubleshooting Tools of the Module	87
	Agilent Lab Advisor Software	111
	Other Lab Advisor Functions	112
7	Error Information	113
	What Are Error Messages	115
	General Error Messages	116
	Detector Error Messages	129
8	Maintenance	144
	Introduction to Maintenance	145
	Safety Information Related to Maintenance	146
	Overview of Maintenance	148
	Cleaning the Module	149
	Remove and Install Doors	150
	Replace the Deuterium Lamp	153
	Replace the Flow Cell / Cuvette Holder	158
	Repairing the Flow Cells	162
	Using the Cuvette Holder	165
	Correcting Leaks	168
	Replace Leak Handling System Parts	170
	Replace the Module Firmware	171
9	Parts and Materials for Maintenance	173
	Overview of Maintenance Parts	174
	Standard Flow Cell 10 mm / 14 μ L	176

Micro Flow Cell 3 mm / 2 μ L 178
Semi-micro Flow Cell (Parts) 180
High Pressure Flow Cell (Parts) 182
Bio Standard Flow Cell 183
Bio Micro Flow Cell 185
Cuvette Holder (Parts) 187
Accessory Kit 188

10 Identifying Cables 189

Cable Overview 190
Analog Cables 192
Remote Cables 194
BCD Cables 198
CAN/LAN Cables 200
RS-232 Cables 201
USB 202

11 Hardware Information 203

General Hardware Information 204
Module-Specific Hardware Information 218

12 LAN Configuration 222

What You Have to Do First 223
TCP/IP Parameter Configuration 224
Configuration Switch 225
Initialization Mode Selection 226
Dynamic Host Configuration Protocol (DHCP) 228
Manual Configuration 231
PC and User Interface Software Setup 236

13 Appendix 239

General Safety Information 240
Material Information 248
At-a-Glance Details About Agilent Capillaries 254
Waste Electrical and Electronic Equipment (WEEE) Directive 258

Radio Interference	259
RFID Statement	260
Sound Emission	262
Declaration of Conformity for HOX2 Filter	263
Agilent Technologies on Internet	265



In This Book

This manual covers the following Agilent InfinityLab LC Series modules:

- Agilent 1260 Infinity III Variable Wavelength Detector (G7114A)
- Agilent 1290 Infinity III Variable Wavelength Detector (G7114B)

Find information on other Agilent Variable Wavelength Detectors in separate manuals.

This chapter gives an introduction to the module and instrument overview.

Overview of the Module 9

Product Description of the 1260 Infinity III Variable Wavelength Detector (G7114A) 10

Features of the 1260 Infinity III Variable Wavelength Detector (G7114A) 11

Product Description of the 1290 Infinity III Variable Wavelength Detector (G7114B) 12

Features of the 1290 Infinity III Variable Wavelength Detector (G7114B) 13

Operating Principle 14

Hydraulic Path 15

Flow Cell 15

Lamp 16

Source Lens Assembly 17

Entrance Slit Assembly 17

Filter Assembly 17

Mirror Assemblies M1 and M2 18

Grating Assembly 18

Beam Splitter Assembly 18

Photo Diodes Assemblies 19

Photo Diode ADC (Analog-to-Digital Converter) 19

Dual-Wavelength Mode 20

Overview of the Module

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

- Higher data rate up to 120 Hz (G7114A) or 240 Hz for ultra-fast-HPLC (G7114B)
- Deuterium lamp for highest intensity and lowest detection limit over a wavelength range of 190 to 600 nm
- Optional flow-cell cartridges (standard 10 mm, 14 μ L; high pressure 10 mm, 14 μ L; micro 3 mm, 2 μ L; semi-micro 6 mm, 5 μ L), Bio, and Prep Cells are available and can be used depending on the application needs (other types may be introduced later)
- Dual wavelength mode, see [Dual-Wavelength Mode](#) on page 20
- Easy front access to lamp and flow cell for fast replacement
- Electronic identification of flow cell and lamp with RFID (Radio Frequency Identification) tag for unambiguous identification
 - Lamp information: part number, serial number, production date, ignitions, burn time
 - Cell information: part number, serial number, production date, nominal path length, volume, maximum pressure
- Built-in electronic temperature control (ETC) for improved baseline stability
- Built-in holmium oxide filter for fast wavelength accuracy verification

Introduction

Product Description of the 1260 Infinity III Variable Wavelength Detector (G7114A)

Product Description of the 1260 Infinity III Variable Wavelength Detector (G7114A)

The Agilent 1260 Infinity III Variable Wavelength Detector (VWD) is the most sensitive and fastest detector in its class.

Time-programmable wavelength switching provides sensitivity and selectivity for your applications.

More sample information can be acquired in the dual wavelength mode.

Low detector noise ($<\pm 2.5 \mu\text{AU}$) and baseline drift ($<1 \cdot 10^{-4} \text{ AU/h}$) facilitates precise quantification of trace levels components.

High productivity can be achieved with fast analysis at up to 120 Hz data rates.

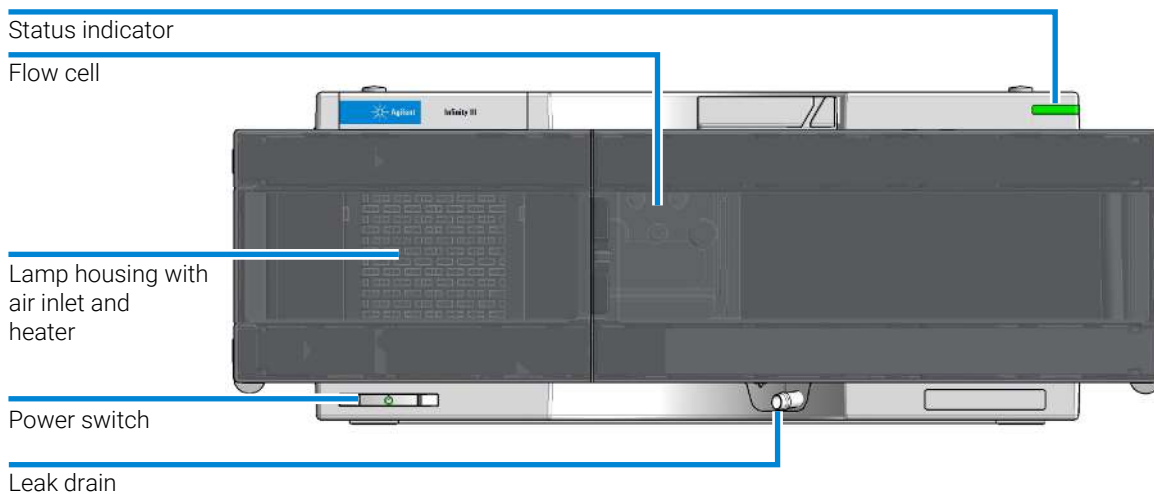


Figure 1: Overview of the G7114A Detector

Features of the 1260 Infinity III Variable Wavelength Detector (G7114A)

- High sensitivity with lowest baseline drift achieved by significant hardware improvements over the most successful predecessor Agilent 1200 Series VWD.
- Electronic temperature control (ETC) - maximum baseline stability and practical sensitivity under fluctuating ambient temperature and humidity conditions.
- Up to 100 % resolution gain in fast LC by 120 Hz data acquisition rate.
- Wide linear range - for reliable, simultaneous quantification of primary compounds, by-products and impurities.
- Stop-flow wavelength scanning for quick wavelength optimization.
- Radio frequency identification (RFID) technology for flow cells and lamps - a new level of data traceability.
- Automatic wavelength verification by built-in holmium oxide filter.
- Extensive diagnostics, error detection and display with Instant Pilot controller and Agilent Lab Advisor software.
- Early maintenance feedback (EMF) for continuous tracking of instrument usage in terms of lamp burn-time with user-settable limits and feedback messages.
- Front access to lamps and flow cells enables fast and convenient maintenance for maximum uptime.

Introduction

Product Description of the 1290 Infinity III Variable Wavelength Detector (G7114B)

Product Description of the 1290 Infinity III Variable Wavelength Detector (G7114B)

The Agilent 1290 Infinity III Variable Wavelength Detector (VWD) is the most sensitive and fastest detector in its class.

Time-programmable wavelength switching provides sensitivity and selectivity for your applications.

More sample information can be acquired in the dual wavelength mode.

Low detector noise ($< \pm 1.5 \mu\text{AU}$) and baseline drift ($< 1 \cdot 10^{-4} \text{ AU/h}$) facilitates precise quantification of trace levels components.

High productivity can be achieved with fast analysis at up to 240 Hz data rates.

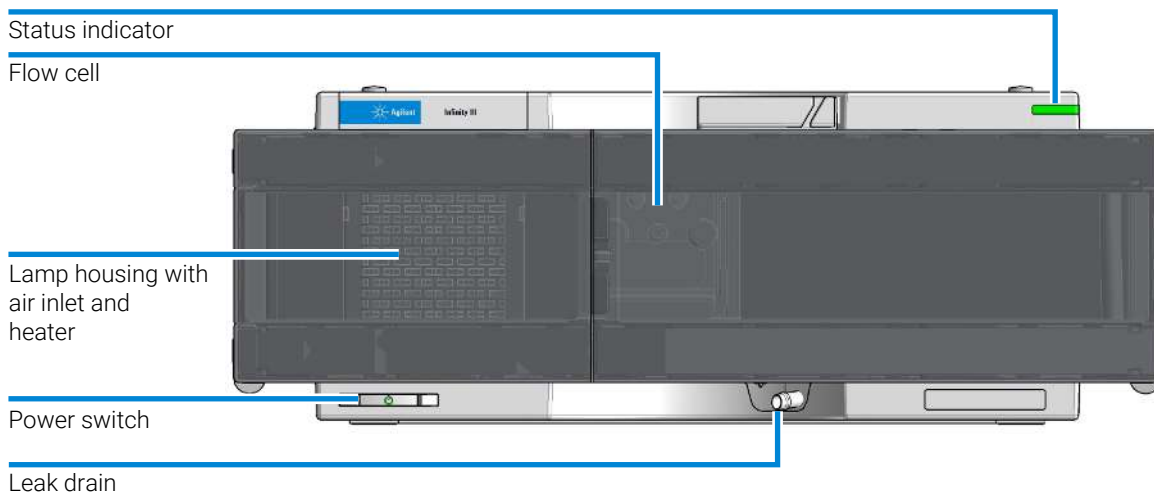


Figure 2: Overview of the G7114B Detector

Features of the 1290 Infinity III Variable Wavelength Detector (G7114B)

- Low baseline noise and drift results in low detection limits for robust quantification of trace level components.
- Dual-wavelength capabilities offer more analyte information per run.
- Up to 100 % resolution gain in fast LC, at up to 240 Hz data acquisition rate.
- Wide linear range (>2.5 AU upper limit) enables simultaneous quantification of primary compounds, by-products and impurities.
- Electronic temperature control (ETC) – for high baseline stability and practical sensitivity under fluctuating ambient temperature and humidity conditions.
- Automatic wavelength verification provided through the use of built-in holmium oxide filter.
- Radio frequency identification (RFID) technology for flow cells and lamps improve data traceability.
- Early maintenance feedback (EMF) for continuous tracking of instrument usage, in terms of lamp burn time, with user-defined limits and message types.
- Extensive diagnostics, error detection and displays provided using Agilent Lab Advisor software.
- Stop-flow wavelength scanning for fast wavelength optimization.

Operating Principle

The optical system of the detector is shown in the figure below. Its radiation source is a deuterium-arc discharge lamp for the ultraviolet (UV) wavelength range from 190 to 600 nm. The light beam from the deuterium lamp passes through a lens, a filter assembly, an entrance slit, a spherical mirror (M1), a grating, a second spherical mirror (M2), a beam splitter, and finally through a flow cell to the sample diode. The beam through the flow cell is absorbed depending on the solutions in the cell, in which UV absorption takes place, and the intensity is converted to an electrical signal by means of the sample photodiode. Part of the light is directed to the reference photodiode by the beam splitter to obtain a reference signal for compensation of intensity fluctuation of the light source. A slit in front of the reference photodiode cuts out light of the sample bandwidth. Wavelength selection is made by rotating the grating, which is driven directly by a stepper motor. This configuration allows fast change of the wavelength. The cutoff filter is moved into the lightpath above 370 nm to reduce higher order light.

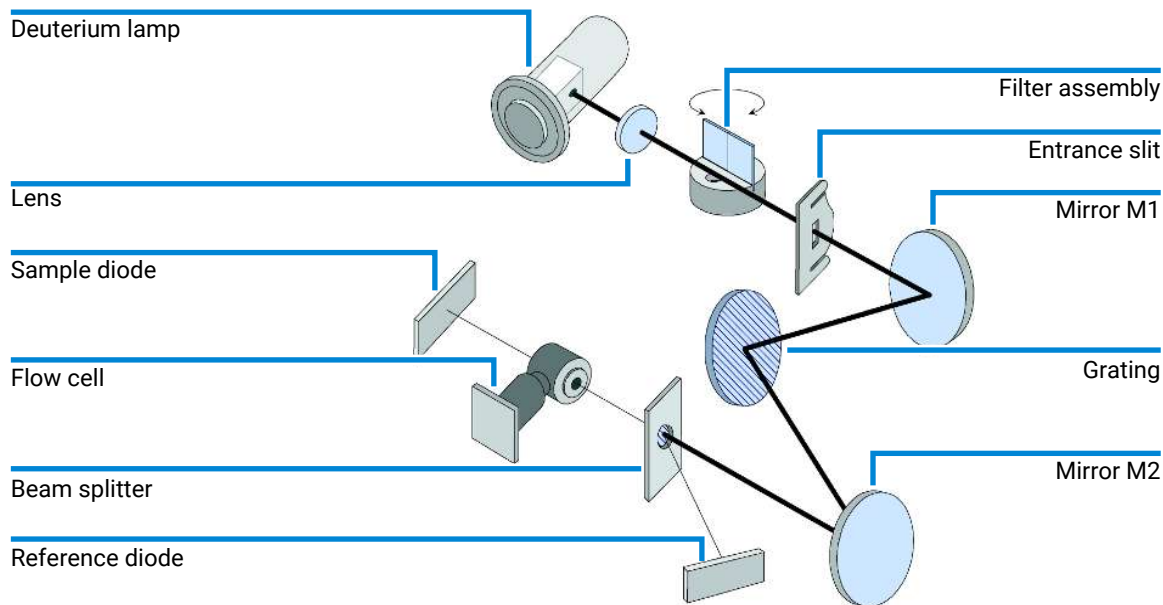


Figure 3: Optical Path of the Variable Wavelength Detector

Hydraulic Path

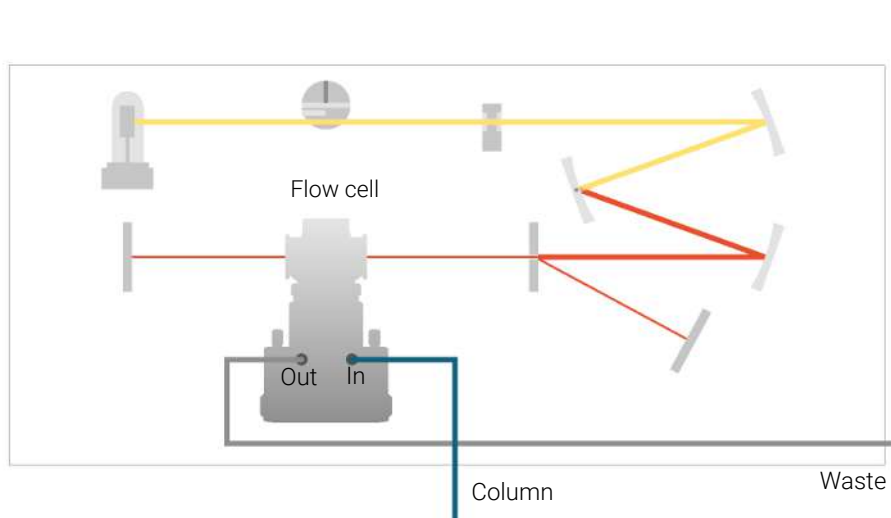


Figure 4: Hydraulic path

Flow Cell

A variety of flow-cell cartridges can be inserted using the same quick and simple mounting system.

The flow cells have an integrated RFID tag that contains the flow cell specific information (e.g. part number, cell volume, path length, ...). A RFID tag reader reads out this information and transfers it to the user interface.

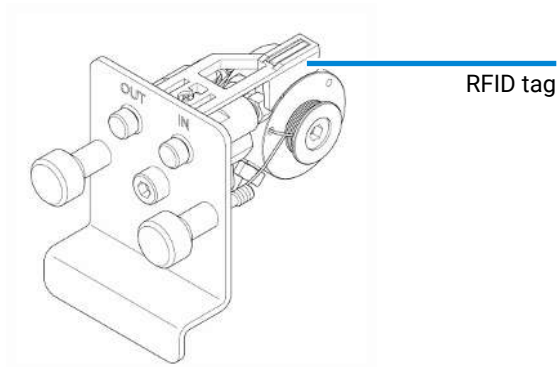


Figure 5: Flow Cell with RFID tag

Table 1: Flow cell data

	Standard	Semi-micro	Micro	High Pressure	Bio Standard	Bio Micro
Maximum pressure (bar (MPa))	40 (4)	40 (4)	120 (12)	400 (40)		
Path length (mm)	10 (conical)	6 (conical)	3 (conical)	10 (conical)	10 (conical)	3 (conical)
Volume (μL)	14	5	2	14	14	2
Inlet i.d. (mm)	0.25	0.17	0.12	0.25	14	0.12
Inlet length (mm)	750	250	310	750	750	310
Outlet i.d. (mm)	0.30	0.17	0.17	0.17	0.30	310
Outlet length (mm)	120	120	120	120	120	310
Total volume (μL)	60.77	14.49	14.00	60.77	60.77	310
Materials in contact with solvent	SST, quartz, PTFE, PEEK	SST, quartz, PTFE	SST, quartz, PTFE	SST, quartz, Kapton	MP35N, sapphire	MP35N, sapphire

Lamp

The light source for the UV wavelength range is a deuterium lamp. As a result of plasma discharge in a low pressure deuterium gas, the lamp emits light over the 190 – 600 nm wavelength range.

Introduction

Operating Principle

The lamp has an integrated RFID tag that contains the lamp specific information (e.g. part number, burn time, ...). An RFID tag reader reads out this information and transfers it to the user interface.

Source Lens Assembly

The source lens receives the light from the deuterium lamp and focuses it onto the entrance slit.

Entrance Slit Assembly

The entrance slit assembly has an exchangeable slit. The standard one has a 1-mm slit. For replacement and calibration purposes to optimize the alignment, a slit with a hole is needed.

Filter Assembly

The filter assembly is electromechanically actuated. During wavelength calibrations it moves into the light path.

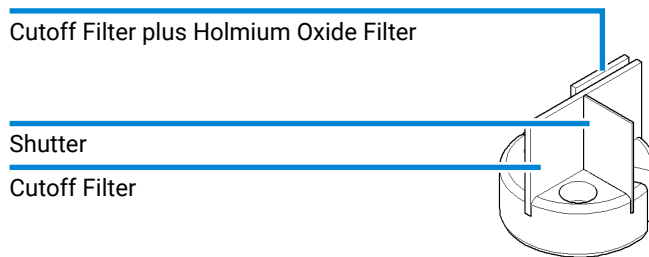


Figure 6: Filter Assembly

Introduction

Operating Principle

The filter assembly has two filters installed and is processor-controlled.

OPEN	nothing in light path at $\lambda < 370$ nm
CUTOFF	cut off filter in light path at $\lambda > 370$ nm
HOLMIUM	holmium oxide filter for wavelength check
SHUTTER	for measurement of dark current of photo diodes

A photo sensor determines the correct position.

Mirror Assemblies M1 and M2

The instrument contains two spherical mirrors (M1 and M2). The beam adjustable is vertically and horizontally. Both mirrors are identical.

Grating Assembly

The grating separates the light beam into all its component wavelengths and reflects the light onto mirror #2.

The stepper motor reference position is determined by a plate fitted onto the motor shaft, interrupting the beam of a photo sensor. The wavelength calibration of the grating is done at the zero order light position and at 656 nm, which is the emission line of the deuterium lamp.

Beam Splitter Assembly

The beam splitter splits the light beam. One part goes directly to the sample diode. The other part of the light beam goes to the reference diode.

Photo Diodes Assemblies

Two photo diode assemblies are installed in the optical unit. The sample diode assembly is located on the left side of the optical unit. The reference diode assembly is located in the front of the optical unit.

Photo Diode ADC (Analog-to-Digital Converter)

The photo diode current is directly converted to digital data direct photo current digitalization. The data is transferred to the detector main board . The photo diode ADC boards are located close to the photo diodes.

Dual-Wavelength Mode

The detector provides a Dual-Wavelength mode that offers additional operation functions.

Features:

- 200 ms acquisition time per data point
 - 5 Hz data rate distributed to two channels
 - 2.5 Hz data rate for each channel,
- delta wavelength max. 150 nm,
- scans during Dual-Wavelength mode are possible,
- the second order filter is disabled when one wavelength is < 370 nm.

Timetable:

- Wavelength settings are timetable programmable (depends if enough time for implementation is available),
- switching from Single-Wavelength mode to Dual-Wavelength mode is NOT timetable programmable,
- filter settings are not timetable programmable.



2

Site Requirements and Specifications

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

Site Requirements 22

Power Consideration 22

Power Cords 23

Bench Space 24

Environment 24

Specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A) 25

Specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B) 28

Specification Conditions 31

Site Requirements

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

Power Consideration

The detector power supply has wide ranging capabilities, see Physical specifications. It accepts any line voltage in the above mentioned range. Consequently, there is no voltage selector in the rear of the detector. There are also no externally accessible fuses, because automatic electronic fuses are implemented in the power supply.

WARNING

Hazard of electrical shock or damage of your instrumentation can result, if the devices are connected to a line voltage higher than specified.

- Connect your instrument to the specified line voltage only.

WARNING

Inaccessible power plug.

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

- Make sure the power connector of the instrument can be easily reached and unplugged.
 - Provide sufficient space behind the power socket of the instrument to unplug the cable.
-

WARNING**Electrical shock hazard**

The module is partially energized when switched off, as long as the power cord is plugged in.

The cover protects users from personal injuries, for example electrical shock.

- Do not open the cover.
 - Do not operate the instrument and disconnect the power cable in case the cover has any signs of damage.
 - Contact Agilent for support and request an instrument repair service.
-

Power Cords

Country-specific power cords are available for the module. The female end of all power cords is identical. It plugs into the power-input socket at the rear. The male end of each power cord is different and designed to match the wall socket of a particular country or region.

Agilent makes sure that your instrument is shipped with the power cord that is suitable for your particular country or region.

WARNING**Unintended use of power cords**

Using power cords for unintended purposes can lead to personal injury or damage of electronic equipment.

- Never use a power cord other than the one that Agilent shipped with this instrument.
 - Never use the power cords that Agilent Technologies supplies with this instrument for any other equipment.
 - Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.
-

WARNING**Absence of ground connection**

The absence of ground connection can lead to electric shock or short circuit.

- Never operate your instrumentation from a power outlet that has no ground connection.
-

WARNING**Electrical shock hazard**

Solvents may damage electrical cables.

- Prevent electrical cables from getting in contact with solvents.
 - Exchange electrical cables after contact with solvents.
-

Bench Space

The detector dimensions and weight (see Physical specifications) allows you to place the detector on almost any desk or laboratory bench. It needs an additional 2.5 cm (1.0 inch) of space on either side and approximately 8 cm (3.1 inch) in the rear for air circulation and electric connections.

If the bench should carry an Agilent 1200 Infinity Series system, make sure that the bench is designed to bear the weight of all modules.

The detector should be operated in a horizontal position.

Environment

Your detector will work within specifications at ambient temperatures and relative humidity as described in Physical specifications.

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.
-

Specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A)

Table 2: Physical specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A)

Type	Specification	Comments
Weight	11 kg (24.3 lbs)	
Dimensions (height x width x depth)	140 x 396 x 436 mm (5.5 x 15.6 x 17.2 inches)	
Line voltage	100–240 V~, ±10%	Wide-ranging capability
Line frequency	50 or 60 Hz, ±5%	
Power consumption	80 VA, 70 W	
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 3000 m (9842 ft)	
Safety standards: IEC, EN, CSA, UL	Overvoltage category II, Pollution degree 2	For indoor use only
ISM classification	ISM Group 1 Class B	According to CISPR 11

Table 3: Performance specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A)

Type	Specification
Detection type	Double-beam photometer
Designed for use with Agilent InfinityLab Assist	Intuitive User Interface, Automated Workflows, Predictive Maintenance & Assisted Troubleshooting
Light source	Deuterium lamp
Number of signals	Single and dual wavelength detection
Maximum data rate	120 Hz (single wavelength detection) 2.5 Hz (dual wavelength detection)

Site Requirements and Specifications

Specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A)

Type	Specification
Short term signal noise (ASTM)	< $\pm 0.25 \cdot 10^{-5}$ AU, at 230 nm (single wavelength detection) < $\pm 0.80 \cdot 10^{-5}$ AU, at 230 nm and 254 nm (dual wavelength detection)
Drift	< $1 \cdot 10^{-4}$ AU/h, at 230 nm
Linear absorbance range	> 2.5 AU upper limit
Wavelength range	190 – 600 nm
Wavelength accuracy	± 1 nm, self-calibration with deuterium lines, verification with holmium oxide filter
Wavelength precision	< ± 0.1 nm
Slit width	6.5 nm typical over whole wavelength range
Time programmable	Wavelength, polarity, peak width, lamp on/off
Flow cells	<i>Standard:</i> 14 μ L volume, 10 mm cell path length and 40 bar (580 psi) pressure maximum <i>Micro:</i> 2 μ L volume, 3 mm cell path length and 120 bar (1760 psi) pressure maximum <i>Semi-micro:</i> 5 μ L volume, 6 mm cell path length and 40 bar (588 psi) pressure maximum <i>Standard Bio:</i> 14 μ L volume, 10 mm cell path length and 40 bar (580 psi) pressure maximum <i>Micro Bio:</i> 2 μ L volume, 3 mm cell path length and 120 bar (1760 psi) pressure maximum <i>Preparative SST:</i> 4 μ L volume, 3 mm cell path length and 120 bar (1760 psi) pressure maximum <i>Preparative Quartz:</i> 4 μ L volume, 3 mm cell path length and 50 bar (725 psi) pressure maximum <i>Preparative Quartz:</i> 0.3 mm cell path length and 50 bar (725 psi) pressure maximum <i>Preparative Quartz:</i> 0.06 mm cell path length and 50 bar (725 psi) pressure maximum <i>High pressure:</i> 14 μ L volume, 10 mm cell path length and 400 bar (5800 psi) pressure maximum
Spectral tools	Stop-flow wavelength scan
Analog output	Recorder/Integrator 100 mV or 1 V, 1 output
Instrument Control	LC and CE Drivers A.02.14 or above Instrument Control Framework (ICF) A.02.04 or above Lab Advisor B.02.08 or above InfinityLab Assist (G7180A) with firmware D.07.40 or above Agilent Instant Pilot (G4208A) B.02.19 or above For details about supported software versions refer to the compatibility matrix of your version of the LC and CE Drivers

Site Requirements and Specifications
Specifications of the 1260 Infinity III Variable Wavelength Detector (G7114A)

Type	Specification
Communication	LAN, Controller Area Network (CAN), USB Extended Remote Interface: ready, start, stop and shut-down signals
GLP	Early maintenance feedback (EMF) for continuous tracking of instrument usage in terms of lamp burn time with user settable limits and feedback messages. Electronic records of maintenance and errors. RFID for electronic records of flow cell and UV lamp conditions (path length, volume, product number, serial number, test passed, and usage). Verification of wavelength accuracy with built-in holmium oxide filter.
Safety and maintenance	Extensive diagnostics, error detection and display through Agilent Instant Pilot and Agilent Lab Advisor software. Leak detection, safe leak handling, leak output signal for shutdown of pumping system. Low voltages in major maintenance areas. Tracking of flow cells and lamps with RFID (radio frequency identification) tags.
Housing	All materials recyclable.
Others	Second generation of Electronic temperature control (ETC) for the complete optical unit.

Specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B)

Table 4: Physical specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B)

Type	Specification	Comments
Weight	11 kg (24.3 lbs)	
Dimensions (height x width x depth)	140 x 396 x 436 mm (5.5 x 15.6 x 17.2 inches)	
Line voltage	100–240 V~, ±10%	Wide-ranging capability
Line frequency	50 or 60 Hz, ±5%	
Power consumption	80 VA, 70 W	
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 3000 m (9842 ft)	
Safety standards: IEC, EN, CSA, UL	Overvoltage category II, Pollution degree 2	For indoor use only
ISM classification	ISM Group 1 Class B	According to CISPR 11

Table 5: Performance specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B)

Type	Specification
Detection type	Double-beam photometer
Designed for use with Agilent InfinityLab Assist	Intuitive User Interface, Automated Workflows, Predictive Maintenance & Assisted Troubleshooting
Light source	Deuterium lamp
Number of signals	Single and dual wavelength detection
Maximum data rate	240 Hz (single wavelength detection) 2.5 Hz (dual wavelength detection)

Site Requirements and Specifications

Specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B)

Type	Specification
Short term signal noise (ASTM)	< $\pm 0.15 \cdot 10^{-5}$ AU, at 230 nm (single wavelength detection) < $\pm 0.80 \cdot 10^{-5}$ AU, at 230 nm and 254 nm (dual wavelength detection)
Drift	< $1 \cdot 10^{-4}$ AU/h, at 230 nm
Linear absorbance range	> 2.5 AU upper limit
Wavelength range	190 – 600 nm
Wavelength accuracy	± 1 nm, self-calibration with deuterium lines, verification with holmium oxide filter
Wavelength precision	< ± 0.1 nm
Slit width	6.5 nm typical over whole wavelength range
Time programmable	Wavelength, polarity, peak width, lamp on/off
Flow cells	<i>Standard</i> : 14 μ L volume, 10 mm cell path length and 40 bar (580 psi) pressure maximum <i>Micro</i> : 2 μ L volume, 3 mm cell path length and 120 bar (1760 psi) pressure maximum <i>Semi-micro</i> : 5 μ L volume, 6 mm cell path length and 40 bar (580 psi) pressure maximum <i>Standard Bio</i> : 14 μ L volume, 10 mm cell path length and 40 bar (580 psi) pressure maximum <i>Micro Bio</i> : 2 μ L volume, 3 mm cell path length and 120 bar (1760 psi) pressure maximum <i>Preparative SST</i> : 4 μ L volume, 3 mm cell path length and 120 bar (1760 psi) pressure maximum <i>Preparative Quartz</i> : 4 μ L volume, 3 mm cell path length and 50 bar (725 psi) pressure maximum <i>Preparative Quartz</i> : 0.3 mm cell path length and 50 bar (725 psi) pressure maximum <i>Preparative Quartz</i> : 0.06 mm cell path length and 50 bar (725 psi) pressure maximum <i>High pressure</i> : 14 μ L volume, 10 mm cell path length and 400 bar (5800 psi) pressure maximum
Spectral tools	Stop-flow wavelength scan
Analog output	Recorder/Integrator 100 mV or 1 V, 1 output
Instrument Control	LC and CE Drivers A.02.11 or above Instrument Control Framework (ICF) A.02.04 or above Lab Advisor B.02.06 or above InfinityLab Assist (G7180A) with firmware D.07.40 or above Agilent Instant Pilot (G4208A) B.02.19 or above For details about supported software versions refer to the compatibility matrix of your version of the LC and CE Drivers

Site Requirements and Specifications

Specifications of the 1290 Infinity III Variable Wavelength Detector (G7114B)

Type	Specification
Communication	LAN, Controller Area Network (CAN), USB Extended Remote Interface: ready, start, stop and shut-down signals
GLP	Early maintenance feedback (EMF) for continuous tracking of instrument usage in terms of lamp burn time with user settable limits and feedback messages. Electronic records of maintenance and errors. RFID for electronics records of flow cell and UV lamp conditions (path length, volume, product number, serial number, test passed, and usage). Verification of wavelength accuracy with built-in holmium oxide filter.
Safety and maintenance	Extensive diagnostics, error detection and display through Agilent Instant Pilot and Agilent Lab Advisor software. Leak detection, safe leak handling, leak output signal for shutdown of pumping system. Low voltages in major maintenance areas. Tracking of flow cells and lamps with RFID (radio frequency identification) tags.
Housing	All materials recyclable.
Others	Second generation of Electronic temperature control (ETC) for the complete optical unit.

Specification Conditions

Following many of the principles outlined in ASTM method E165798.

Reference conditions: Standard flow cell, path length 10 mm, flow 1 mL/min LC-grade methanol.

Noise:

$\pm 1.5 \cdot 10^{-6}$ AU at 230 nm, TC 2 s

RT = 2.2 * TC

Linearity:

Linearity is measured with caffeine at 273 nm.

NOTE

The specification are based on the the standard RFID tag lamp (G1314-60101) and may be not achieved when other lamp types or aged lamps are used.

ASTM drift tests require a temperature change below 2°C/hour (3.6°F/hour) 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.

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



3 Installation

The installation of the module will be done by an Agilent service representative. In this chapter, only installation of user-installable options and accessories are described.

Installing Capillaries 33

Install Capillaries 33

Handling Leak and Waste 38

Drain Connectors Installation 41

Waste Concept 46

Waste Guidance 46

Leak Sensor 47

Connecting Modules and Control Software 48

Instrument Configuration 49

Installing Capillaries

This section provides information on how to install capillaries and fittings.

Install Capillaries

Capillaries and connections depend on which system is installed.

NOTE

As you move to smaller-volume, high-efficiency columns, you will want to use narrow id tubing, as opposed to the wider id tubing used for conventional HPLC instruments.

NOTE

Agilent capillaries are color-coded for quick identification, see [At-a-Glance Details About Agilent Capillaries](#) on page 254.

Table 6: Capillary connections for 1260 Infinity III systems

p/n	From	To
G7120-60007 (Bottle Head Assembly)	Solvent Bottle	Infinity III Pump
5500-1246 (Capillary ST 0.17 mm x 500 mm SI/SI)	Pump	Sampler
5500-1217 (Capillary, ST, 0.17 mm x 900 mm SI/SX)	Pump	Vialsampler with ICC
5500-1246 (Capillary ST 0.17 mm x 500 mm SI/SI)	Multisampler	MCT Valve/Heat Exchanger
5500-1252 (Capillary, ST, 0.17 mm x 400 mm SL/SL)	Vialsampler	MCT Valve/Heat Exchanger
5500-1240 (Capillary ST 0.17 mm x 105 mm SL/SL)	Vialsampler	ICC Heat Exchanger
5500-1250 (Capillary, ST, 0.17 mm x 120 mm SL/SL, long socket)	ICC Heat Exchanger	Column
5500-1193 (InfinityLab Quick Turn Capillary ST 0.17 mm x 105 mm, long socket)	MCT Heat Exchanger	Column
5500-1191 (InfinityLab Quick Turn Capillary ST 0.12 mm x 280 mm, long socket)	Column/MCT Valve	Detector
5062-8535 (Waste accessory kit (Flow Cell to waste))	VWD	Waste
5062-2462 (Tube PTFE 0.7 mm x 5 m, 1.6 mm od)	DAD/FLD	Waste
G5664-68712 (Analytical tubing kit 0.25 mm i.d. PTFE-ESD)	Detector	Fraction Collector

Table 7: Capillary connections for 1290 Infinity III systems

p/n	From	To
G7120-60007 (Bottle Head Assembly)	Solvent Bottle	Infinity III Pump
5500-1245 (Capillary ST 0.17 mm x 400 mm SI/SI)	Pump	Sampler

p/n	From	To
5500-1217 (Capillary, ST, 0.17 mm x 900 mm SI/SX)	Pump	Vialsampler with ICC
5500-1157 (Capillary ST 0.12 mm x 500 mm SL/S)	Multisampler	MCT Valve/Heat Exchanger
5500-1251 (Capillary ST 0.12 mm x 400 mm SL/SL)	Vialsampler	MCT Valve/Heat Exchanger
5500-1238 (Capillary ST 0.12 mm x 105 mm SL/SL)	Vialsampler	ICC Heat Exchanger
5500-1249 (Capillary ST 0.12 mm x 120 mm SL/SL, long socket)	ICC Heat Exchanger	Column
5500-1201 (Capillary ST 0.12 mm x 105 mm SL)	MCT Heat Exchanger	Column
5500-1191 (InfinityLab Quick Turn Capillary ST 0.12 mm x 280 mm, long socket)	Column/MCT Valve	Detector
5062-8535 (Waste accessory kit (Flow Cell to waste))	VWD	Waste
5062-2462 (Tube PTFE 0.7 mm x 5 m, 1.6 mm od)	DAD/FLD	Waste
G5664-68712 (Analytical tubing kit 0.25 mm i.d. PTFE-ESD)	Detector	Fraction Collector

Table 8: Capillary connections for 1260 Infinity III Bio-inert LC

p/n	From	To
G7120-60007 (Bottle Head Assembly)	Solvent Bottle	Infinity III Pump
5500-1264 (Capillary Ti 0.17 mm x 500 mm, SL/SLV)	Pump	Multisampler
G5667-81005 (Capillary PK/ST 0.17 mm x 500 mm, RLO/RLO (Bio-inert))	Multisampler	MCT
5067-4741 (ZDV union (Bio-inert))	Capillary	Bio-inert Heat Exchanger
G7116-60041 (Quick Connect Heat Exchanger Bio-inert)		
0890-1763 (Capillary PEEK 0.18 mm x 1.5 m) and 5063-6591 (PEEK Fittings 10/PK)	Column/MCT Valve	Detector
5062-8535 (Waste accessory kit (Flow Cell to waste))	VWD	Waste
5062-2462 (Tube PTFE 0.7 mm x 5 m, 1.6 mm od)	DAD/FLD	Waste
G5664-68712 (Analytical tubing kit 0.25 mm i.d. PTFE-ESD)	Detector	Fraction Collector

Table 9: Capillary connections for 1290 Infinity III Bio LC

p/n	From	To
G7120-60007 (Bottle Head Assembly)	Solvent Bottle	Infinity III Pump

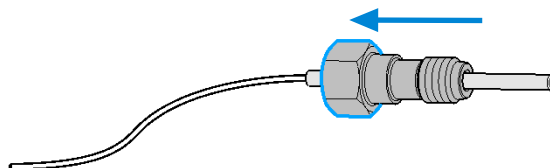
p/n	From	To
5500-1419 (Capillary MP35N 0.17 mm x 500 mm, SI/SI)	Pump	Multisampler
5500-1279 (Capillary MP35N 0.12 mm x 500 mm SI/SI)	Multisampler	MCT
5500-1578 (Quick Connect Capillary MP35N 0.12 mm x 105 mm)	MCT Heat Exchanger	Column
5500-1596 (Quick Turn Capillary MP35N 0.12 mm x 280 mm)	Column/MCT Valve	Detector (DAD)
5500-1598 (Quick Turn Capillary MP35N 0.12 mm x 500 mm)	Column/MCT Valve	Detector (VWD)
5062-8535 (Waste accessory kit (Flow Cell to waste))	VWD	Waste
5062-2462 (Tube PTFE 0.7 mm x 5 m, 1.6 mm od)	DAD/FLD	Waste
G5664-68712 (Analytical tubing kit 0.25 mm i.d. PTFE-ESD)	Detector	Fraction Collector

For correct installation of capillary connections it's important to choose the correct fittings, see Syntax for Capillary Description.

- 1 Select a nut that is long enough for the fitting you'll be using.



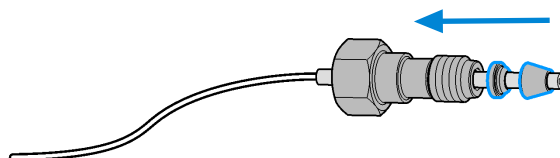
- 2 Slide the nut over the end of the tubing or capillary.



Installation

Installing Capillaries

- 3 Carefully slide the ferrule components on after the nut and then finger-tighten the assembly while ensuring that the tubing is completely seated in the bottom of the end fitting.

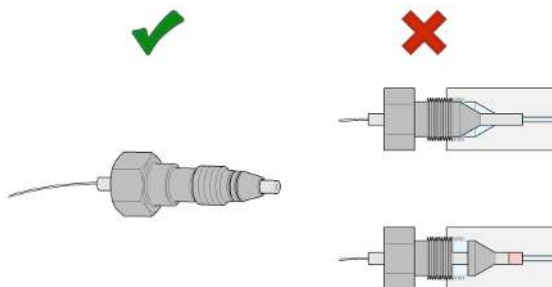


- 4 Use a stable port installed to the module to gently tighten the fitting facing to the module. Or use the column to tighten the fitting facing to the column. This measure forces the ferrule to seat onto the tubing or capillary.

NOTE

Do not overtighten. Over-tightening will shorten the lifetime of the fitting.

- 5 Loosen the nut and verify that the ferrule is correctly positioned on the tubing or capillary.



NOTE

The first time that the Swagelok fitting is used on a column or an injection valve, the position of the ferrule is permanently set. If changing from a column or an injection valve to another, the fitting may leak or decrease the quality of the separation by contributing to band broadening.

For Bio and Bio-Inert Systems, the Swagelok instructions do not apply.

Handling Leak and Waste

The Agilent InfinityLab LC Series has been designed for safe leak and waste handling. It is important that all security concepts are understood and instructions are carefully followed.

The solvent cabinet is designed to store a maximum volume of 8 L solvent. The maximum volume for an individual bottle stored in the solvent cabinet should not exceed 2 L. For details, see the usage guideline for the Agilent Infinity III Solvent Cabinets (a printed copy of the guideline has been shipped with the solvent cabinet, electronic copies are available on the Internet).

All leak plane outlets are situated in a consistent position so that all Infinity and Infinity II/III modules can be stacked on top of each other. Waste tubes are guided through a channel on the right hand side of the instrument, keeping the front access clear from tubes.

The leak plane provides leak management by catching all internal liquid leaks, guiding them to the leak sensor for leak detection, and passing them on to the next module below, if the leak sensor fails. The leak sensor in the leak plane stops the running system as soon as the leak detection level is reached.

Solvent and condensate is guided through the waste channel into the waste container:

- from the detector's flow cell outlet
- from the Multisampler needle wash port
- from the Sample Thermostat (condensate)
- from the pump's Seal Wash Sensor (if applicable)
- from the pump's Purge Valve or Multipurpose Valve

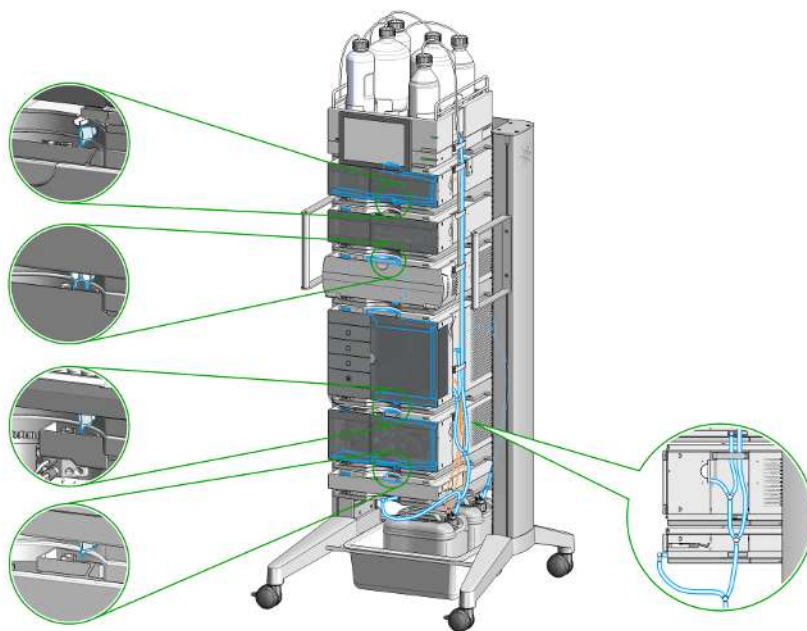


Figure 7: Infinity III Leak Waste Concept (Flex Bench installation)

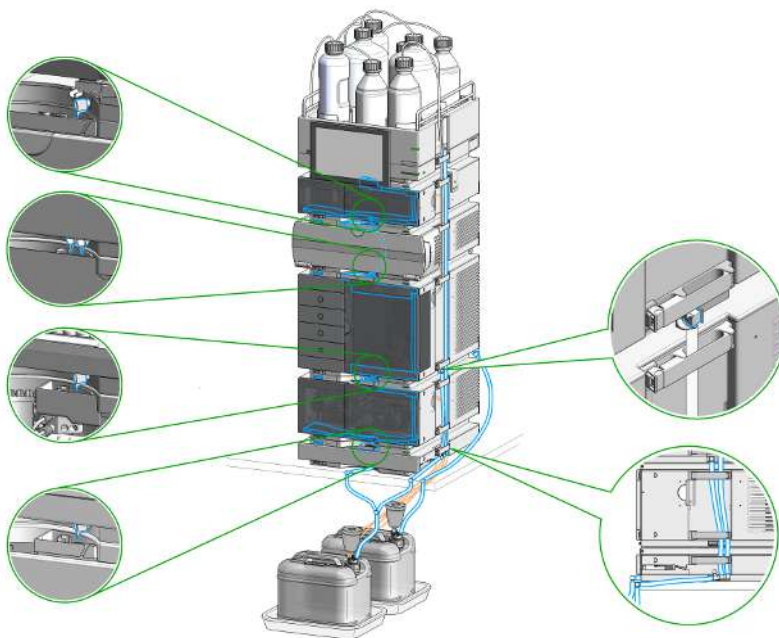


Figure 8: Infinity III Single Stack Leak Waste Concept (bench installation)

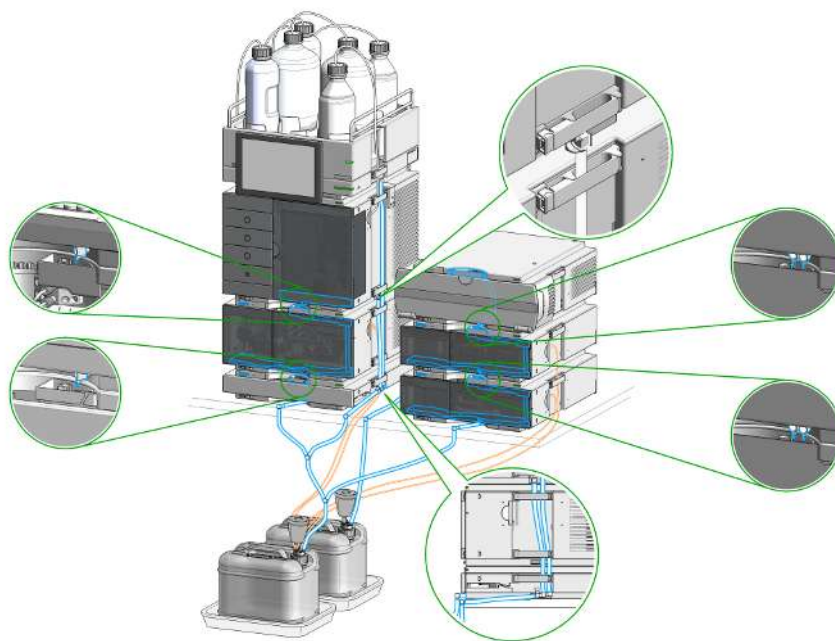


Figure 9: Infinity III Two Stack Leak Waste Concept (bench installation)

The waste tube connected to the leak plane outlet on each of the bottom instruments guides the solvent to a suitable waste container.

Drain Connectors Installation

Drain Connectors have been developed to improve leak drainage for low flow leaks of high viscosity solvents (for example, isopropanol) in Agilent InfinityLab LC Series Systems. Install these parts to modules where they are missing (usually preinstalled).

- Make sure that dripping adapters are correctly installed on each module in the LC stack, excluding lowest module.
- Remove the dripping adapter if it is appeared to be installed on the lowest module in the LC stack and connect waste tube instead.
- Consider 5004-0000 (Drain Connectors Kit) if drain adaptor is missing on some module(s).

For illustration, see [Handling Leak and Waste](#) on page 38.

Parts required

Qty.	p/n	Description
	 5004-0000	Drain Connectors Kit

Content of Drain Connectors Kit (p/n 5004-0000)

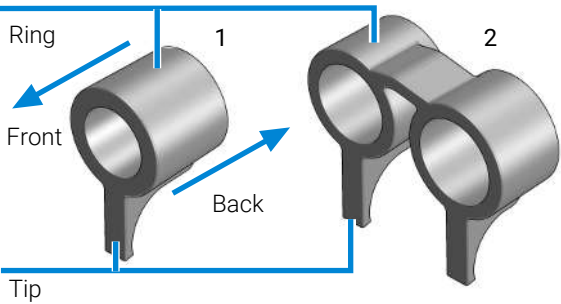


Figure 10: Overview of Drain Connectors: Single (left) and Double (right)



Qty.	p/n	Description
Parts can be ordered only as a complete kit.		
3	 5043-1834	Single Drain Connector ID3.0-Long
1	 5043-1836	Double Drain Connector-Long

Table 10: Compatibility of drain connectors and modules

Drain Connector Type	Compatible Module	Compatible Module Type
Double	G7116A/B	Column Compartment
Single	G7114A/B	Detector
	G7115A	
	G7117A/B/C	
	G7121A/B	
	G7162A/B	
	G7165A	
	G7129A/B/C	Sampler
	G7167A/B/C	
	G5668A	
	G7137A	
	G7157A	
	G4767A	
	G7122A	Degasser
	G7104A/C	
	G7110B	Pump
	G7111A/B	
	G7112B	
	G7120A	
	G7131A/C	
	G7132A	
	G5654A	
	G4782A	

Preparations

- Leak drains of LC modules are clean and free of salt or solvent residuals.

NOTE

Do not install drain connectors on the bottom modules of the stack. Drain outlet of the bottom module has to be connected via waste tubing to a suitable waste container (see Leak and Waste Handling in the manual for a respective module).

NOTE

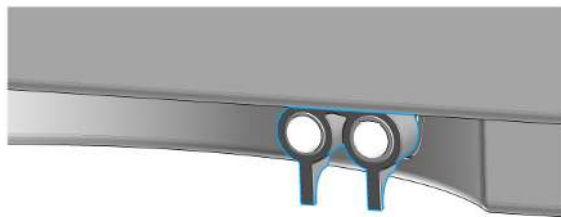
In case of incorrect installation, drain connectors cannot fully perform the intended function.

NOTE

It is not required to power off the HPLC stack to install Single and Double Drain Connectors. The installation of the connectors does not affect the analysis performed during the installation.

**Install the Double Drain Connector on the leak drain of the
1260 Infinity III Multicolumn Thermostat (G7116A)/
1290 Infinity III Multicolumn Thermostat (G7116B)**

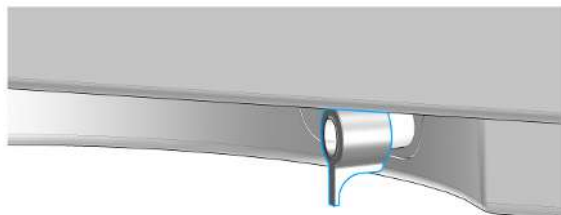
- 1 Align the rings with the leak drain outlets of the module, press slightly with the fingers, and slide the connector along the leak drain outlets until it is aligned with the front of the leak drain.

**Install Single Drain Connectors on other modules in the LC stack**

Installation

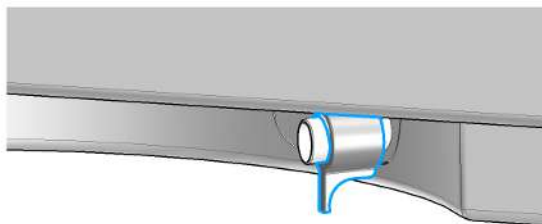
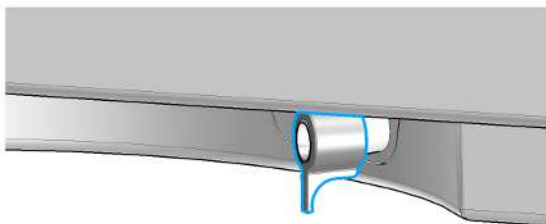
Handling Leak and Waste

- 1 Align the ring with the leak drain outlet of the module, press slightly with the fingers, and slide the connector along the leak drain outlet until it is aligned with the front of the leak drain.



Make sure that the following requirements are covered:

- The tip of the drain connector points straight down.
- The leak drain outlets and the drain connectors are aligned properly.



Waste Concept

Agilent recommends using the 5043-1221 (6 L waste can with 1 Stay Safe cap GL45 with 4 ports) for optimal and safe waste disposal. If you decide to use your own waste solution, make sure that the tubes don't immerse in the liquid.



Waste Guidance



NOTE

The waste drainage must go straight into the waste containers. The waste flow must not be restricted at bends or joints.

Leak Sensor

CAUTION

Solvent incompatibility

The solvent DMF (dimethylformamide) leads to corrosion of the leak sensor. The material of the leak sensor, PVDF (polyvinylidene fluoride), is incompatible with DMF.

- Do not use DMF as mobile phase.
- Check the leak sensor regularly for corrosion.

Connecting Modules and Control Software

WARNING**Use of unsupplied cables**

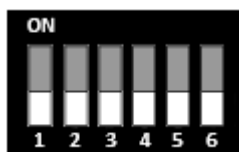
Using cables not supplied by Agilent Technologies can lead to damage of the electronic components or personal injury.

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

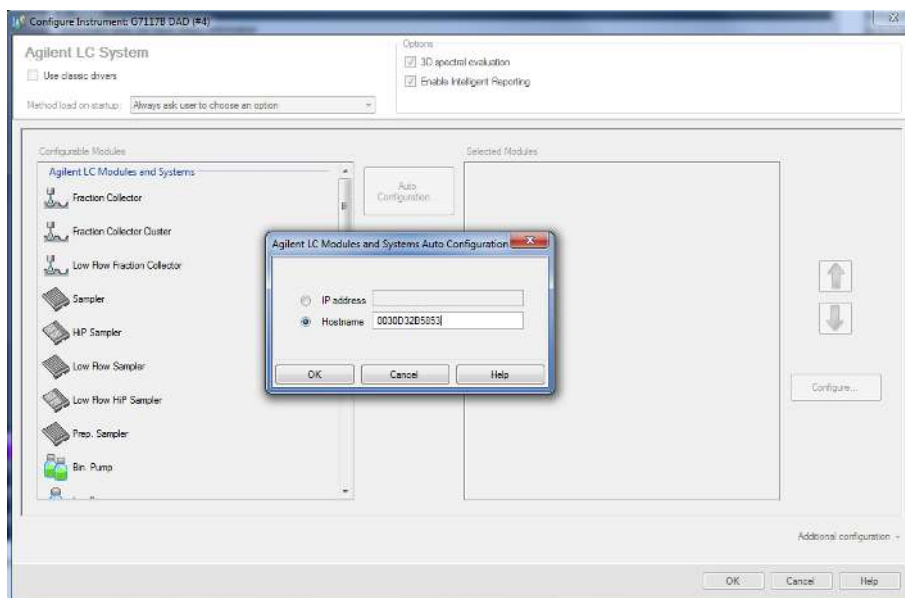
Instrument Configuration

Example shows an instrument configuration with a Diode Array Detector.

- 1 Set the switches of the Configuration switch at the rear of the module:
 - a All switches DOWN: module uses the default IP address 192.168.254.11.



- b Switch 4 UP and others DOWN: module uses DHCP.
 - c Switch 5 UP and others DOWN: modules uses STORED address.
- 2 Enter the setup information (MAC ¹ / IP address and/or Instrument Name).
 - a Agilent OpenLab ChemStation (Configure Instrument):



¹ MAC address can only be used in DHCP DIP-switch configuration.

Installation

Instrument Configuration


b Lab Advisor (Instrument Overview - Add Instrument):

System Properties

System

System Name: Description:

Instruments

Instrument Name:	Instrument Address:	Instrument Type:
 <input type="text" value="G7117B"/>	<input type="text" value="0030D32B5853"/> <input type="text" value="192.168.254.11"/>	<input type="text" value="Agilent LC/CE"/>

Reconnect: ☐



4 Using the Module

This chapter provides information on how to use the module.

General Information 52

Turn On/Off 52

Status Indicators 54

Preparation of the System 56

Prepare a Run 56

Prime and Purge the System 63

Preparing the Detector 65

Preparing the Module 66

The Detector User Interface 66

Detector Control Settings 68

Method Parameter Settings 69

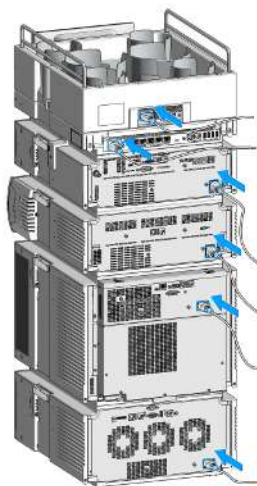
Scanning with the VWD 72

General Information

Turn On/Off

This procedure exemplarily shows an arbitrary LC stack configuration.

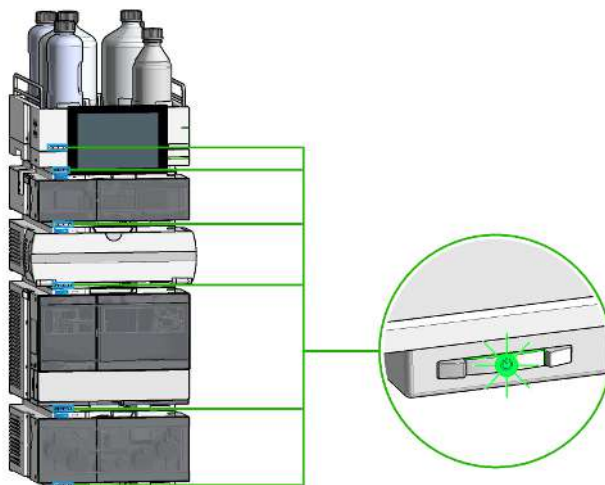
1



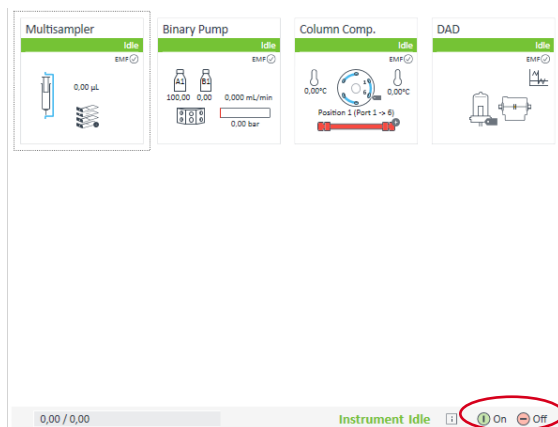
Using the Module

General Information

2 On/Off switch: On



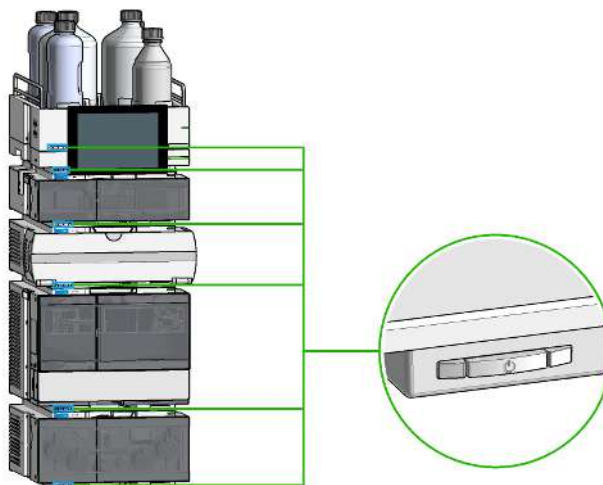
3 Turn instrument On/Off with the control software.



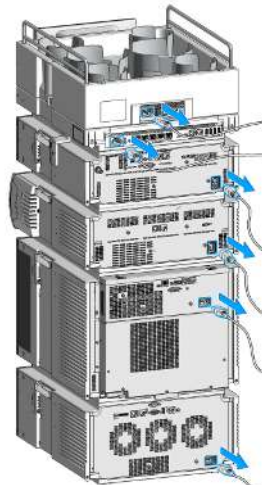
Using the Module

General Information

4 On/Off switch: Off



5



Status Indicators

The module status indicator indicates one of six possible module conditions.

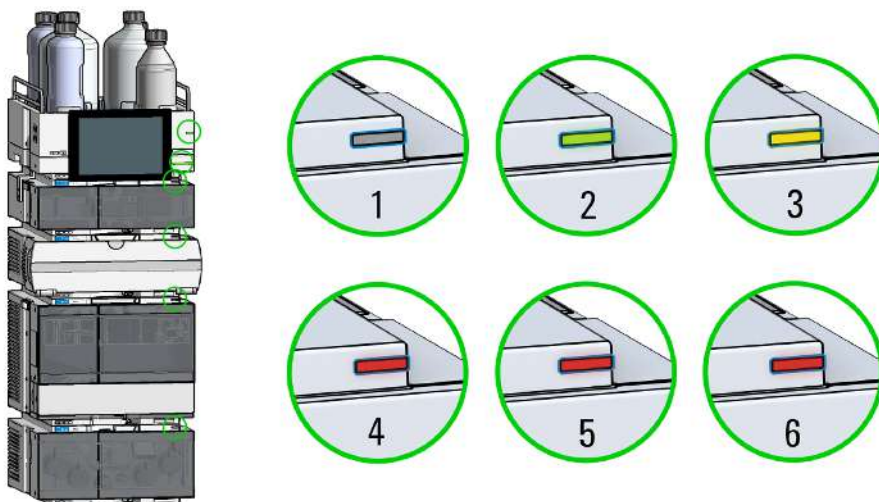


Figure 11: Arbitrary LC stack configuration (example)

1	Idle
2	Run mode
3	Not-ready. Waiting for a specific pre-run condition to be reached or completed.
4	Error mode - interrupts the analysis and requires attention (for example, a leak or defective internal components).
5	Resident mode (blinking) - for example, during update of main firmware.
6	Bootloader mode (fast blinking). Try to re-boot the module or try a cold-start. Then try a firmware update.

InfinityLab Assist Hub Status Indicator

The Assist Hub status indicator displays the status of the entire system. If a module in the system is not ready (yellow), the Assist Hub status indicator also shows not ready (yellow). The same applies for the module conditions **Idle**, **Run mode**, and **Error mode**.

Preparation of the System

Prepare a Run

This procedure exemplarily shows how to prepare a run. Parameters as shown in the screenshots may vary, depending on the system installed.

WARNING

Toxic, flammable and hazardous solvents, samples and reagents

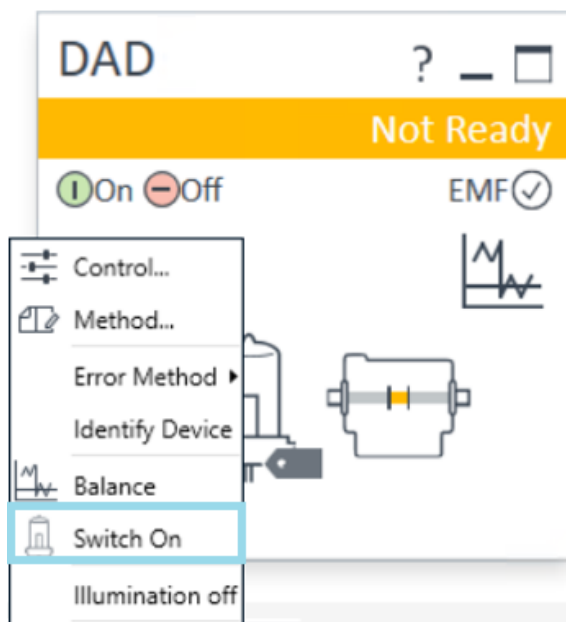
The handling of solvents, samples and reagents can hold health and safety risks.

- 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.
- Do not use solvents with an auto-ignition temperature below 200 °C (392 °F). Do not use solvents with a boiling point below 56 °C (133 °F).
- Avoid high vapor concentrations. Keep the solvent temperature at least 40 °C (72 °F) below the boiling point of the solvent used. This includes the solvent temperature in the sample compartment. For the solvents methanol and ethanol keep the solvent temperature at least 25 °C (45 °F) below the boiling point.
- Do not operate the instrument in an explosive atmosphere.
- Do not use solvents of ignition Class IIC according IEC 60079-20-1 (for example, carbon disulfide).
- Reduce the volume of substances to the minimum required for the analysis.
- Never exceed the maximum permissible volume of solvents (8 L) in the solvent cabinet. Do not use bottles that exceed the maximum permissible volume as specified in the usage guideline for solvent cabinet.
- Ground the waste container.
- Regularly check the filling level of the waste container. The residual free volume in the waste container must be large enough to collect the waste liquid.
- To achieve maximal safety, regularly check the tubing for correct installation.

Using the Module

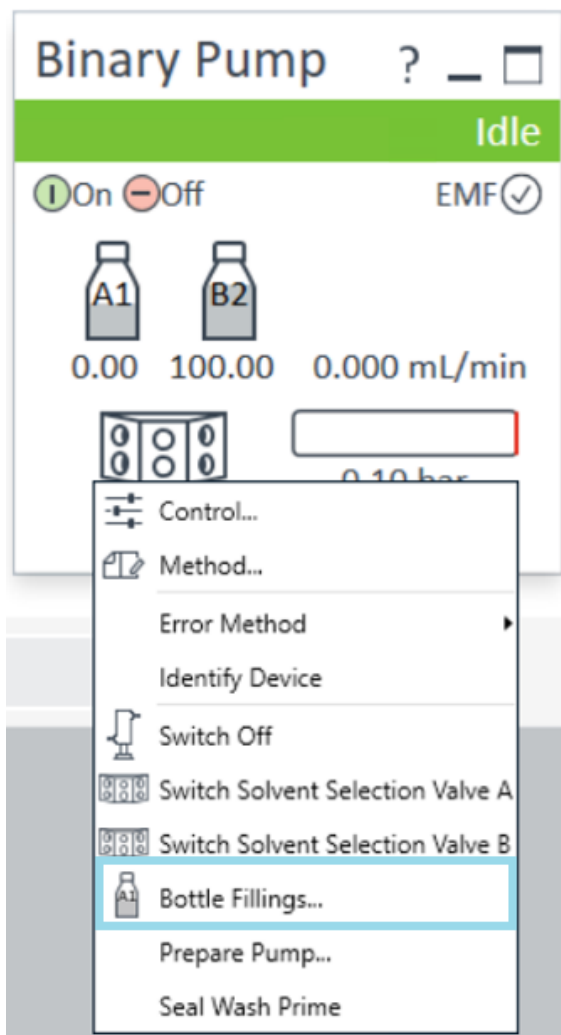
Preparation of the System

- 1 Switch on the detector.



- 2 Fill the solvent bottles with adequate solvents for your application.
- 3 Place solvent tubings with bottle head assemblies into the solvent bottles.
- 4 Place solvent bottles into the solvent cabinet.

- 5 Solvent bottle filling dialog (in the software).



Using the Module

Preparation of the System

Bottle Fillings

Solvent Bottle

Fillings

	Actual Volume		Total Volume	
A1	0.22	liter	1.00	liter
A2	0.29	liter	1.00	liter
B1	0.16	liter	1.00	liter
B2	0.49	liter	1.00	liter

Actions

☐ Prevent analysis if level falls below liter

☐ Turn pump off if running out of solvent

Waste Bottle

Filling

	Actual Volume		Total Volume	
Waste bottle:	0.00	liter	0.00	liter

Actions

☐ Prevent analysis if level raises above liter

☐ Turn pump off if waste volume has reached maximum limit

Ok Cancel Help

6 Purge the pump.

NOTE

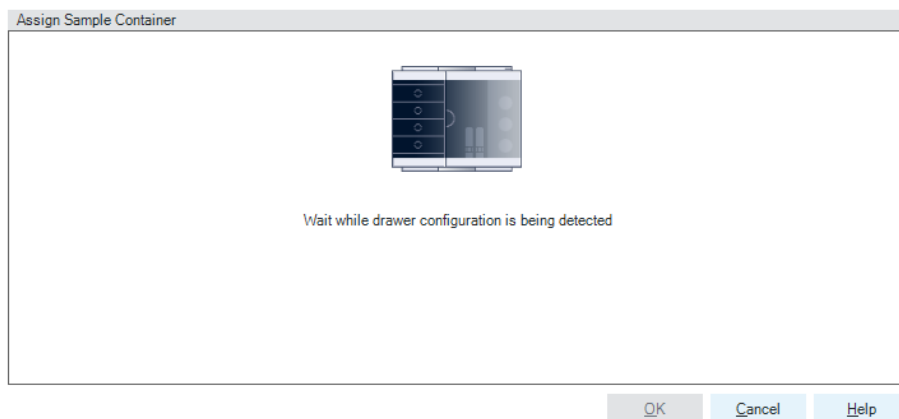
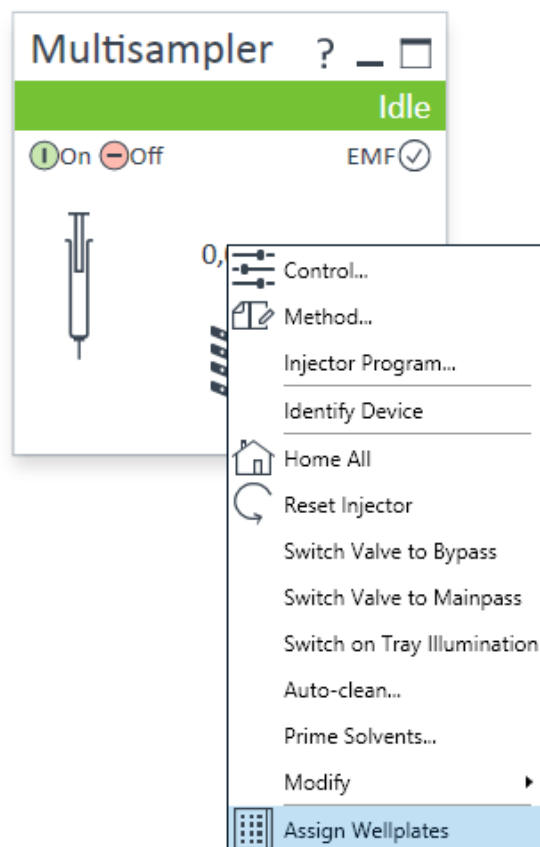
For details on priming and purging, refer to the technical note *Best Practices for Using an Agilent LC System Technical Note (InfinityLab-BestPractice-en-SD-29000194.pdf, SD-29000194)*.

7 Change solvent type if necessary.

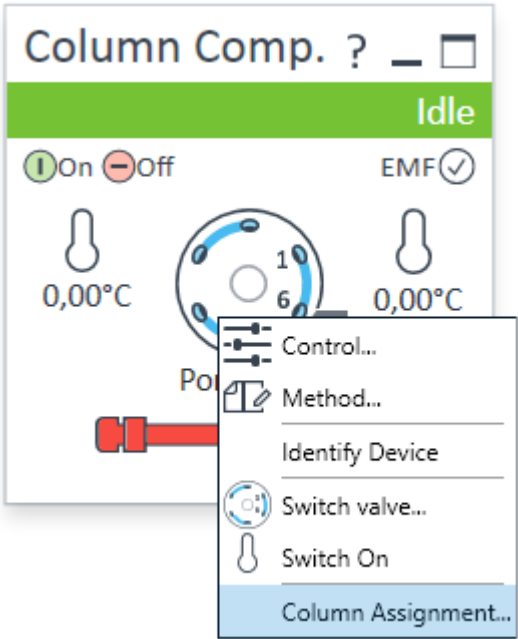
Using the Module

Preparation of the System

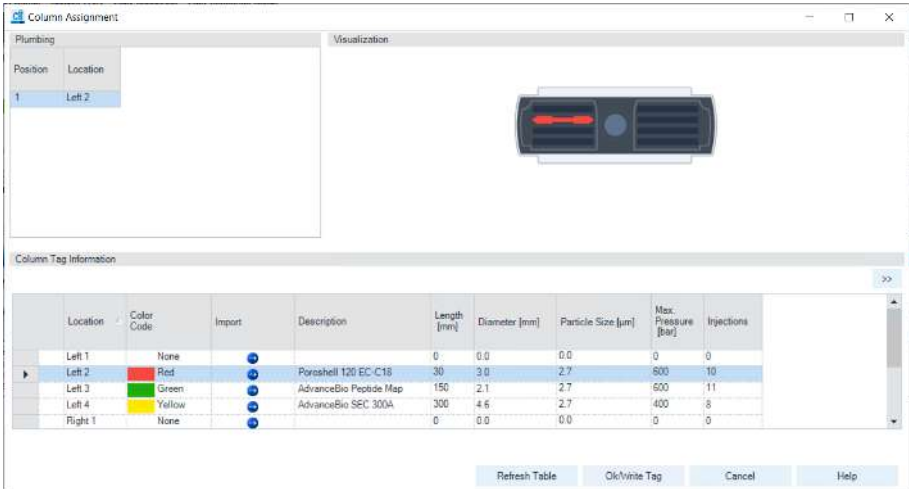
- 8 Choose the tray format of the sampler.



9 Add a new column.



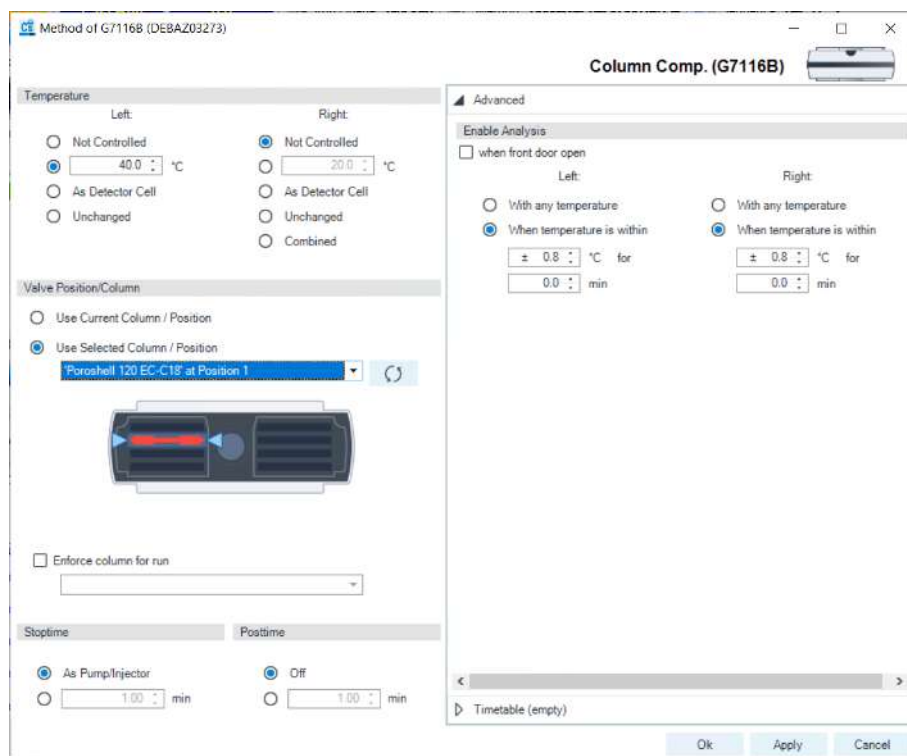
10 Enter the column information.



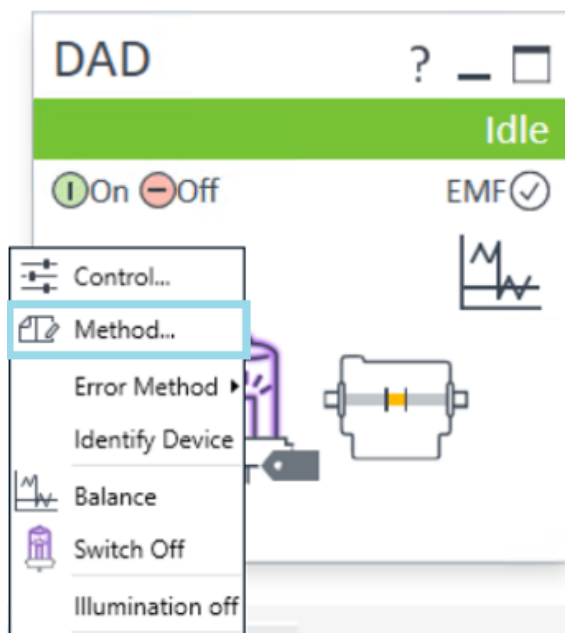
Using the Module

Preparation of the System

11 Select the column in the Method settings of the column compartment.



12 Set the detector parameters according to the needs of your method.



Prime and Purge the System

When the solvents have been exchanged or the pumping system has been turned off for a certain time (for example, overnight) oxygen will re-diffuse into the solvent channel between the solvent reservoir, vacuum degasser (when available in the system) and the pump. Solvents containing volatile ingredients will slightly lose these. Therefore priming of the pumping system is required before starting an application.

Table 11: Choice of priming solvents for different purposes

Activity	Solvent	Comments
After an installation	Isopropanol	Best solvent to flush air out of the system
When switching between reverse phase and normal phase (both times)	Isopropanol	Best solvent to flush air out of the system
After an installation	Ethanol or Methanol	Alternative to Isopropanol (second choice) if no Isopropanol is available
To clean the system when using buffers	Bidistilled water	Best solvent to re-dissolve buffer crystals
After a solvent change	Bidistilled water	Best solvent to re-dissolve buffer crystals
After the installation of normal phase seals (P/N 0905-1420)	Hexane + 5% Isopropanol	Good wetting properties

NOTE

The pump should never be used for priming empty tubings (never let the pump run dry). Use a syringe to draw enough solvent for completely filling the tubings to the pump inlet before continuing to prime with the pump.

- 1 Open the purge valve of your pump (by turning it counterclockwise) and set flow rate to 3 – 5 mL/min.
- 2 Flush all tubes with at least 30 mL of solvent.
- 3 Set flow to required value of your application and close the purge valve.

NOTE

Pump for approximately 10 minutes before starting your application.

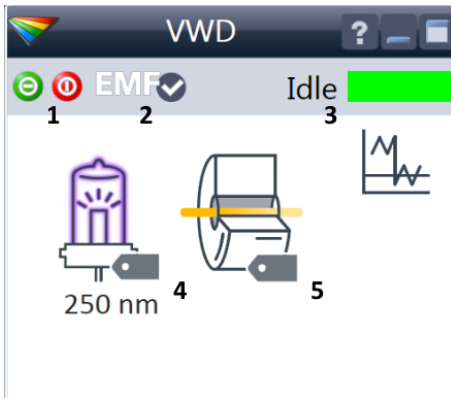
Preparing the Detector

For best performance of the detector

- Let the lamp warm-up and stabilize for at least one hour (initial turn on of the module requires a longer time depending on the environment and the application needs); refer to [Specification Conditions](#) on page 31.
- For high sensitivity measurements, a stable environment is required; refer to [Environment](#) on page 24. Prevent drafts from air condition systems.
- Do not work with removed/open front panels/doors. When the system includes a G1316 TCC (typically located below the detector) and its front panel is removed while the TCC is set to high temperatures, the up-streaming air could influence the stability of the detector baseline.

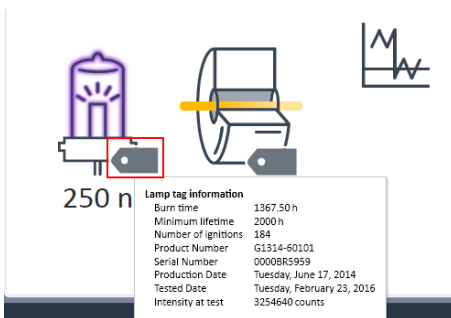
Preparing the Module

The Detector User Interface



Within the detector GUI, there are active areas. If you move the mouse cursor across the icons the cursor will change.

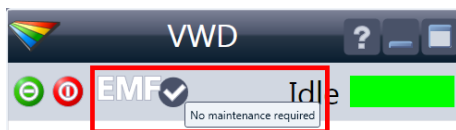
1. Lamp: turn on and off of UV-lamp
2. EMF status
3. Detector status
4. Lamp status (on/off) and information (RFID tag)
5. Flow Cell information (RFID tag)



RFID tag information is displayed when moving with the mouse cursor on to the tag attached to the flow cell or lamp. The information provides flow cell and lamp related information like

- Part number
- Production date
- Serial number

and other details.



EMF Status shows Run / Ready / Error state and “Not Ready text” or “Error text”

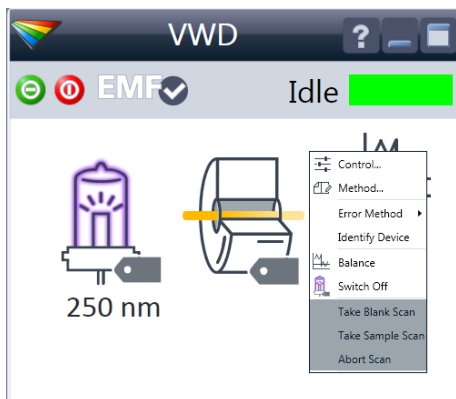
- Offline (gray)
- Ok. No Maintenance required (green)
- EMF warning. Maintenance might be required (yellow)
- EMF warning. Maintenance required (red)

Important: The EMF settings can be accessed via Agilent Lab Advisor. The limit(s) can be changed. Based on the limit, the User Interface displays the above status.



Module Status shows Run / Ready / Error state and “Not Ready text” or “Error text”

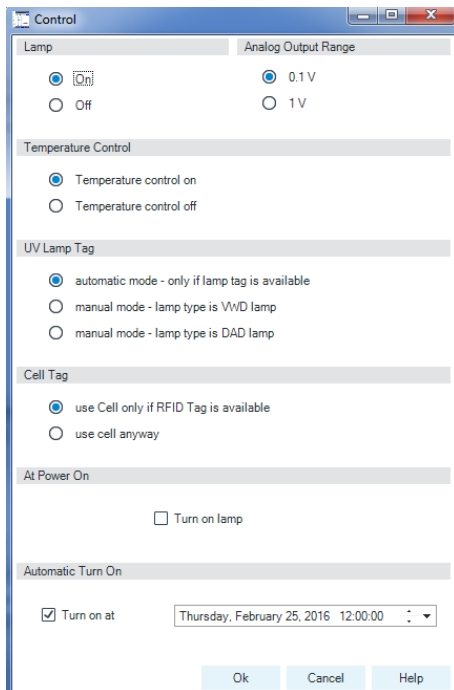
- Error (red)
- Not ready (yellow)
- Ready (green)
- Pre run, Post run (purple)
- Run (blue)
- Idle (green)
- Offline (dark gray)
- Standby (light gray)



A right-click into the Active Area will open a menu to

- Show the Control Interface (special module settings)
- Show the Method interface (similar as via menu **Instrument > Setup Instrument Method**)
- Set Error Method
- Identify Module (Status LED will blink)
- Perform a Balance
- Switch the UV-lamp on/off (same as click on button “Make Device Ready/Turn device off (standby)”))
- Take / Abort Scans (during flow off)

Detector Control Settings



The figure shows the default settings.

- **Lamps:** can be turned ON/OFF.
- **Analog Output Range:** can be set to either 100 mV or 1 V full scale, for additional settings see “Analog” (under [Table 12](#) on page 70).
- **Temperature Control:** The optical unit is kept on constant temperature (some degrees above ambient) and improves the baseline stability in unstable environments, see also [Environment](#) on page 24.
If the flow cell temperature is critical for your chromatography or your environment is stable, you may set the Temperature Control to off. This will lower the optical unit and flow cell temperature by some degree.
- **UV Lamp Tag**
 - **Automatic** detects a lamp with RFID tag. If no RFID tag lamp is used, “UV lamp not ready” is displayed and it cannot be ignited. A compatible mode has to be selected based on the used lamp; see Non-RFID-tag lamp information below.
 - **Manual** (by PN) uses the selected “heating” mode. This mode can also be used when the RFID tag of the standard lamp (G1314-60101) is not recognized (defect RFID tag or reader).
 - **Non-RFID-tag lamp:** In case a non-RFID-tag lamp is used, the user interface will show this when selecting a compatible mode. You may operate the detector outside of the guaranteed specification. The correct selection is important for optimal performance and lifetime.
- **Cell Tag:** Automatic mode for Agilent flow cells with RFID tags. If no RFID tag cell is used, detector icon will become gray (cell tag not ready) and analysis is disabled.
- **At Power On:** automatic lamp-on at power on.
- **Automatic Turn On:** automatic detector power on.

Method Parameter Settings

These settings are available via **Menu > Instrument > Set up Instrument Method** or via right click into the module's active area (does not show the **Instrument Curves** tab).

Setup Method

VWD Instrument Curves

VWD (G7114B)

Signal

☐ Enable Dual-Wavelength

Acquire Wavelength

Signal A: ☒ nm

Peakwidth:

Stoptime **Posttime**

☒ As Pump/Injector ☐ Off

☐ min ☐ min

Advanced

Analog Output

Zero Offset: %

Attenuation: mAU

Signal Polarity **Autobalance**

☒ Positive (+) ☒ Prerun

☐ Negative (-) ☐ Postrun

Miscellaneous

☒ Lamp on required for acquisition

Scan Range: to nm

Step: nm

Timetable (empty)

☐ Show timetable graph

OK Apply Cancel Help

Figure 12: Method parameter settings

NOTE

For additional help and support: Highlight the desired cell and press **F1**. A help screen will open with additional information and documentation about the topic.

Table 12: Method Parameter Settings

Signal

☐ Enable Dual-Wavelength

Acquire

Wavelength

Signal A: ☒ nm

Peakwidth:

> 0.1 min (2 s resp. time) (5 Hz)

< 0.00078 min (0.0078 s resp. time) (240 Hz)

> 0.00078 min (0.016 s resp. time) (240 Hz)

> 0.0016 min (0.031 s resp. time) (240 Hz)

> 0.0031 min (0.063 s resp. time) (160 Hz)

> 0.0063 min (0.13 s resp. time) (80 Hz)

> 0.013 min (0.25 s resp. time) (40 Hz)

> 0.025 min (0.5 s resp. time) (20 Hz)

> 0.05 min (1 s resp. time) (10 Hz)

> 0.1 min (2 s resp. time) (5 Hz)

> 0.2 min (4 s resp. time) (2.5 Hz)

> 0.4 min (8 s resp. time) (1.25 Hz)

Stoptime

☒ As Pump/Inje

☐

min

Signal

Wavelength

Single Wavelength (190 – 600 nm, step 1)

Dual Wavelength Mode enables the multi-wavelength mode with two wavelengths.

Peakwidth (Responsetime, Data Rate)

Peakwidth enables you to select the peak width (response time) for your analysis. The peak width is defined as the width of a peak, in minutes, at half the peak height. Set the peak width to the narrowest expected peak in your chromatogram. The peak width sets the optimum response time for your detector. The peak detector ignores any peaks that are considerably narrower, or wider, than the peak width setting. The response time is the time between 10 % and 90 % of the output signal in response to an input step function. When the All spectrum storage option is selected, then spectra are acquired continuously depending on the setting of the peak width. The time specified by the peak width is used as a factor in the acquisition of spectra. The acquisition time for one spectrum is slightly less than the peak width divided by 8 , which is the acquisition time.

Limits: When you set the peak width (in minutes), the corresponding response time is set automatically and the appropriate data rate for signal and spectra acquisition is selected.

NOTE: The G7114A VWD has a data rate of up to 120 Hz. The G7114B VWD has a data rate of up to 240 Hz.

G7114B Peakwidth settings up to 240 Hz

Signal

☐ Enable Dual-Wavelength

Acquire

Wavelength

Signal A: ☒ nm

Peakwidth:

> 0.1 min (2 s resp. time) (5 Hz)

< 0.0016 min (0.016 s resp. time) (120 Hz)

> 0.0016 min (0.031 s resp. time) (120 Hz)

> 0.0031 min (0.063 s resp. time) (120 Hz)

> 0.0063 min (0.13 s resp. time) (80 Hz)

> 0.013 min (0.25 s resp. time) (40 Hz)

> 0.025 min (0.5 s resp. time) (20 Hz)

> 0.05 min (1 s resp. time) (10 Hz)

> 0.1 min (2 s resp. time) (5 Hz)

> 0.2 min (4 s resp. time) (2.5 Hz)

> 0.4 min (8 s resp. time) (1.25 Hz)

Stoptime

☒ As Pump/Inje

☐

min

G7114A Peakwidth settings up to 120 Hz

Dual-Wavelength Settings

☒ Enable Dual-Wavelength

Wavelength

Signal A: nm

Signal B: nm

Peakwidth:

< 0.4 min (4 s resp. time) (2.5 Hz)

G7114B Dual Wavelength Settings

Stoptime	Posttime
<input checked="" type="radio"/> As Pump/Injector <input type="radio"/> 1.00 min	<input checked="" type="radio"/> Off <input type="radio"/> 1.00 min

Stoptime/Posttime

The stoptime is the time where either the complete system stops (**As Pump/Injector**) or the module (if different from system stop time). The data collection is stopped at this time. A posttime period can be used to allow module's items to equilibrate (e.g. after gradient change or temperature change).

Analog Output
Zero Offset: 5 % Attenuation: 1000 mAU

Analog Output

The range can be set to either 100 mV or 1 V full scale, see [Table](#) on page 68.

- **Zero Offset:** 1 – 99 % in steps of 1 % (5 % equal to 50 mV).
- **Attenuation:** 0.98 – 4000 mAU at discrete values for either 100 mV or 1 V full scale.

Signal Polarity	Autobalance
<input checked="" type="radio"/> Positive (+) <input type="radio"/> Negative (-)	<input checked="" type="checkbox"/> Prerun <input type="checkbox"/> Postrun

Signal Polarity

Can be switched to negative (if required).

Autobalance

Defines, whether a balance is performed prior to a run and/or after a run has finished.

Miscellaneous
<input checked="" type="checkbox"/> Lamp on required for acquisition Scan Range: 190 to 400 nm Step: 2 nm

Miscellaneous

Lamp on required for acquisition: If unchecked, the lamp will be turned off after the analysis has finished. Note that the lamp on requires at least one hour warm-up time, see [Warm up of the Detector](#) on page 82.

Scan Range / Step: Stop-Flow scan range / step. Access to the scan feature is only possible during run. See [Scanning with the VWD](#) on page 72.

▲ Timetable (1/100 events)

Time [min]	Function	Parameter
▶ 0.00	Balance	Autobalancing
	Balance	
	Change Signal	
	Take Scan	

Timetable

You may set up time events to change functions with their parameters over the run time. Add lines as required.

Time Limits: 0.00 to 99999.00 min in steps of 0.01 min.

Via the buttons in the bottom area, time table lines can be added, removed, cut copied, pasted or completely cleared.

Based on the chosen function, a certain parameter can be selected.

Add	Remove	Clear all
Cut	Copy	Paste

Setup Method

WWD

Instrument Curves

VWD:

☐ Signal without Reference (mAU)

☐ Reference only (mAU)

☐ Signal B without Reference (mAU)

☐ Reference B only (mAU)

☐ Board Temperature (°C)

☐ UV Lamp Anode Voltage (V)

☐ Sample Intensity (counts)

☐ Reference Intensity (counts)

☐ Optical Unit Temperature (°C)

☐ Show timetable graph

OK

Apply

Cancel

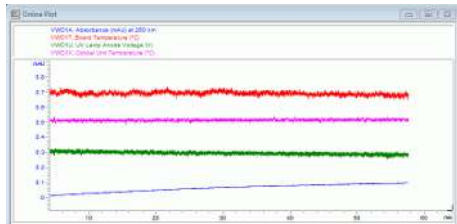
Help

Instrument Curves

The detector has several signals (internal temperatures, voltages of lamps) that can be used for diagnosing problems. These can be baseline problems deriving from deuterium lamps wander / drift problems due to temperature changes.

These signals can be used in addition to the normal baseline signal to determine whether correlation to temperature or voltage/current of the lamp.

These signals are available via the Agilent ChemStation Online Plot/Data Signal and/or Agilent Lab Advisor Software.



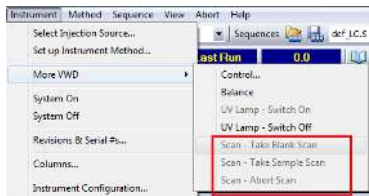
Scanning with the VWD

NOTE

Access to the scan feature is only possible during run with stopped flow. The spectrum is taken during a stop-flow condition only while the peak is kept in the flow cell.

- 1 Set up a run.
- 2 Start a run.
- 3 While running on the baseline, take a **Blank Scan**. A background scan is stored in the memory.

Table 13: Blank scan



- Step 1: Blank Scan: scan of the background (solvent) is stored in the memory.
- Step 2: Sample Scan: scan of the peak of interest is taken while the peak stays in the flow cell (stop-flow condition).
- Online Spectrum: Sample Scan minus Blank Scan.

Here the functions are inactive (grayed out). Will be active in run mode.

- 4 When the peak of interest enters the flow cell, stop the flow (set flow rate to zero or open the purge valve) and wait a few moments to stabilize the concentration.

NOTE

Turning off the pump would stop the run and no access to the sample scan is possible.

Using the Module

Preparing the Module

- 5 Open the Online Spectra window (**View > Online Spectra > VWD**) and change the absorbance and wavelength range according your needs.

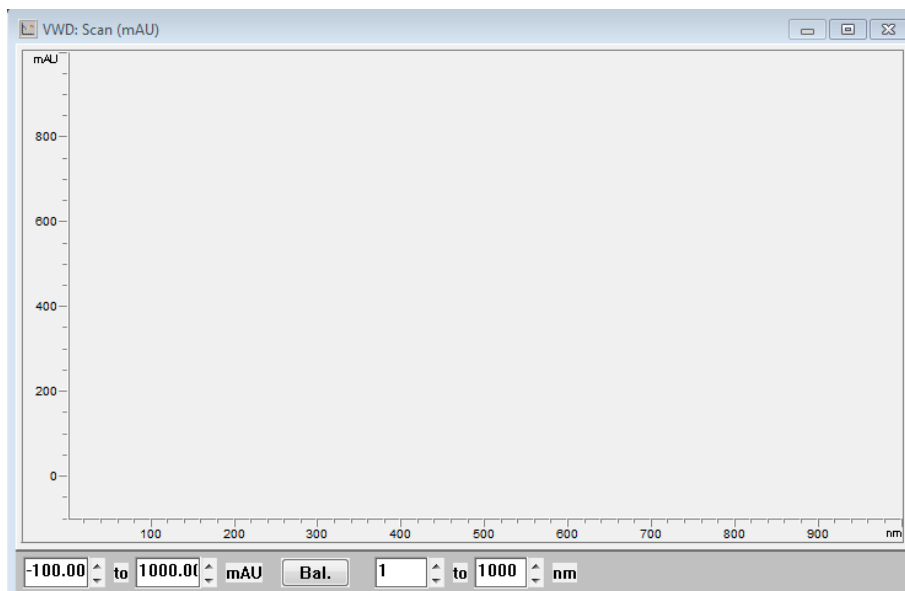


Figure 13: Online Spectra Window

- 6 Select **Sample Scan**. A sample scan is taken in the range defined under **Miscellaneous** in [Table 12](#) on page 70 and displays the result (Sample Scan minus Blank Scan).



5

Optimizing the Performance of the Module

This chapter provides information on how to optimize the module.

Introduction 76

Match the Flow Cell to the Column 77

Set the Detector Parameters 81

Warm up of the Detector 82

Introduction

The detector has a variety of parameters that can be used to optimize performance.

The information below will guide you on how to get the best detector performance. Follow these rules as a start for new applications. It gives a rule-of-thumb for optimizing the detector parameters.

Match the Flow Cell to the Column

The tables below recommend the flow cell that matches 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.

Standard HPLC Applications

Column length	Typical peak width	Recommended flow cell				
<= 5 cm	0.025 min	Micro flow cell				High Pressure flow cell
10 cm	0.05 min		Semimicro flow cell			
20 cm	0.1 min			Standard flow cell		
>= 40 cm	0.2 min					
	Typical flow rate	0.05-0.2 ml/min	0.2- 0.4 ml/min	0.4- 0.8 ml/min	1-2 ml/min	0.01- 5 ml/min
	Internal column diameter	1.0 mm	2.1mm	3.0 mm	4.6 mm	

Figure 14: Choosing a Flow Cell (Standard HPLC Applications)

Ultra Fast Separation With RRLC Systems

Column ID	2.1 mm	3.0 mm	4.6 mm
Configuration	No damper No mixer ++	Damper Mixer	Damper Mixer
Flow cell	2 µl, 3 mm	5 µl, 6 mm +	14 µl, 10 mm +

Figure 15: Choosing a Flow Cell for G7114B (for ultra fast separation with RRLC systems)

- (+) For ultra fast analysis with step gradients the micro flow cell (2 µL, 3 mm) gives the best performance
- (++) In high resolution analysis time is not the highest priority. Higher delay volumes are accepted. Therefore we recommend to use the damper plus mixer for a highest signal to noise.
- If longer columns (> 50 mm) for higher resolution are used, then the next larger flow cell is the preferred choice for higher sensitivity.

Flow Cell Path Length

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

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

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 [mol/L]	is the concentration of the absorbing species,
d [m]	is the path length of the cell used for the measurement.

Therefore, flow cells with longer path lengths yield higher signals. Although noise usually increases little with increasing path length, there is a gain in the signal-to-noise ratio. For example, in [Figure 16](#) on page 79 the noise increased by less than 10 % but a 70 % increase in signal intensity was achieved by increasing the path length from 6 mm to 10 mm.

When increasing the path length, the cell volume usually increases – in the example from 5 – 14 µL. Typically, this causes more peak dispersion. As demonstrated, this did not affect the resolution in the gradient separation in the example that is shown below.

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).

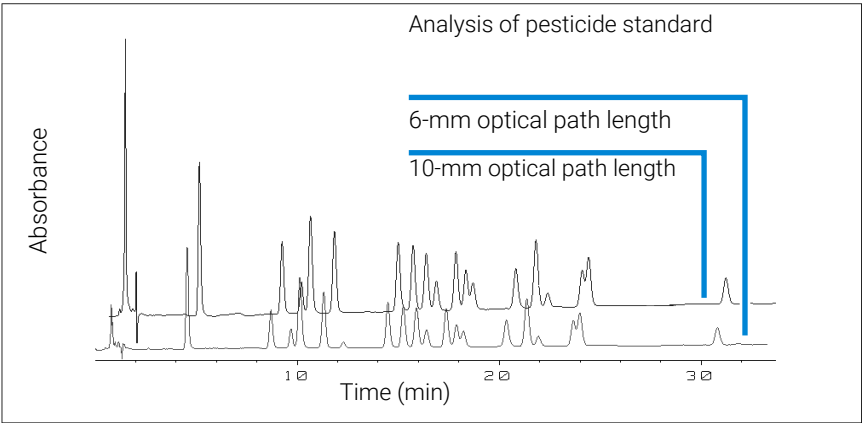


Figure 16: Influence of Cell Path Length on Signal Height

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

Please find below the details of the flow cells:

Table 14: Correction factors for Agilent VWD flow cells

Part number	Path length (actual)	Correction factor
G1314-60186 (Standard flow cell 10 mm, 14 µL, 40 bar)	10.15 ± 0.19 mm	10/10.15
G1314-60183 (Semi-micro flow cell 6 mm, 5 µL)	6.10 ± 0.19 mm	6/6.10
G1314-60187 (Micro flow cell 3 mm, 2 µL, 120 bar)	2.80 ± 0.19 mm	3/2.8
G1314-60182 (High pressure flow cell 10 mm, 14 µL, 400 bar)	10.00 ± 0.19 mm	10/10
G1314-60188 (Bio standard flow cell VWD, 10 mm, Cell Vol. 14 µl, Sapphire, MP35N)	10.15 ± 0.19 mm	10/10.15
G1314-60189 (Bio micro flow cell VWD, 3 mm, Cell Vol. 2 µl, Sapphire, MP35N)	2.80 ± 0.19 mm	3/2.8

NOTE

However you have to be aware that there is additional tolerance of gasket thickness and its compression ratio which is supposed to be very small in comparison with the machining tolerance.

Set the Detector Parameters

- 1 Set peakwidth as close as possible to the width (at half height) of a narrow peak of interest.
- 2 Choose the sample wavelength.
 - at a longer wavelength than the cut-off wavelength of the mobile phase,
 - at a wavelength where the analytes have strong absorptivity if you want to get the lowest possible detection limit,
 - at a wavelength with moderate absorptivity if you work with high concentrations, and
 - preferably where the spectrum is flat for better linearity.
- 3 Consider to use time-programming to further optimization.

Warm up of the Detector

Give the optical unit enough time to warm-up and stabilize (> 60 minutes). The detector is temperature controlled. After turn-on of the detector, it goes through a cycle of different states:

- 0 to 0.5 minutes the heater control is OFF and the heater element runs at 0 % duty cycle.
- 0.5 to 1 minutes the heater control is OFF and the heater element runs at 66% duty cycle. This first minute is used as self-test of the heater functionality.
- 1 to 30 minutes the heater control is OFF and the heater element runs at 40% duty cycle.
- After 30 minutes the heater control is ON and is working with optimized parameters to get the optical unit into the optimal temperature window stabilized.

This cycle starts

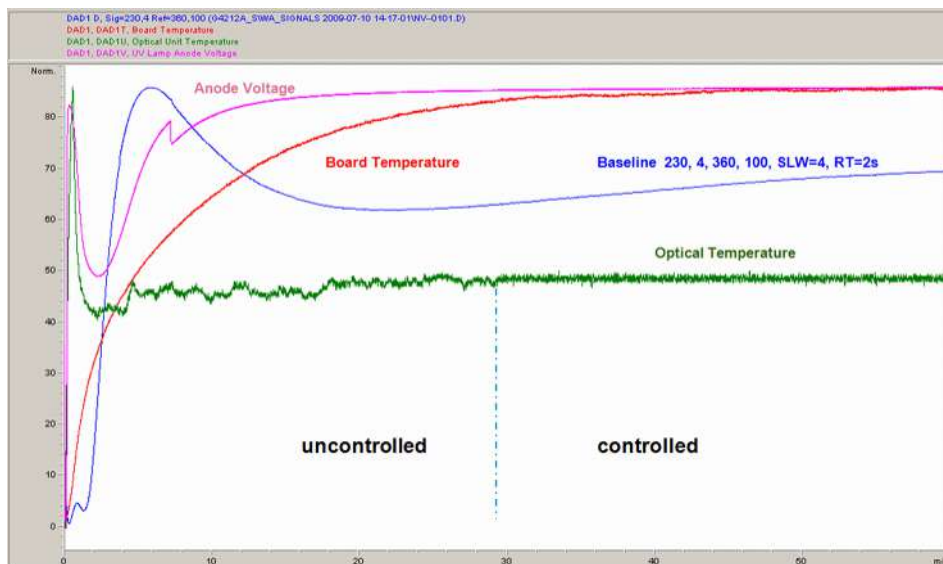
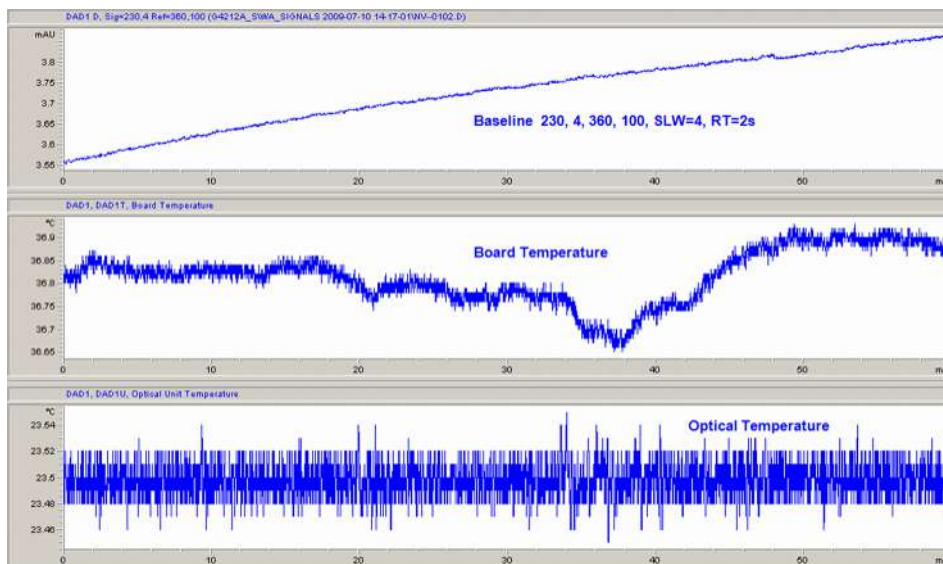
- when the detector is turned off/on
- when the lamp is turned off/on

to ensure that the temperature control operates in a defined control range.

NOTE

The times to stabilize the baseline may vary from instrument to instrument and depends on the environment. The example below was done under stable environmental conditions.

The figures below show the first two hours of a detector warm-up phase. The lamp was turned on immediately after turn on of the detector.

Figure 17: Detector Warm-up – 1st hourFigure 18: Detector Warm-up – 2nd hour

6

Diagnostics and Troubleshooting

This chapter gives an overview of the maintenance, troubleshooting, and diagnostic features available.

Diagnostic Features 85

User Interfaces 85

Troubleshooting With HPLC Advisor 85

Overview of Available Tests and Tools 86

Tests and Calibrations in Agilent Lab Advisor 86

Maintenance and Troubleshooting Tools of the Module 87

Introduction 87

Available Tests versus Interfaces 90

Intensity Test 91

Cell Test 93

Wavelength Calibration 95

ASTM Drift and Noise Test 97

Quick Noise Test 98

Dark Current Test 100

Holmium Oxide Test 103

D/A Converter Test 105

Using the Built-In Test Chromatogram 108

Agilent Lab Advisor Software 111

Other Lab Advisor Functions 112

EMFs - Early Maintenance Feature 112

Diagnostic Features

This section gives an overview of the diagnostic features available.

User Interfaces



InfinityLab Assist

InfinityLab Assist provides you with assisted troubleshooting and maintenance at your instrument.

If the system in use supports the InfinityLab Assist, follow the instructions provided. Else, the preferred solution is to use Agilent Lab Advisor Software.

- Depending on the user interface, the available tests and the screens/reports may vary.
- The preferred tool for troubleshooting and diagnostics should be Agilent Lab Advisor Software, see [Agilent Lab Advisor Software](#) on page 111.
- Screenshots used within these procedures are based on the Agilent Lab Advisor Software.

Troubleshooting With HPLC Advisor

Baseline, Peak Shape, Pressure, Retention related issues, can be solved using the HPLC Advisor App. For more information, see Troubleshooting Reversed-Phase Chromatographic Techniques With HPLC Advisor.

If using an InfinityLab Assist, navigate to **Health > Troubleshooting** to help solve baseline, peak shape, pressure, and retention related issues.

Overview of Available Tests and Tools

Tests and Calibrations in Agilent Lab Advisor

Use the tests and diagnostic features provided in the Agilent Lab Advisor software to check if your module is working correctly.

For further details, refer to the Agilent Lab Advisor software help files.

Maintenance and Troubleshooting Tools of the Module

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

Introduction

All tests are described based on the Agilent Lab Advisor Software B.02.08. Other user interfaces may not provide any test or just a few.

Table 15: Interfaces and available test functions

Interface	Comment	Available Function
Agilent Lab Advisor	For functions, refer to Function Overview Lab Advisor • Table 16 on page 87	Available functions depend on Product Level (Basic – Advanced – FSE)
Agilent ChemStation	No tests available Adding of temperature/lamp signals to chromatographic signals possible	<ul style="list-style-type: none">• Temperature main board• Temperature optical unit• Lamp anode voltage

Table 16: Function Overview Lab Advisor Basic/Advanced (G7114A/G7114B)

Function	Product Level	
Tests		
- ASTM Drift and Noise Test	Basic	Advanced
- Cell Test	Basic	Advanced
- D/A Converter Test	Basic	Advanced
- Dark Current Test	Basic	Advanced
- Filter/Grating Motor Test	Basic	Advanced
- Holmium Oxide Test	Basic	Advanced
- Intensity Test	Basic	Advanced

Function	Product Level	
- Quick Noise Test	Basic	Advanced
Calibrations		
- Wavelength Calibration	Basic	Advanced
Tools		
- Diagnostic Buffers	Basic	Advanced
- Board Check and Change		
- Module Info	Basic	Advanced
- Firmware Declustering		
- Test Chromatogram	Basic	Advanced
- Spectral Scan	Basic	Advanced
Controls		
- Advanced Method Parameters		
- Analog Output 1 Attenuation		Advanced
- D2 lamp required		Advanced
- Analog Output 1 Offset [% Full Scale]		Advanced
- Configuration		
- Remote Pulse Duration [s] *	Basic	Advanced
- Analog Output 1 Range		Advanced
- Control		
- Balance Detector		Advanced
- UV Lamp	Basic	Advanced
- Conversions		
- Transfer to Resident FW		
- Transfer to Main FW *		
- G7114B allows G1314E and G1314F	Basic	Advanced
- G7114A allows NONE	Basic	Advanced
- Method Parameters		
- Set Data Rate [Hz]		Advanced
- Set Wavelength [nm]		Advanced
- Module Information		

Function	Product Level	
- Firmware Version		
- Identify Module	Basic	Advanced
- Special Commands		
- Lamp tag required	Basic	Advanced
- Cell tag required	Basic	Advanced
- Clear Error	Basic	Advanced
- Detector Reset	Basic	Advanced
- Forced Cold Start		
Actuals		
- Signal A [mAU]		Advanced
- Sample Signal		
- Reference Signal		
Statemachines		
- UV Lamp	Basic	Advanced
Signals		
- Signal A [mAU]		Advanced
- Sample Signal [mAU]		Advanced
- Reference Signal [mAU]		Advanced
- Board Temperature [°C]		Advanced
- Lamp Voltage [V]		Advanced
EMF Counters		
- Accumulated UV Lamp On-Time	Basic	Advanced
- Number of UV Lamp Ignitions	Basic	Advanced

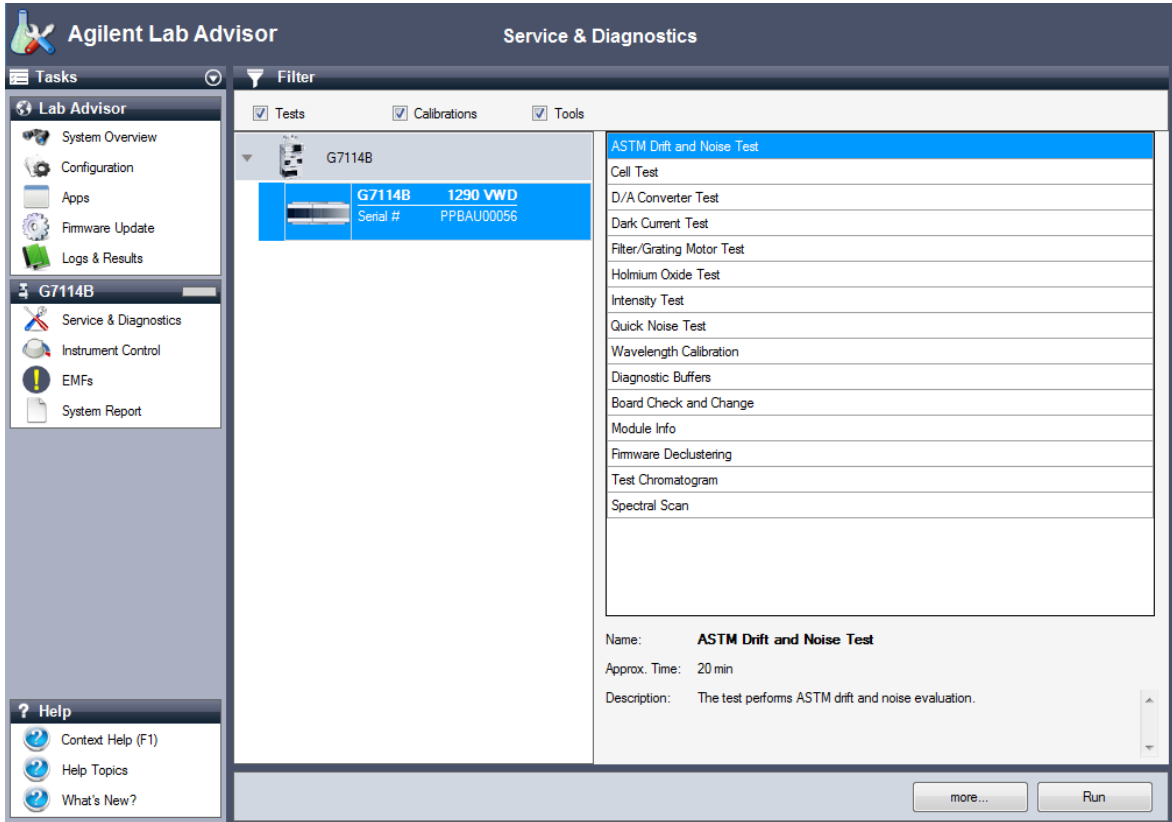


Figure 19: The Lab Advisor shows the available test

Available Tests versus Interfaces

NOTE

Depending on the used interface, the available tests and the screens/reports may vary.

Preferred tool should be the Agilent Lab Advisor, see [Agilent Lab Advisor Software](#) on page 111.

Agilent Lab Advisor B.02.08 or later is required.

The Instant Pilot (G4208A) supports the G7114B with B.02.19 and the G7114A with B.02.20 or later.

Diagnostics and Troubleshooting

Maintenance and Troubleshooting Tools of the Module

- Preferred tool should be the Agilent Lab Advisor software, see [Agilent Lab Advisor Software](#) on page 111
- Screenshots used within these procedures are based on the Agilent Lab Advisor software.

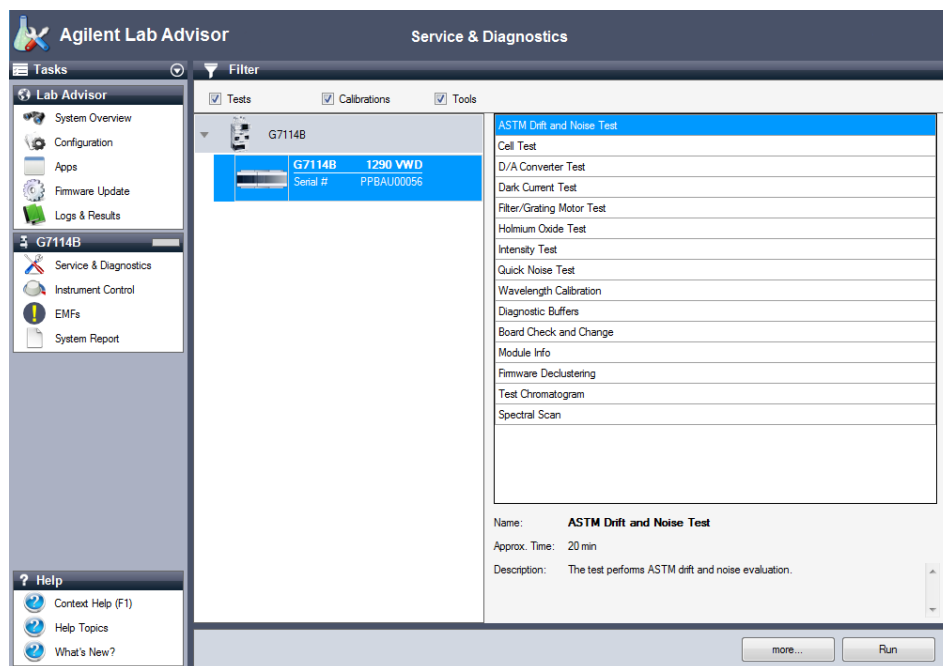


Figure 20: The Lab Advisor shows the available test

Intensity Test

The **Intensity Test** measures the intensity of the deuterium lamp over the full VWD wavelength range (190 - 600 nm). The test can be used to determine the performance of the lamp, and to check for dirty or contaminated flow cell windows. When the test is started, 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 characteristics. Therefore, intensity spectra will differ slightly between instruments. The figure below shows a typical intensity test spectrum.

The Intensity Test is available in Agilent Lab Advisor (preferred tool).

Intensity Test evaluation

The Agilent Lab Advisor and the Instant Pilot evaluate three values automatically and display the limits for each value, the average, the minimum and the maximum of all data points and **passed** or **failed** for each value.

- 1 Run the Intensity-Test with Agilent Lab Advisor (for further information see Online-Help of user interface).

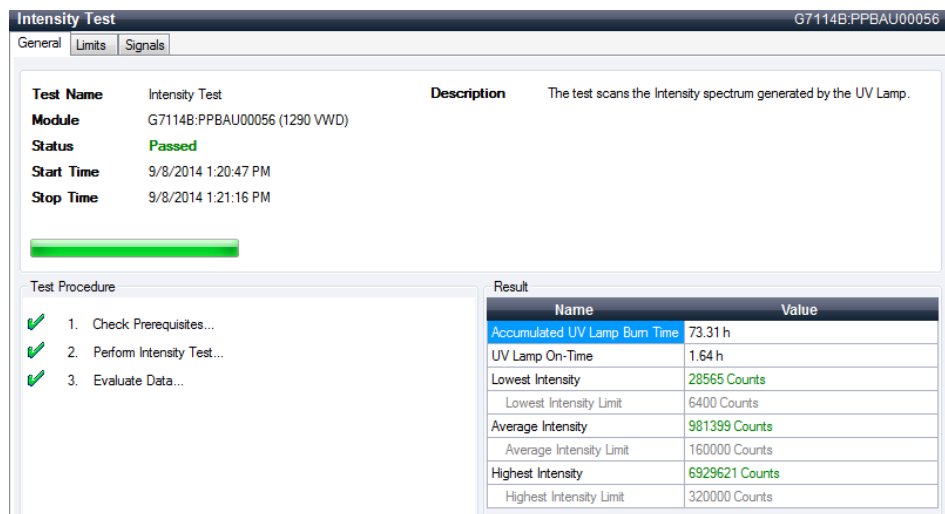


Figure 21: Intensity Test – Results

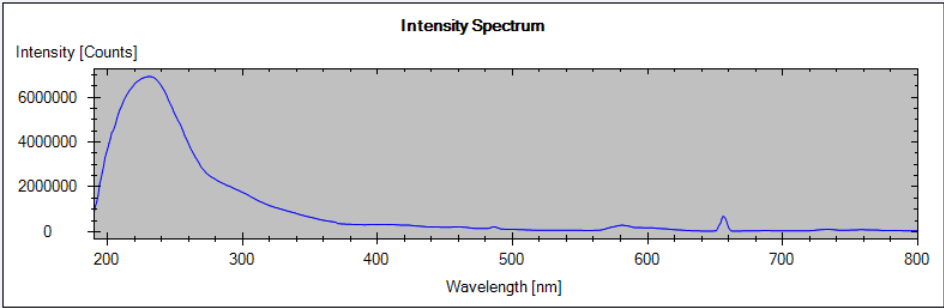


Figure 22: Intensity Test – Signal

Intensity Test Failed

Probable cause		Suggested actions
1	Empty flow cell	• Ensure the flow cell is filled with water.
2	Flow cell windows dirty	• Repeat the test with the flow cell window removed. If the test passes, exchange the flow cell windows.
3	Optics defect	• Please contact your Agilent service representative.
4	Defective lamp or optics.	• Exchange the lamp.

Cell Test

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.

Limits: No real limit. The reason is that it depends on the position/alignment of the reference side (beam splitter – reference slit – reference diode). Therefore the reference side value can be higher/smaller than the sample side value.

With a clean cell the counts for sample and reference (photocurrent) are in the same range. If the sample side shows much lower values than the reference side the flow cell might have a problem.

Preparations

- Flush the flow cell with a flow of 1 mL/min for at least 10 minutes.
- 1 Run the Cell-Test with Agilent Lab Advisor (for further information see Online-Help of user interface).

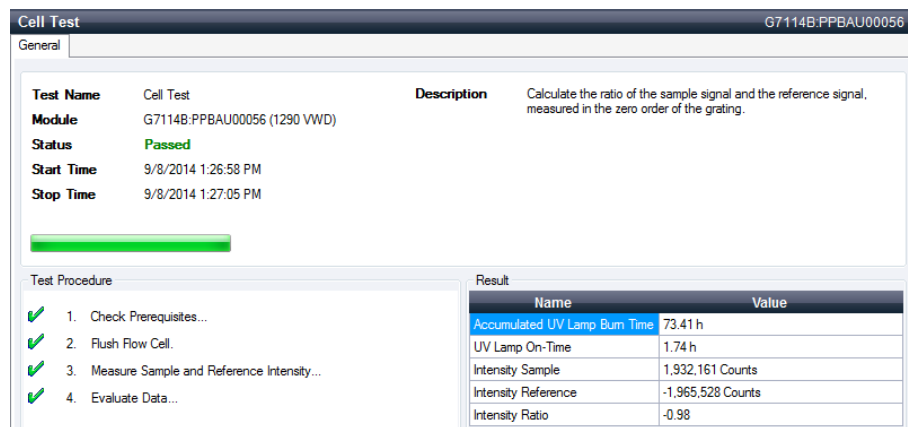


Figure 23: Test – Results

Cell test failed

Probable cause		Suggested actions
1	Cell contaminated	<ul style="list-style-type: none">• Flush flow cell
2	Cell windows are contaminated	<ul style="list-style-type: none">• Clean/replace cell windows
3	Mechanical problem	<ul style="list-style-type: none">• Check cell position

Wavelength Calibration

Wavelength Calibration of the detector is done using the zero-order position and 656 nm emission line position of the deuterium lamp. The calibration procedure involves two 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.

In addition to the zero-order and 656 nm (alpha-emission line) calibration, the beta-emission line at 486 nm and the three holmium lines are used for the complete wavelength calibration process. These holmium lines are at 360.8 nm, 418.5 nm and 536.4 nm.

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.

The Wavelength Verification/Calibration is available in Agilent Lab Advisor (preferred tool).

When

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.

NOTE

If the detector was repaired (opened covers), the wavelength calibration can be done 10 minutes after lamp on. A final wavelength calibration should be repeated after complete warm-up of the detector.

- 1 Run the Wavelength Calibration with the Agilent Lab Advisor (for further information see Online-Help of user interface).

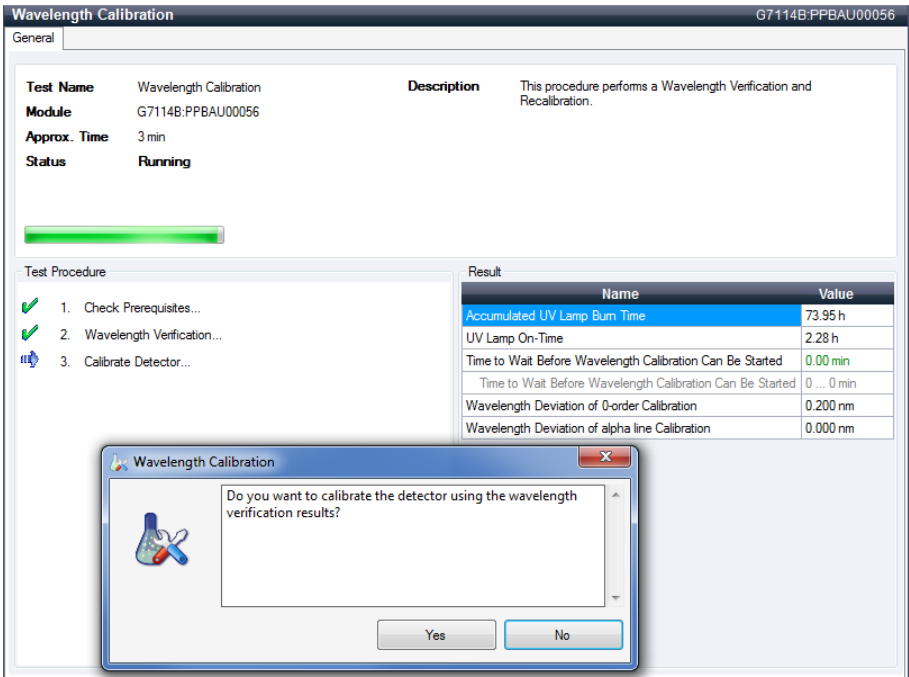


Figure 24: Wavelength Calibration - Results

If you select **No**, the test is aborted.
If you select **Yes**, the re-calibration is performed (the offset is corrected).

ASTM Drift and Noise Test

The **ASTM Drift and Noise Test** determines the detector noise over a period of 20 minutes. The test is done with HPLC-grade water flowing through the flow cell at 1 mL/min. On completion of the test, the noise result is displayed automatically.

- 1 Run the ASTM Drift and Noise Test with Agilent Lab Advisor (for further information see Online- Help of user interface).

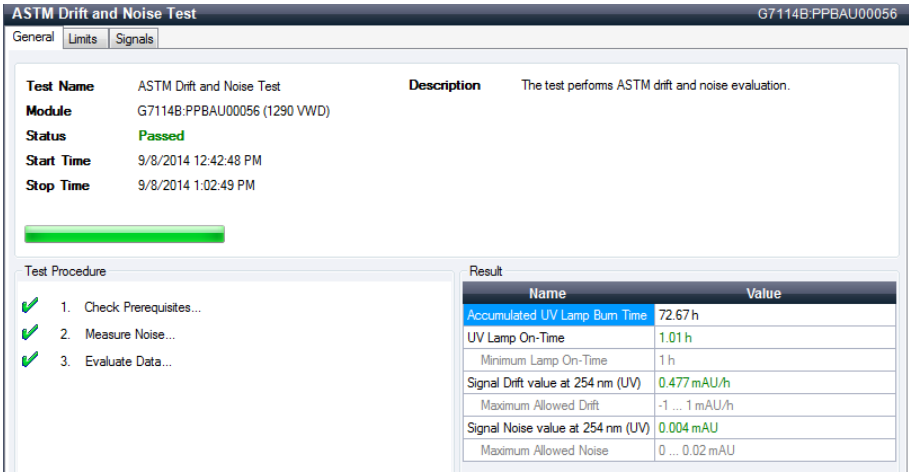


Figure 25: ASTM Drift and Noise Test – Results

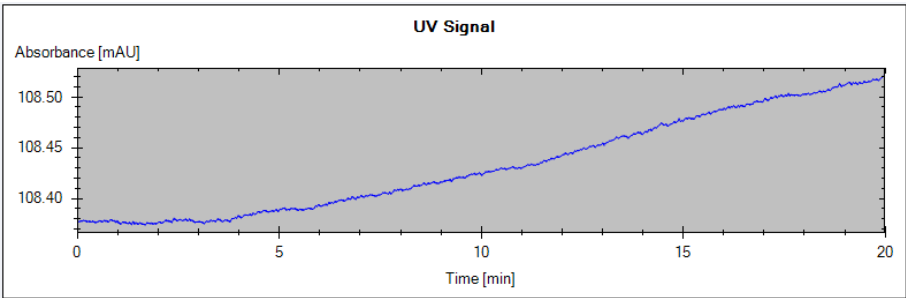


Figure 26: Drift and Noise Test – Signal

Quick Noise Test

The noise test measures the noise of the detector, with HPLC-grade water flowing through the flow cell at 1 mL/min, in one minute intervals over a total of 5 minutes.

The noise of the detector is calculated by using the maximum amplitude for all random variations of the detector signal of frequencies greater than one cycle per hour. The noise is determined for 5 one minute intervals and is based on the accumulated peak-to-peak noise for the intervals. At least seven data points per cycles are used in the calculation.

The cycles in the noise determination are not overlapping.

In order to obtain reliable results, the lamp should be turned on for at least 10 minutes prior to measurement.

- 1 Run the Quick Noise Test with Agilent Lab Advisor (for further information see Online-Help of user interface).

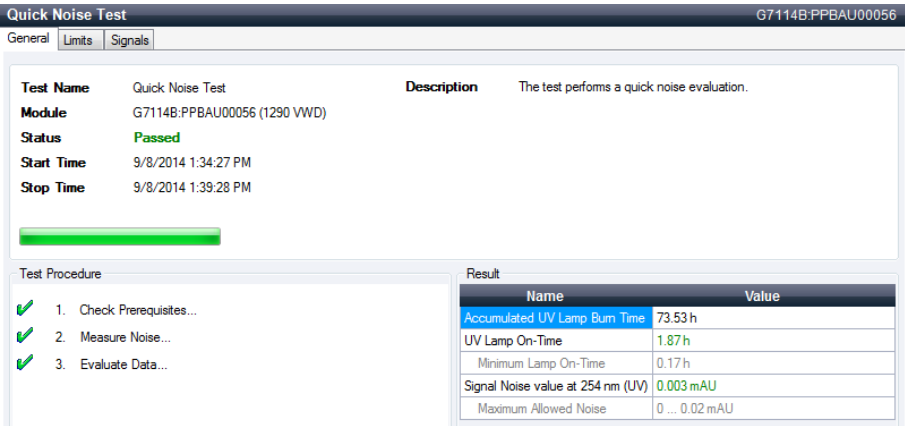


Figure 27: Quick Noise Test – Results

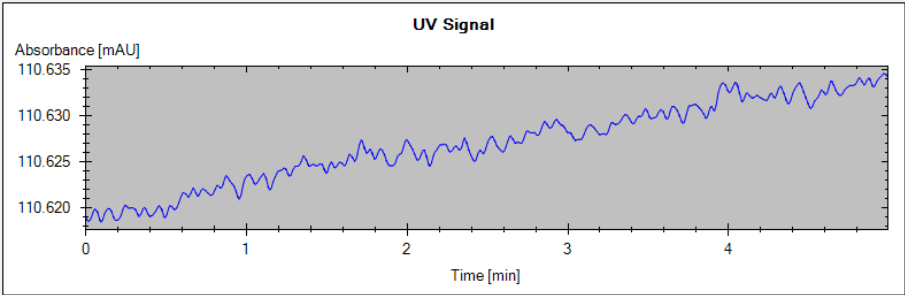


Figure 28: Quick Noise Test – Signal

Dark Current Test

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 shutter is moved into the light path. Next, the leakage current from both diodes is measured.

- 1 Run the Dark Current Test with the Agilent Lab Advisor (for further information see Online-Help of user interface).

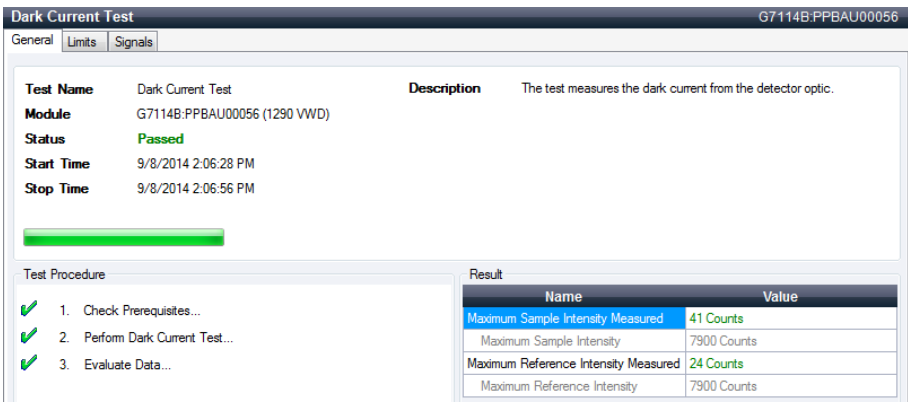


Figure 29: Dark Current Test - Results

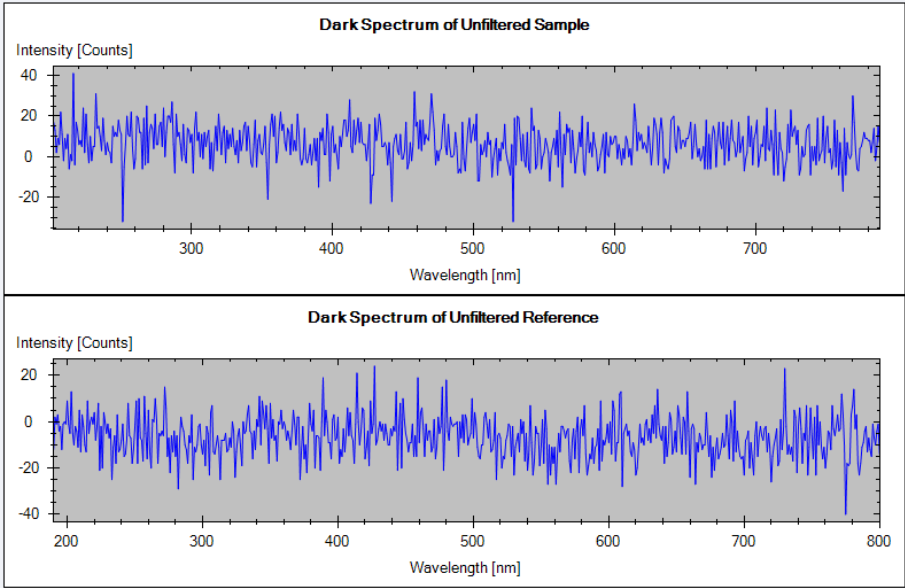


Figure 30: Dark Current Test - Signals

Dark Current Test Failed

Probable cause		Suggested actions
1	Defective sample or reference diode.	• Please contact your Agilent service representative.
2	Defective sample or reference ADC board.	• Please contact your Agilent service representative.
3	Defective mainboard.	• Please contact your Agilent service representative.

Holmium Oxide Test

This **Holmium Oxide Test** verifies the calibration of the detector against the three wavelength maxima of the built-in holmium oxide filter. The test displays the difference between the expected and measured maxima. The figure below shows a holmium test spectrum.

The Holmium Oxide Test is available in Agilent Lab Advisor (preferred tool).

The test uses the following holmium maxima:

- 360.8 nm
- 418.5 nm
- 536.4 nm

NOTE

See also [Declaration of Conformity for HOX2 Filter](#) on page 263.

When

- after recalibration,
- as part of the Operational Qualification/Performance Verification procedure, or
- after flow cell maintenance or repair.

Interpreting the results

The test is passed successfully when all three wavelengths are within ± 1 nm of the expected value. This indicates the detector is calibrated correctly.

- 1 Run the Holmium Oxide Test with the Agilent Lab Advisor (for further information see Online- Help of user interface).

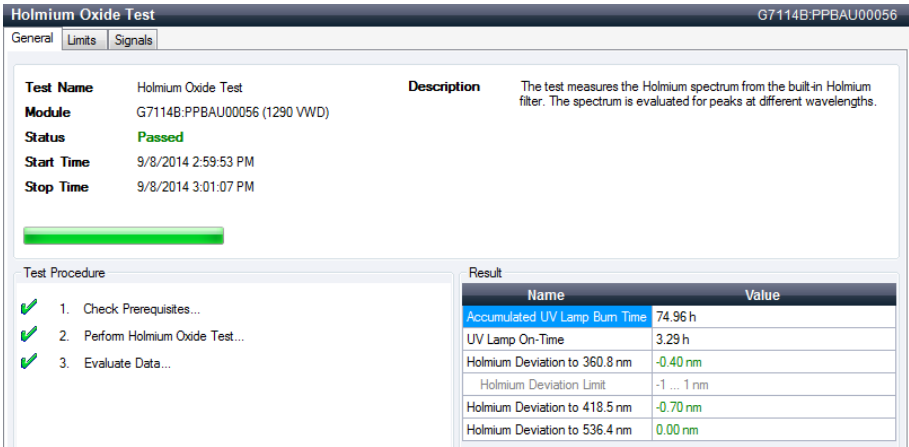


Figure 31: Holmium Oxide Test - Results

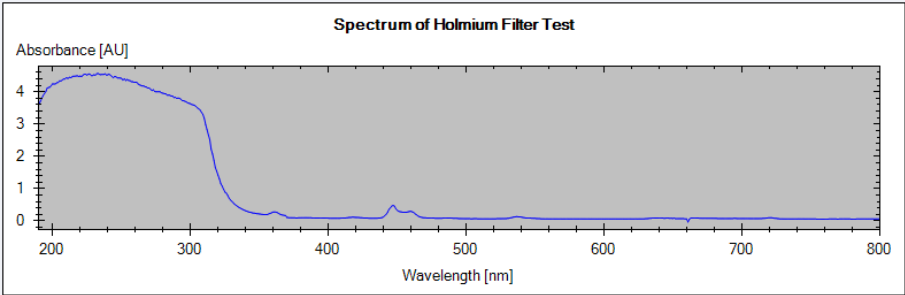


Figure 32: Holmium Oxide Test - Signals

Holmium Oxide Test Failed

Probable cause		Suggested actions
1	Detector not calibrated.	<ul style="list-style-type: none">Recalibrate the detector.
2	Dirty or defective flow cell.	<ul style="list-style-type: none">Repeat the test with the flow cell removed. If the test is OK, exchange the flow cell components.
3	Dirty or defective holmium oxide filter.	<ul style="list-style-type: none">Run the holmium oxide filter test. If the test fails, contact your Agilent service representative.
4	Optical misalignment.	<ul style="list-style-type: none">Please contact your Agilent service representative.

D/A Converter 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 **D/A Converter 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 μ V 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 Lab Advisor

- 1 Run the D/A Converter Test with the Agilent Lab Advisor (for further information see Online- Help of user interface).

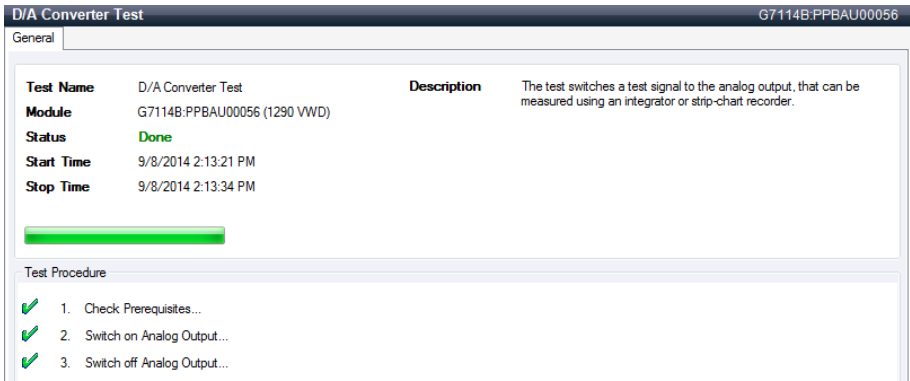


Figure 33: Converter Test - Results

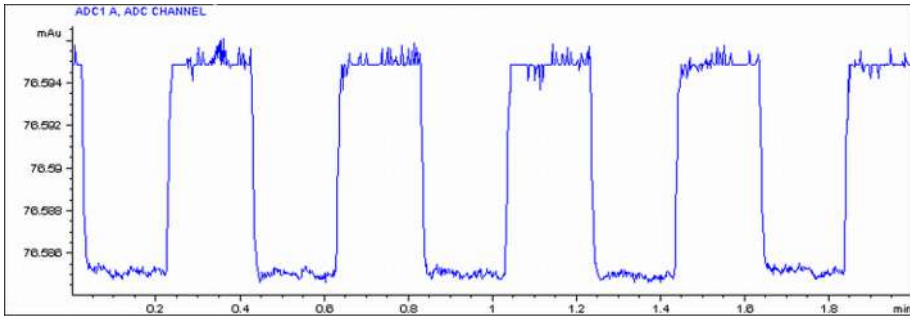


Figure 34: D/A Converter (DAC) Test – Example of Integrator Plot

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`

D/A Converter Test failed

D/A Converter Test evaluation

The noise on the step should be less than 3 μ V.

Probable cause		Suggested actions
1	Bad cable or grounding problem between detector and external device.	<ul style="list-style-type: none">• Check or replace the cable.
2	Defective detector main board.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

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.

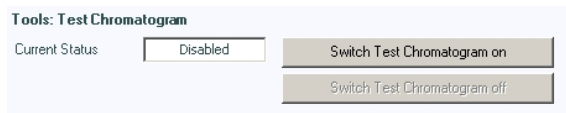
NOTE

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

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

Procedure using the Agilent Lab Advisor

- 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.

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

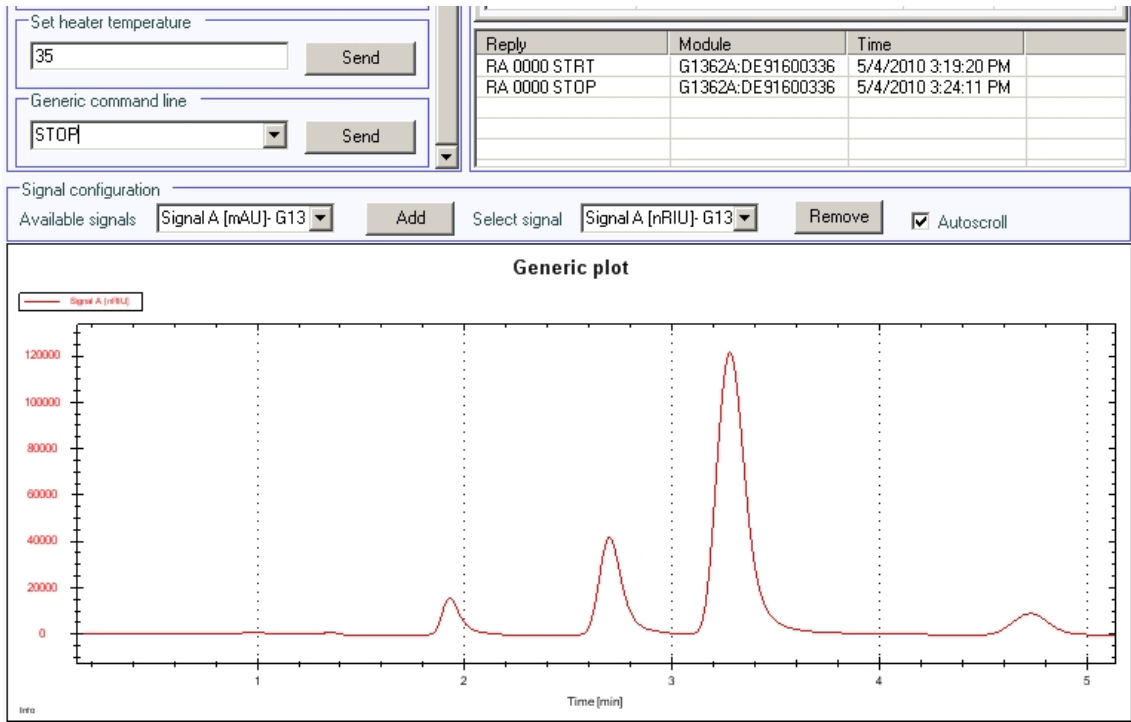


Figure 35: Test Chromatogram with Agilent Lab Advisor

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.

Agilent Lab Advisor Software

The Agilent Lab Advisor Software (basic license, shipped with an Agilent LC pump) is a standalone product that can be used with or without a chromatographic data system. Agilent Lab Advisor helps to manage the lab for high-quality chromatographic results by providing a detailed system overview of all connected analytical instruments with instrument status, Early Maintenance Feedback counters (EMF), instrument configuration information, and diagnostic tests. With the push of a button, a detailed diagnostic report can be generated. Upon request, the user can send this report to Agilent for a significantly improved troubleshooting and repair process.

The Agilent Lab Advisor software is available in two versions:

- Lab Advisor Basic
- Lab Advisor Advanced

Lab Advisor Basic is included with every Agilent 1200 Infinity Series and Agilent InfinityLab LC Series instrument.

The Lab Advisor Advanced features can be unlocked by purchasing a license key, and include real-time monitoring of instrument actuals, all various instrument signals, and state machines. In addition, all diagnostic test results, calibration results, and acquired signal data can be uploaded to a shared network folder. The Review Client included in Lab Advisor Advanced makes it possible to load and examine the uploaded data no matter on which instrument it was generated. This makes Data Sharing an ideal tool for internal support groups and users who want to track the instrument history of their analytical systems.

The optional Agilent Maintenance Wizard Add-on provides an easy-to-use, step-by-step multimedia guide for performing preventive maintenance on Agilent 1200 Infinity LC Series instrument.

The tests and diagnostic features that are provided by the Agilent Lab Advisor software may differ from the descriptions in this manual. For details, refer to the Agilent Lab Advisor software help files.

Other Lab Advisor Functions

EMFs - Early Maintenance Feature

The EMFs screen allows you to view and manage the EMF counters for all modules in all systems.

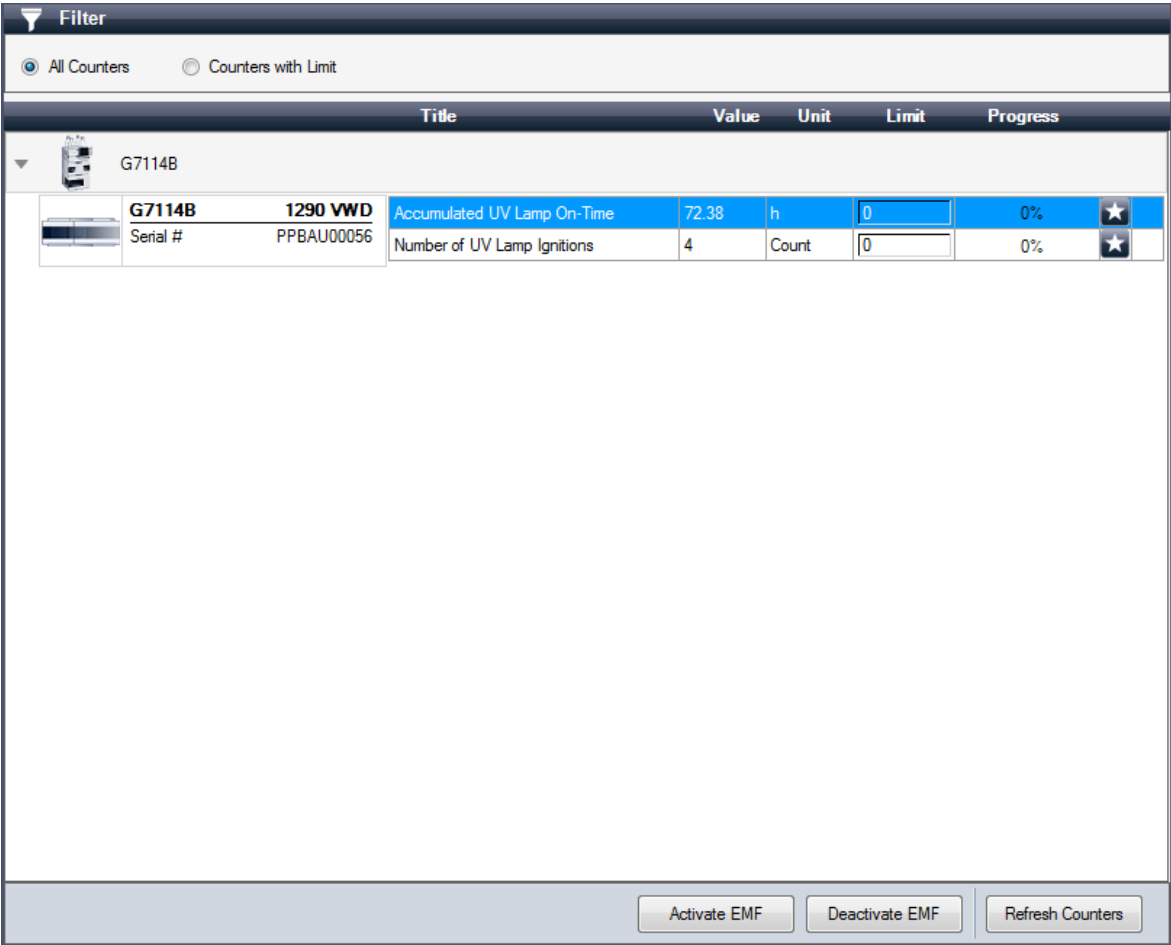


Figure 36: EMFs

This chapter describes the meaning of detector error messages, and provides information on probable causes and suggested actions how to recover from error conditions.

What Are Error Messages 115

General Error Messages 116

- Timeout 116
- Shutdown 116
- Remote Timeout 117
- Lost CAN Partner 118
- Leak 119
- Leak Sensor Open 120
- Leak Sensor Short 121
- Compensation Sensor Open 122
- Compensation Sensor Short 123
- Fan Failed 124
- Open Cover 125
- Cover Violation 126
- ERI Messages 127

Detector Error Messages 129

- UV Lamp: No Current 129
- UV Lamp: No Voltage 129
- Lamp Ignition Failed 130
- No Heater Current 131
- Wavelength Calibration Setting Failed 132
- Wavelength Holmium Check Failed 133
- Grating or Filter Motor Errors 134
- Wavelength Test Failed 135
- Cutoff Filter Doesn't Decrease the Light Intensity at 250 nm 136
- ADC Hardware Error 137

Illegal Temperature Value from Sensor on Main Board 138
Illegal Temperature Value from Sensor at Air Inlet 139
Heater at Fan Assembly Failed 140
Heater Power at Limit 141
Cover Violation 142

What Are Error Messages

Error messages are displayed in the user interface when an electronic, mechanical, or hydraulic (flow path) failure occurs that 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).

If using the InfinityLab Assist, instrument errors will generate a notification. To view the probable causes and recommended actions for this error, click on **Help** button displayed on the notification.

General Error Messages

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

Timeout

Error ID: 62

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.

Shutdown

Error ID: 63

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.	<ul style="list-style-type: none">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.	<ul style="list-style-type: none">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.	<ul style="list-style-type: none">Check external instruments for a shut-down condition.
4	The degasser failed to generate sufficient vacuum for solvent degassing.	<ul style="list-style-type: none">Check the vacuum degasser for an error condition. Refer to the Service Manual for the degasser or the pump that has the degasser built-in.Check the external vacuum degasser module (if installed) for an error condition. Refer to the <i>Service Manual</i> for the degasser or the pump that has the degasser built-in.

Remote Timeout

Error ID: 70

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.	<ul style="list-style-type: none">• Ensure the instrument showing the not-ready condition is installed correctly, and is set up correctly for analysis.
2	Defective remote cable.	<ul style="list-style-type: none">• Exchange the remote cable.
3	Defective components in the instrument showing the not-ready condition.	<ul style="list-style-type: none">• Check the instrument for defects (refer to the instrument’s documentation).

Lost CAN Partner

Error ID: 71

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.

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.	<ul style="list-style-type: none">• Exchange the CAN cable.
3	Defective mainboard in another module.	<ul style="list-style-type: none">• Switch off the system. Restart the system, and determine which module or modules are not recognized by the system.

Leak

Error ID: 64

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 mainboard.

Probable cause		Suggested actions
1	Loose fittings.	• Ensure all fittings are tight.
2	Broken capillary.	• Exchange defective capillaries.
3	Leaking flow cell.	• Exchange flow cell components.

Leak Sensor Open

Error ID: 83

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 on/off switch 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.
4	On/Off switch assembly defective.	• Please contact your Agilent service representative.

Leak Sensor Short

Error ID: 82

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.
3	On/Off switch assembly defective.	• Please contact your Agilent service representative.
4	Cable or contact problem.	• Please contact your Agilent service representative.

Compensation Sensor Open

Error ID: 81

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 on/off switch board and the mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
2	Defective on/off switch assembly.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Compensation Sensor Short

Error ID: 80

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	Defective on/off switch assembly.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
2	Loose connection between the on/off switch board and the mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Fan Failed

Error ID: 68

The cooling fan in the module has failed.

The hall sensor on the fan shaft is used by the mainboard 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.

This limit is given by 2 revolutions/second for longer than 5 seconds.

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 mainboard.	• Please contact your Agilent service representative.

Open Cover

Error ID: 205

The top foam has been removed.

The sensor on the main board detects when the top foam is in place. If the foam is removed, the fan is switched off, and the error message is generated.

Probable cause		Suggested actions
1	The top foam was removed during operation.	• Please contact your Agilent service representative.
2	Foam not activating the sensor.	• Please contact your Agilent service representative.
3	Defective sensor or main board.	• Please contact your Agilent service representative.

Cover Violation

Error ID: 7461

The top foam has been removed.

The sensor on the main board detects when the top foam is in place. If the foam is removed while the lamps are on (or if an attempt is made to switch on for example the lamps with the foam removed), the lamps are switched off, and the error message is generated.

Probable cause		Suggested actions
1	The top foam was removed during operation.	• Please contact your Agilent service representative.
2	Foam not activating the sensor.	• Please contact your Agilent service representative.

ERI Messages

Error ID: 11120, 11121

The ERI (Enhanced Remote Interface) provides two error events related to over current situations on the +5 V and +24 V lines.

Probable cause		Suggested actions
1	The load on the ERI is too high.	<ul style="list-style-type: none">• Reduce the load.

Detector Error Messages

These errors are detector specific.

UV Lamp: No Current

Error ID: 7450

The lamp anode current is missing. 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.	<ul style="list-style-type: none">• Ensure the lamp connector is seated firmly.• Ensure the lamp is connected.
2	Top foam removed while lamp is on.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
3	Defective or non-Agilent lamp.	<ul style="list-style-type: none">• Exchange the lamp.
4	Defective mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
5	Defective power supply.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

UV Lamp: No Voltage

Error ID: 7451

The 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 or non-Agilent lamp.	• Exchange the lamp.
2	Defective power supply.	• Please contact your Agilent service representative.
3	Defective mainboard.	• Please contact your Agilent service representative.

Lamp 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 too hot. Hot gas discharge lamps may not ignite as easily as cold lamps.	<ul style="list-style-type: none">Switch off the lamp and allow it to cool down for at least 15 minutes.
2	Lamp disconnected.	<ul style="list-style-type: none">Ensure the lamp connector is seated firmly.Ensure the lamp is connected.
3	Defective or non-Agilent lamp.	<ul style="list-style-type: none">Exchange the lamp.
4	Defective power supply.	<ul style="list-style-type: none">Please contact your Agilent service representative.
5	Defective mainboard.	<ul style="list-style-type: none">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.	<ul style="list-style-type: none">• Ensure the lamp connector is seated firmly.• Ensure the lamp is connected.
2	Ignition started without the top foam in place.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
3	Fan not running (permitting lamp on).	<ul style="list-style-type: none">• Please contact your Agilent service representative.
4	Defective mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
5	Defective or non-Agilent lamp.	<ul style="list-style-type: none">• Exchange the lamp.
6	Defective power supply.	<ul style="list-style-type: none">• 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.	<ul style="list-style-type: none">• Switch on the lamp.
2	Incorrect flow cell installation.	<ul style="list-style-type: none">• Ensure the flow cell is installed correctly.
3	Flow cell contamination or air bubbles.	<ul style="list-style-type: none">• Clean/replace flow cell windows or remove air bubbles.
4	Intensity too low.	<ul style="list-style-type: none">• Replace lamp.
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.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
7	Defective mainboard.	<ul style="list-style-type: none">• 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 etc.).• 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.

Grating or Filter Motor Errors

Error ID: 7800, 7801, 7802, 7803, 7804, 7805, 7806, 7808, 7809, 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		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.

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.	<ul style="list-style-type: none">Recalibrate the detector.

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.

ADC Hardware Error

Error ID: 7830

A/D-Converter hardware is defective.

Probable cause		Suggested actions
1	A/D-Converter hardware is defective.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

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.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
2	Detector is exposed to illegal ambient conditions.	<ul style="list-style-type: none">• 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">• Please contact your Agilent service representative.
2	Detector is exposed to illegal ambient conditions.	<ul style="list-style-type: none">• Verify that the ambient conditions are within the allowed range.

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.	<ul style="list-style-type: none">• Wait until temperature control equilibrates.

Cover Violation

Error ID: 7461

The top foam has been removed.

The sensor on the main board detects when the top foam is in place. If the foam is removed while the lamps are on (or if an attempt is made to switch on for example the lamps with the foam removed), the lamps are switched off, and the error message is generated.

Probable cause		Suggested actions
1	The top foam was removed during operation.	• Please contact your Agilent service representative.
2	Foam not activating the sensor.	• Please contact your Agilent service representative.

8 Maintenance

This chapter provides general information on maintenance of the module.

Introduction to Maintenance 145

Safety Information Related to Maintenance 146

Overview of Maintenance 148

Cleaning the Module 149

Remove and Install Doors 150

Replace the Deuterium Lamp 153

Replace the Flow Cell / Cuvette Holder 158

Agilent InfinityLab LC Series Variable Wavelength Detectors User Manual 000

Repairing the Flow Cells 162

Using the Cuvette Holder 165

Correcting Leaks 168

Replace Leak Handling System Parts 170

Replace the Module Firmware 171

Introduction to Maintenance

The module is designed for easy maintenance. Maintenance can be done from the front with module in place in the system.

NOTE

There are no serviceable parts inside.
Do not open the module.

Safety Information Related to Maintenance

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.

WARNING**Fire and damage to the module****Wrong fuses**

- Make sure that only fuses with the required rated current and of the specified type (super-fast, fast, time delay etc) are used for replacement.
- The use of repaired fuses and the short-circuiting of fuse-holders must be avoided.

WARNING**Personal injury or damage to the product**

Agilent is not responsible for any damages caused, in whole or in part, by improper use of the products, unauthorized alterations, adjustments or modifications to the products, failure to comply with procedures in Agilent product user guides, or use of the products in violation of applicable laws, rules or regulations.

- Use your Agilent products only in the manner described in the Agilent product user guides.

WARNING**Electrical shock**

Repair work at the module can lead to personal injuries, e.g. shock hazard, when the cover is opened.

- Do not remove the cover of the module.
- Only certified persons are authorized to carry out repairs inside the module.

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

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.

Overview of Maintenance

The following pages describe maintenance (simple repairs) of the detector that can be carried out without opening the main cover.

Table 17: Simple Repairs

Procedures	Typical Frequency	Notes
Deuterium lamp exchange	If noise and/or drift exceeds your application limits or lamp does not ignite.	A wavelength calibration test and an intensity test should be performed after replacement.
Flow cell exchange	If application requires a different flow cell type.	A wavelength calibration test should be performed after replacement.
Cleaning 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.
Leak sensor drying	If leak has occurred.	Check for leaks.
Leak handling system replacement	If broken or corroded.	Check for leaks.

Cleaning the Module

To keep the module case clean, use a soft cloth slightly dampened with water, or a solution of water and mild detergent. Avoid using organic solvents for cleaning purposes. They can cause damage to plastic parts.

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.**

NOTE

A solution of 70 % isopropanol and 30 % water might be used if the surface of the module needs to be disinfected.

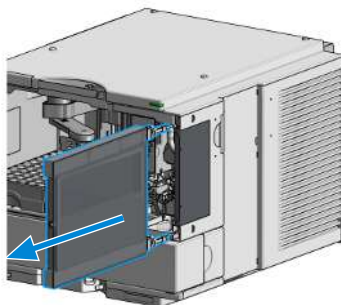
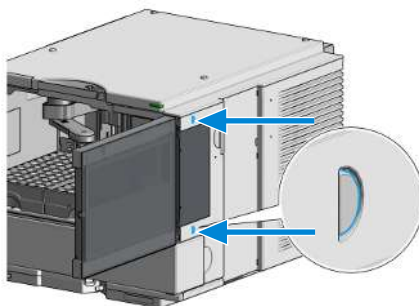
Remove and Install Doors

When	• The instrument doors or the hinges are broken.		
Tools required	Qty.	p/n	Description
	1	 5023-3138	Reversible Screwdriver + Blade 1,0 x 5,5
Parts required (Infinity III)	Qty.	p/n	Description
		 5004-3140	Door Kit Infinity III 140mm
Parts required (Infinity II)	Qty.	p/n	Description
		 5004-0140	Door Kit Infinity II 140mm
Preparations	• Finish any pending acquisition job.		

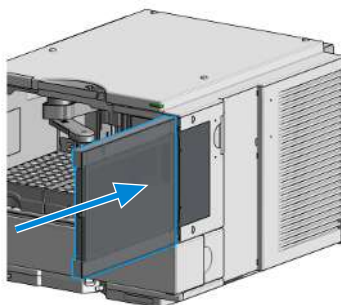
NOTE

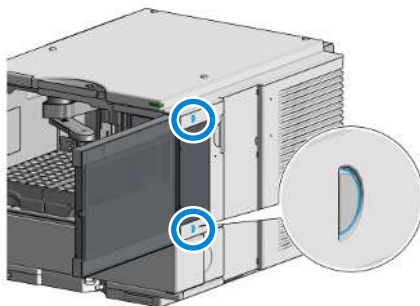
The figures shown in this procedure exemplarily show the Infinity III Vialsampler module. The principle of how to remove and/or install doors works in the same way for all Infinity III modules.

- 1 Press the release buttons and pull the front door out.




- 2 For the Installation of the front door, insert the hinges into their guides and push the door in until the release buttons click into their final position.





Replace the Deuterium Lamp

When	• If noise or drift exceeds application limits or lamp does not ignite.		
Tools required	Qty.	p/n	Description
	1		Screwdriver, Pozidriv #1 PT3
Parts required	Qty.	p/n	Description
	1	 G1314-60101	Deuterium lamp (with RFID tag)
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.

WARNING

Injury by sharp metal edges

- Be careful when touching the RFI sheet metal at the rear of the fan. There are sharp edges.

CAUTION

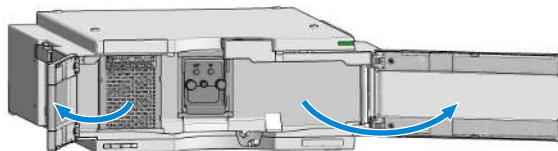
Electronic boards and components are sensitive to electrostatic discharge (ESD).

- To prevent accidental electrostatic discharge when coming into contact with components inside the instrument, touch one of the metal housing panels at the front of the instrument.

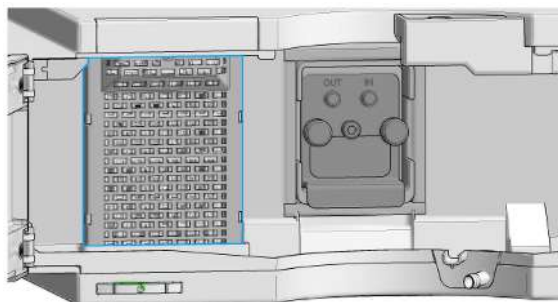
Maintenance

Replace the Deuterium Lamp

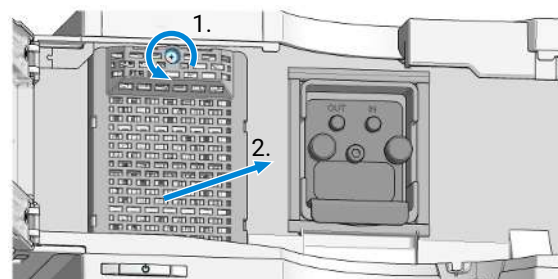
- 1 Open the doors.



- 2 Locate the heater fan cover.

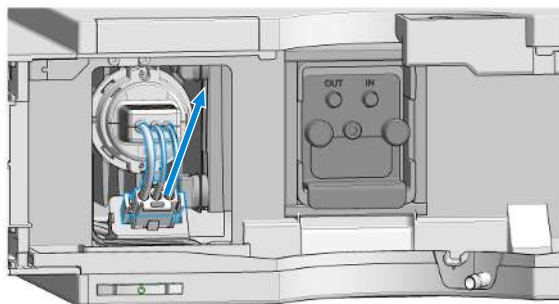


- 3 Unscrew the heater (1.) and remove it (2.).

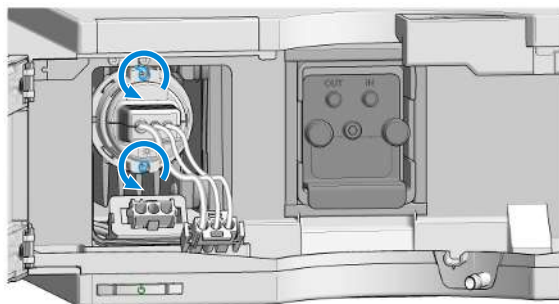


Maintenance**Replace the Deuterium Lamp**

- 4 Unplug the lamp connector.



- 5 Unscrew the two lamp screws (Pozidriv).



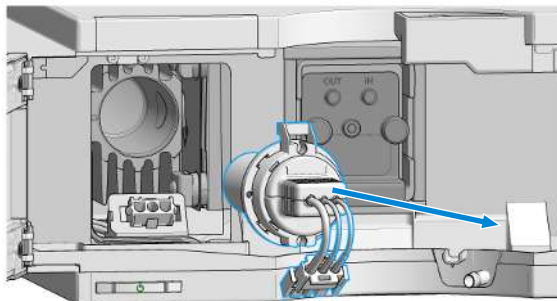
WARNING

Injury by touching hot lamp

If the detector has been in use, the lamp may be hot.

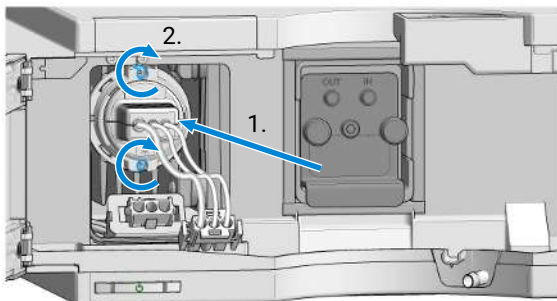
— If so, wait for lamp to cool down.

6 Remove the lamp and place it on a clean place.

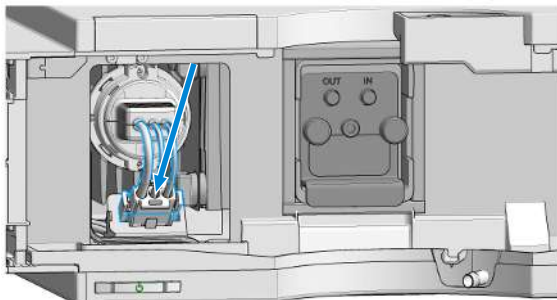
**NOTE**

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.).



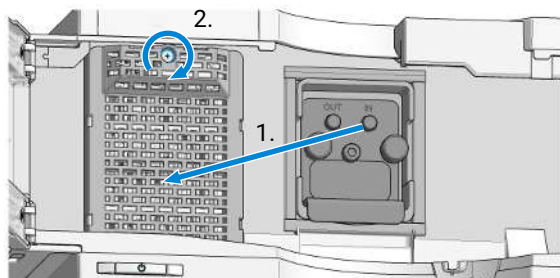
8 Reconnect the connector.



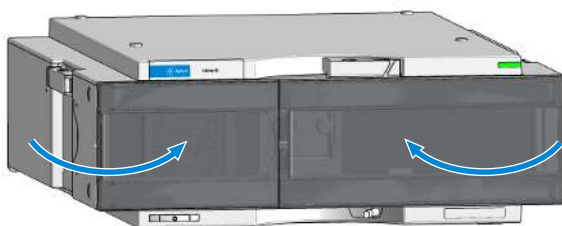
Maintenance

Replace the Deuterium Lamp

- 9 Replace the heater (1.) and fix its screw (2.).



- 10 Close the doors.




- 11 Reset the lamp counter as described in the User Interface documentation (required for non-RFID tag lamps only).
- 12 Turn the lamp ON and give the lamp more than 10 minutes to warm-up.
- 13 Perform a Wavelength Re-calibration after lamp warm-up.

NOTE

After lamp on, the detector requires a warm-up time of 60 min. No measurements should be performed during this time.

Replace the Flow Cell / Cuvette Holder

When • If application needs a different type of flow cell or the flow cell needs repair.

Tools required	Qty.	p/n	Description
	1		Wrench, 1/4 inch, or
	1	 5043-0915	Fitting mounting tool

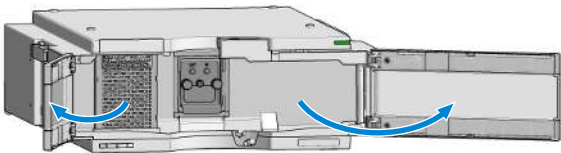
Parts required	Qty.	p/n	Description
			Flow cell

For flow cell details see:

- [Standard Flow Cell 10 mm / 14 µL](#) on page 176
- [Micro Flow Cell 3 mm / 2 µL](#) on page 178
- [Semi-micro Flow Cell \(Parts\)](#) on page 180
- [High Pressure Flow Cell \(Parts\)](#) on page 182
- [Bio Standard Flow Cell](#) on page 183
- [Bio Micro Flow Cell](#) on page 185

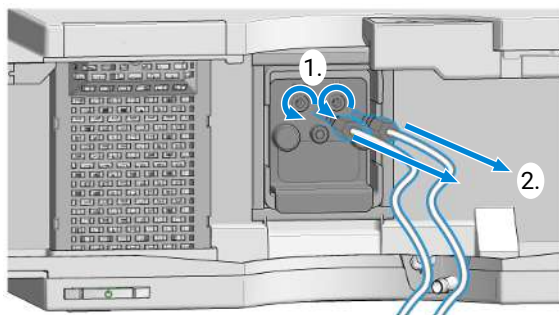
Preparations • Turn the lamp OFF.

- 1 Open the doors.

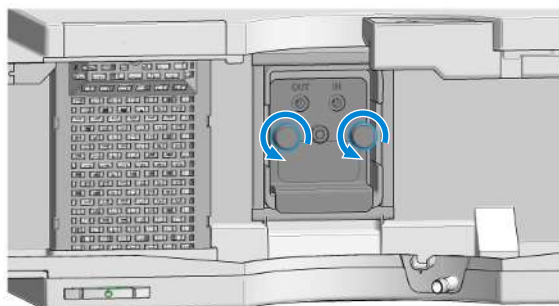


Replace the Flow Cell / Cuvette Holder

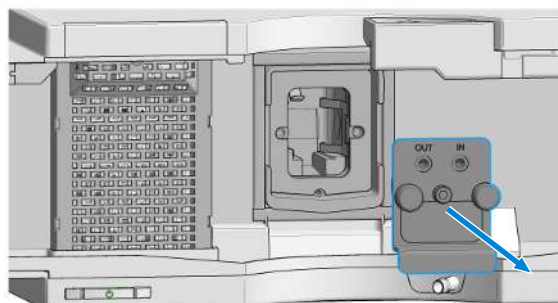
- 2 Disconnect the inlet capillary and outlet tubing from the flow cell.



- 3 Unscrew the two thumb screws.

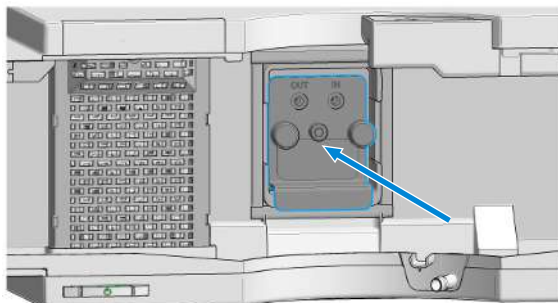


- 4 Pull the flow cell out of its location.

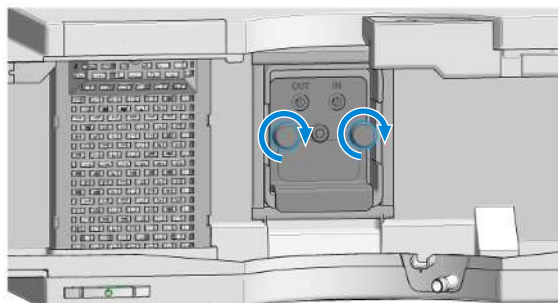


Replace the Flow Cell / Cuvette Holder

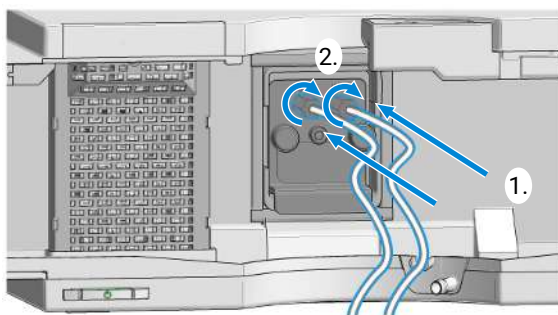
- 5 Insert the flow cell into its location and press it in in the center of the flow cell.



- 6 Fix the two thumb screws parallel and tight.



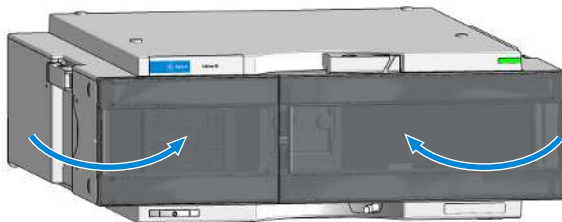
- 7 Reconnect the inlet capillary and the outlet tubing.



Maintenance

Replace the Flow Cell / Cuvette Holder

- 8 Close the doors.



- 9 Configure the flow cell. For further information see Agilent Information Center or the online help of your CDS.
- 10 Perform a **Wavelength Verification-Calibration** to check the correct position of the flow cell/cuvette holder.

Documents about this

- 📄 Agilent InfinityLab LC Series Variable Wavelength Detectors User Manual (<https://www.agilent.com/cs/library/usermanuals/public/G7114-VWD-UseMa-en-SD-29000240.pdf>)

Repairing the Flow Cells

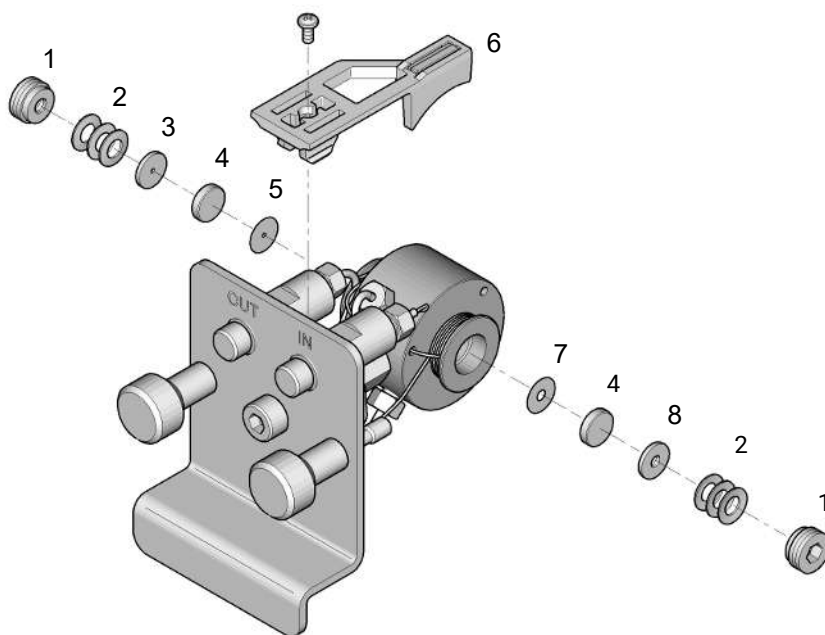
Parts required	Qty.	p/n	Description
	1		Flow cell

For details on flow cells see:

- [Standard Flow Cell 10 mm / 14 µL](#) on page 176
- [Micro Flow Cell 3 mm / 2 µL](#) on page 178
- [Semi-micro Flow Cell \(Parts\)](#) on page 180
- [High Pressure Flow Cell \(Parts\)](#) on page 182
- [Bio Standard Flow Cell](#) on page 183
- [Bio Micro Flow Cell](#) on page 185

NOTE

The shown cell parts will differ depending upon the flow cell type. For detailed parts schematics, refer to above mentioned pages.



- | | | | |
|---|------------------------|---|-----------------|
| 1 | Cell screw | 2 | Conical springs |
| 3 | Ring #1 PEEK | 4 | Window Quartz |
| 5 | Gasket #1 (small hole) | 6 | RFID tag |
| 7 | Gasket #2 (large hole) | 8 | Ring #2 PEEK |

1 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.

CAUTION

Scratched window surfaces by tweezers

Window surfaces can easily be scratched by using tweezers for removing the windows.

— Do not use tweezers to remove windows

- c Use adhesive tape to remove the peek ring, the window and the gasket.
- d Repeat step a through step c for the other window (keep the parts separate - otherwise they could be mixed!).

2 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 it with a piece of lint-free cloth.

NOTE

Always use new gaskets.

3 Reassembling the Flow Cell

- a Hold the flow cell cassette horizontally and place gasket in position. Ensure both cell holes can be seen through the holes of gasket.

NOTE

The semi-micro #1 and #2 gaskets (items 6 and 7, [Semi-micro Flow Cell \(Parts\)](#) on page 180) look very similar. Do not mix them up.

- b Place the window on gasket.
- c Place the peek ring on the window.
- d Insert the conical springs. Make sure the conical springs point towards the window. Otherwise tightening the cell screw might break the window.

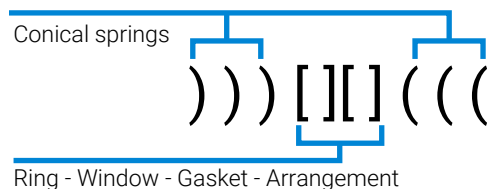


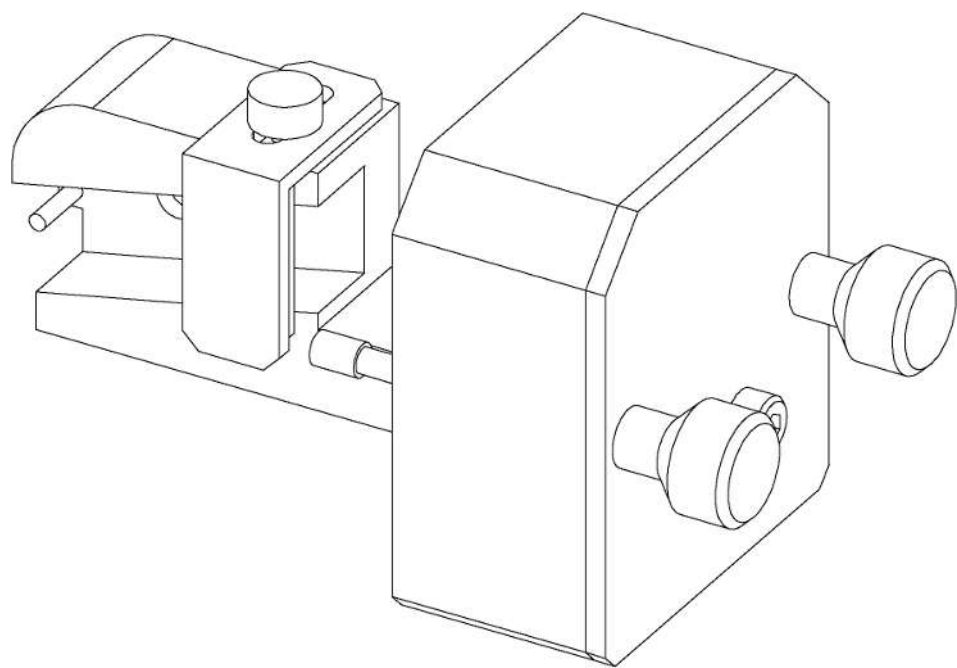
Figure 37: Orientation of conical springs


- e Screw the cell screw into the flow cell and tighten the screw.
- 4 Repeat the procedure for the other cell side.
 - 5 Reconnect the capillaries.
 - 6 Perform a leak test. If OK, insert the flow cell.
 - 7 Perform [Wavelength Calibration](#) on page 95 to check the correct positioning of the flow cell.
 - 8 Replace the front cover.

Using the Cuvette Holder

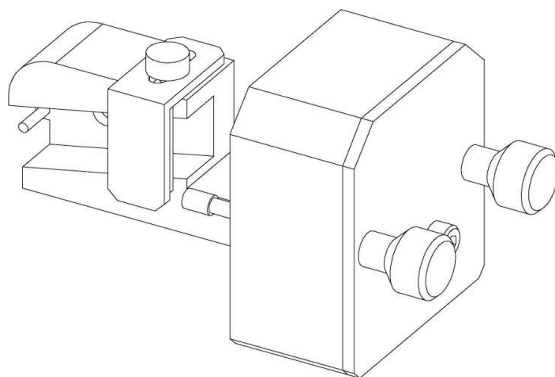
This cuvette holder can be placed instead of a flow cell in the variable wavelength detector. Standard cuvettes with standards in it, for example, National Institute of Standards & Technology (NIST) holmium oxide solution standard, can be fixed in it.

This can be used for wavelength verifications.

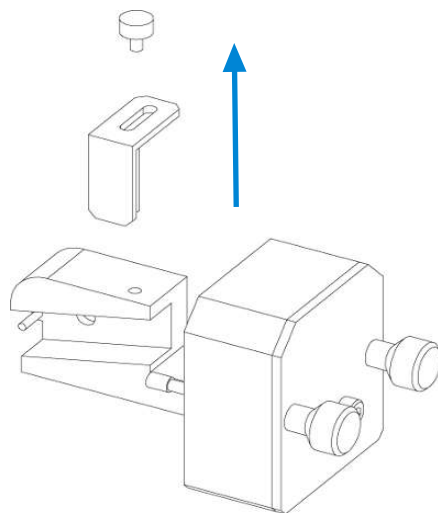


When		<ul style="list-style-type: none">• If your own standard should be used to checkout the instrument.	
Parts required	Qty.	p/n	Description
	1	 G1314-60200	Cuvette Holder
	1		Cuvette with the “standard”, e.g. NIST certified holmium oxide sample
Preparations		<ul style="list-style-type: none">• Remove the normal flow cell.• Have cuvette with standard available.	

- 1 Locate the cuvette holder on the desk.



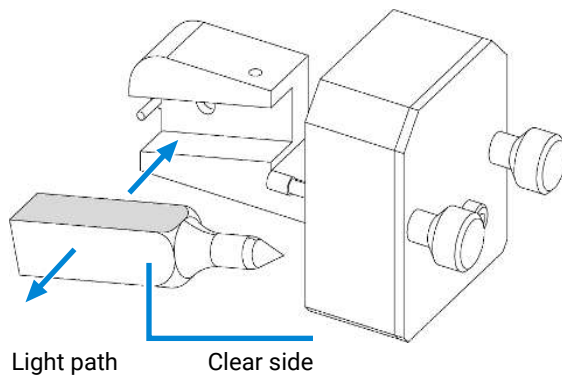
- 2 Unscrew the bracket.



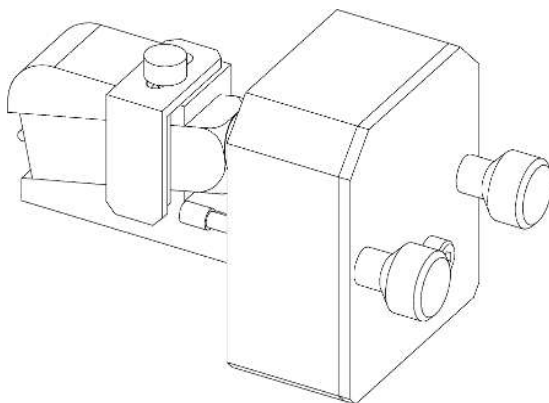
Maintenance

Using the Cuvette Holder

- 3 Insert the cuvette with the sample into the holder. The clear side of the cuvette must be visible.



- 4 Replace the bracket and fix the cuvette.



- 5 Install the cuvette holder in the instrument.
- 6 Perform your Wavelength Verification/Calibration (see [Wavelength Calibration](#) on page 95) to check the correct position of the cuvette holder.

Correcting Leaks

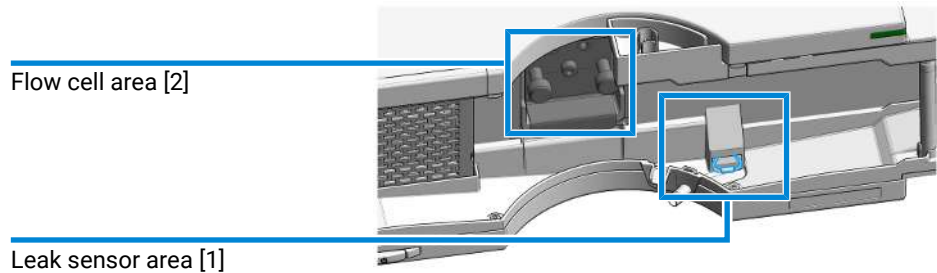
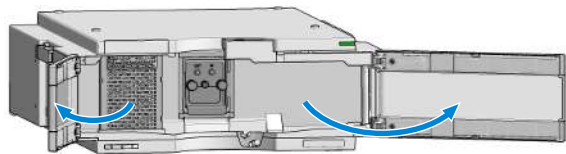


Figure 38: Correcting leaks

When • If a leakage has occurred in the flow cell area or at the capillary connections.

Tools required	Qty.	p/n	Description
	1		Tissue
	1		Wrench, 1/4 inch for capillary connections

- 1 Open the doors.

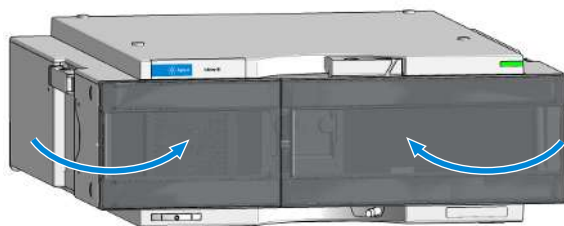


- 2 Use tissue to dry the leak sensor area [1].


Maintenance

Correcting Leaks

- 3 Observe the capillary connections and the flow cell area [2] for leaks and correct, if required.
- 4 Close the doors.



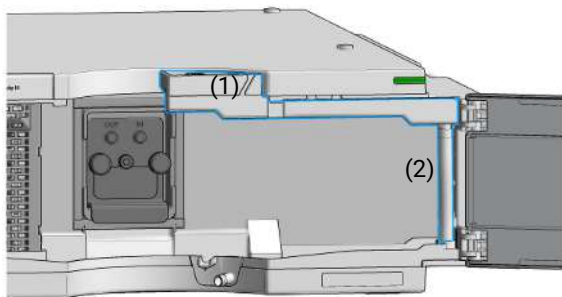
Replace Leak Handling System Parts

Parts required	Qty.	p/n	Description
	1	 5043-0856	Leak Adapter
	1	 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm

Preparations

- Open or remove the doors

- 1 Locate the Leak Adapter [1] and Tubing [2]



- 2 Press the Leak Adapter [1] down and remove it together with the tubing.
- 3 Install the Leak Adapter by pressing it into the Main Cover.
- 4 Insert the Tubing [2] (ca. 85 mm required for replacement) between Leak Adapter outlet and Leak Pan.
- 5 Insert/close the doors.

Replace the Module Firmware

When	Install a newer firmware <ul style="list-style-type: none">• It fixes known problems of older versions, or• It introduces new features, or• It ensures keeping all systems at the same (validated) revision		
When	Install an older firmware <ul style="list-style-type: none">• It ensures keeping all systems at the same (validated) revision, or• It ensures compatibility after adding a new module to the system, or• A third-party control software requires a special version		
Software required	<ul style="list-style-type: none">• Agilent Lab Advisor software		
Tools required	Qty.	p/n	Description
	1		Firmware, tools and documentation from Agilent web site
Preparations	<ul style="list-style-type: none">• Read update documentation provided with the Firmware Update Tool. <p>To upgrade/downgrade the module’s firmware carry out the following steps:</p> <ol style="list-style-type: none">1 Download the required module firmware, the latest FW Update Tool and the documentation from the Agilent web. https://www.agilent.com/en-us/firmwareDownload?whid=697612 For loading the firmware into the module follow the instructions in the documentation.		

Module Specific Information

Table 18: Module Specific Information (G7114A/B)

	G7114B	G7114A
Initial firmware	D.06.70	D.07.01
Compatibility with 1100 / 1200 series modules	When using the G7114A/B in a system, all other modules must have firmware from set 6.50 or above (main and resident). Otherwise the communication will not work.	
Conversion / emulation	Possible via Lab Advisor software: <ul style="list-style-type: none">• G7114B: G1314E/F• G7114A: G1314F	

9

Parts and Materials for Maintenance

This chapter provides information on parts for maintenance.

Overview of Maintenance Parts 174

Standard Flow Cell 10 mm / 14 μ L 176

Micro Flow Cell 3 mm / 2 μ L 178

Semi-micro Flow Cell (Parts) 180

High Pressure Flow Cell (Parts) 182

Bio Standard Flow Cell 183

Bio Micro Flow Cell 185

Cuvette Holder (Parts) 187

Accessory Kit 188

Overview of Maintenance Parts

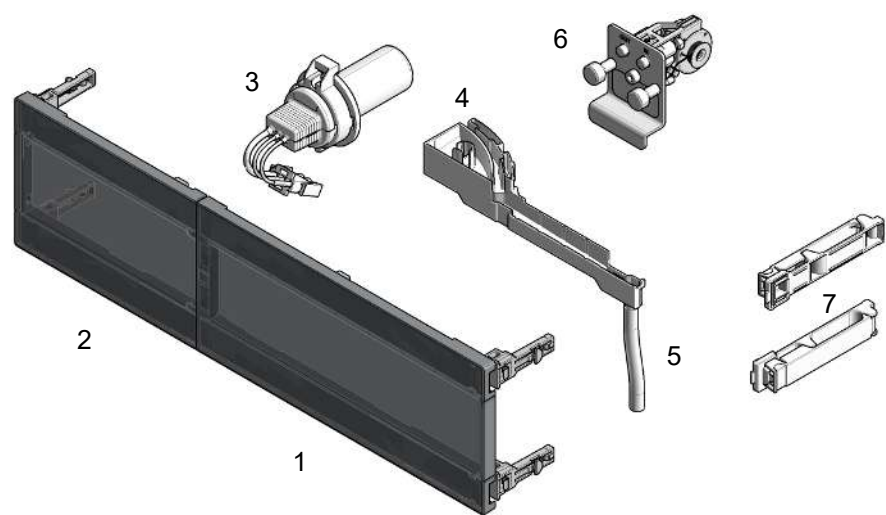




















Figure 39: Maintenance Parts

#	Qty.	p/n	Description
1	1	 5360-0015	Door 140mm right Infinity III (only orderable as part of 5004-3140 Door Kit Infinity III 140mm)
2	1	 5360-0016	Door 140mm left Infinity III (only orderable as part of 5004-3140 Door Kit Infinity III 140mm)
1	1	 5360-0003	Door 140mm right Infinity II (only orderable as part of 5004-0140 Door Kit Infinity II 140mm)
2	1	 5360-0002	Door 140mm left Infinity II (only orderable as part of 5004-0140 Door Kit Infinity II 140mm)
3	1	 G1314-60101	Deuterium lamp (with RFID tag)
4	1	 5043-0856	Leak Adapter
5	1	 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm

#	Qty.	p/n	Description
6	1	 G1314-60186	Standard flow cell 10 mm, 14 µL, 40 bar OR
	1	 G1314-60187	Micro flow cell 3 mm, 2 µL, 120 bar OR
	1	 G1314-60183	Semi-micro flow cell 6 mm, 5 µL OR
	1	 G1314-60182	High pressure flow cell 10 mm, 14 µL, 400 bar OR
	1	 G1314-60023	Prep flow cell 0.06 mm OR
	1	 G1314-60025	Prep flow cell 0.3 mm OR
	1	 G1314-60024	Prep flow cell 3 mm OR
	1	 G1314-60188	Bio standard flow cell VWD, 10 mm, Cell Vol. 14 µl, Sapphire, MP35N OR
	1	 G1314-60189	Bio micro flow cell VWD, 3 mm, Cell Vol. 2 µl, Sapphire, MP35N
	1	 5062-8535	Waste accessory kit (Flow Cell to waste) (Flow Cell to Waste)
7	1	 5043-1013	Tubing Clip

Standard Flow Cell 10 mm / 14 μ L

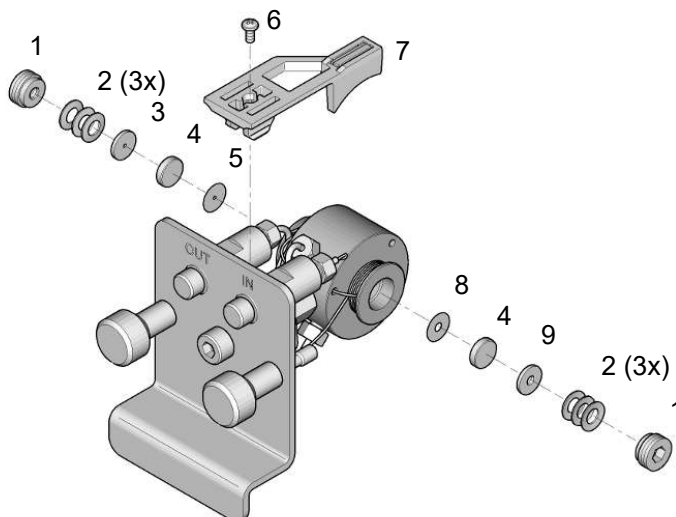


Figure 40: Standard Flow Cell

#	Qty.	p/n	Description
	1	G1314-60186	Standard flow cell 10 mm, 14 μ L, 40 bar
	1	5062-8522	Capillary column - detector PEEK 600 mm lg, 0.17 mm i.d., 1/16 inch o.d.
	1	G1314-65061	Cell Repair Kit, includes 2x Gasket #1, 2x Gasket #2, 2x Window Quartz
1	1	G1314-65062	Cell screw kit
2	1	79853-29100	Conical spring kit, 10/pk
3	1	G1314-65065	Ring #1 kit (OUT large hole, i.d. 2.4 mm) PEEK, 2/pk
4	1	79853-68742	Window quartz kit, 2/pk
5	1	G1314-65063	Gasket #1 kit (OUT large hole, i.d. 2.4 mm) KAPTON, 2/pk
6	1	0515-4780	Screw for Clip, M2.2, 4.5 mm long
7	1	G1314-44010	Clip for RFI ID tag

Parts and Materials for Maintenance

Standard Flow Cell 10 mm / 14 µL

#	Qty.	p/n	Description
8	1	 G1314-65064	Gaskets #2 IN (small hole i.d. 1 mm), KAPTON 10/pk
9	1	 G1314-65066	Ring #2 kit (IN small hole, i.d. 1 mm) PEEK, 2/pk

Micro Flow Cell 3 mm / 2 μ L

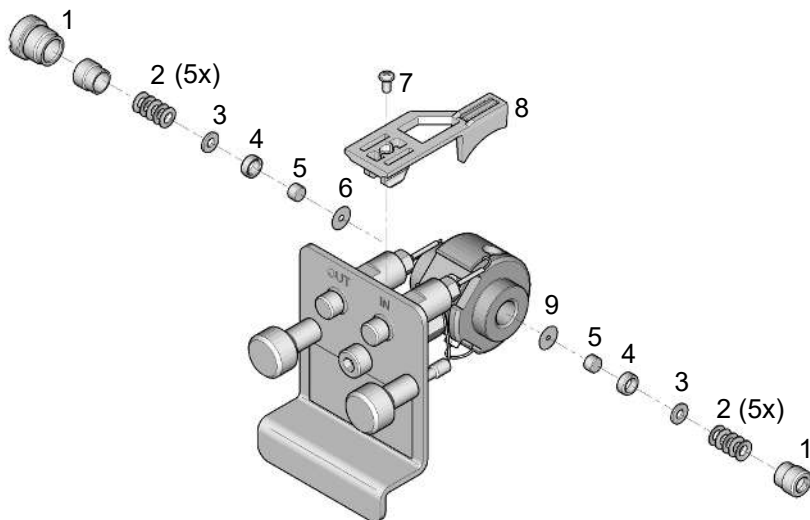


Figure 41: Micro Flow Cell

#	Qty.	p/n	Description
1		G1314-60187	Micro flow cell 3 mm, 2 μ L, 120 bar
1		5021-1823	Capillary column – detector SST 400 mm lg, 0.12 mm i.d.
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		79883-68702	Gasket BACK (PTFE), 1.8 mm hole, outlet side (12/ pk)
7		0515-4780	Screw for Clip, M2.2, 4.5 mm long
8		G1314-44010	Clip for RFI ID tag

Parts and Materials for Maintenance
Micro Flow Cell 3 mm / 2 µL

#	Qty.	p/n	Description
9	1	 G1315-68710	Gasket FRONT (PTFE), 1.3 mm hole, inlet side (12/ pk)
	1	 G1314-87301	Capillary IN (0.12 mm, 310 mm lg)
	1	 G1314-87302	Capillary OUT (0.17 mm, 120 mm lg)
	1	 G1315-68713	Cell repair kit semi-micro, includes window screw kit, Gasket Kit BACK, Gasket Kit FRONT and 4 mm hexagonal wrench
	1	 79883-68703	Window screw kit, includes 2 quartz windows, 2 compression washers, 2 window holders, 2 window screws and 10 washers

Semi-micro Flow Cell (Parts)

NOTE

The semi-micro #1 and #2 gaskets (items 6 and 7) look very similar. Do not mix them up.

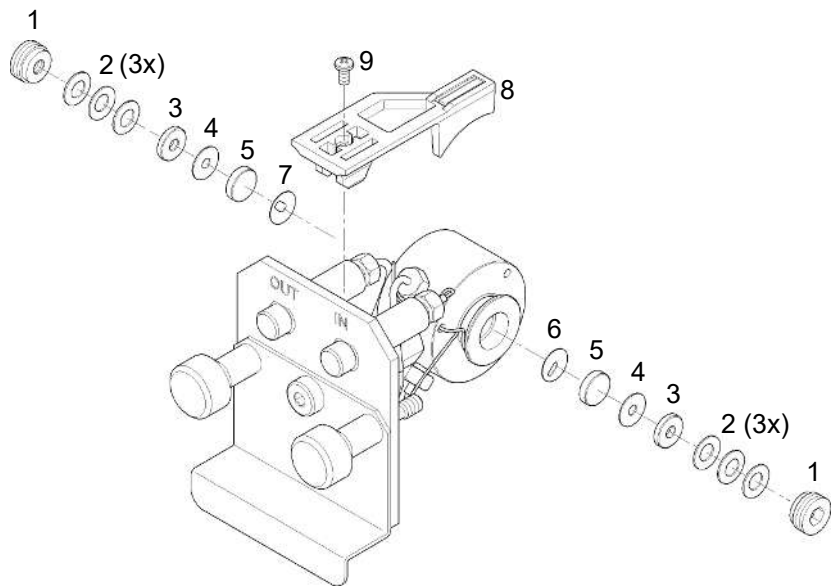





Figure 42: Semi-micro Flow Cell

#	Qty.	p/n	Description
1	1	G1314-60183	Semi-micro flow cell 6 mm, 5 μ L
1	1	5021-1823	Capillary column – detector SST 400 mm lg, 0.12 mm i.d.
1	1	G1314-20047	Cell screw
	1	G1314-65056	Semi-micro cell kit, includes two quartz windows, one gasket #1, one #2 and two PTFE gaskets.
2	1	79853-29100	Conical spring kit, 10/pk
3	1	79853-22500	Ring SST, 2/pk
4	1	79853-68743	PTFE gasket (round hole i.d. 2.5 mm, o.d. 8 mm), (10/pk)

Parts and Materials for Maintenance**Semi-micro Flow Cell (Parts)**

#	Qty.	p/n	Description
5	1	 79853-68742	Window quartz kit, 2/pk
6	1		Semi-micro #1 gasket (long hole 1.5 x 3.5), PTFE
7	1		Semi-micro #2 gasket (long hole 2 x 4), PTFE
8	1	 G1314-44010	Clip for RFI ID tag
9	1	 0515-4780	Screw for Clip, M2.2, 4.5 mm long

High Pressure Flow Cell (Parts)

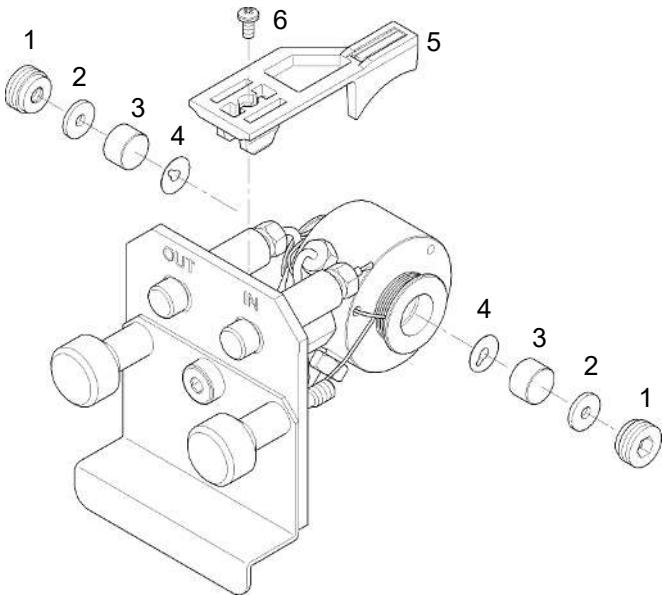


Figure 43: High Pressure Flow Cell

#	Qty.	p/n	Description
	1	G1314-60182	High pressure flow cell 10 mm, 14 μ L, 400 bar
	1	G1315-87311	Capillary ST 0.17 mm x 380 mm S/S
1	1	G1314-20047	Cell screw
	1	G1314-65054	Cell kit Agilent, comprises: two windows, two KAPTON gaskets and two PEEK rings
2	1		Ring PEEK kit
3	1		Window quartz kit
4	1		Gasket kit, KAPTON
5	1	G1314-44010	Clip for RFI ID tag
6	1	0515-4780	Screw for Clip, M2.2, 4.5 mm long

Bio Standard Flow Cell

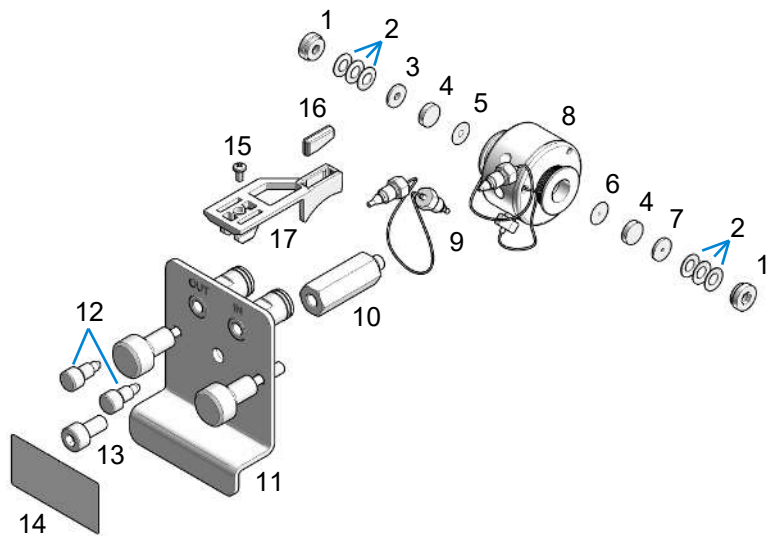









Figure 44: Bio Standard Flow Cell

#	Qty.	p/n	Description
1	2	G1314-20047	Cell screw
2	6	5022-3942	Disc spring
3	1	G1314-20046	Ring (Out)
4	2	5190-0939	Sapphire window VWD
5	1	G1314-40044	Gasket Out Bio
6	1	G1314-40043	Gasket In Bio
7	1	G1314-20045	Ring (In)
8	1	G1314-60054	Cell Block Assembly VWD MP35N Fitting
9	1	5500-1297	Capillary MP35N 0.3 x 120 mm
10	1	G1314-20011	Cell shaft
11	1	G1314-60058	Cell-Plate Assembly Bio

Parts and Materials for Maintenance
Bio Standard Flow Cell

#	Qty.	p/n		Description
12	2		0100-1259	Plug-Screw 1032- Fitting
13	1		0515-1096	SCR-SKT-HD HEX M5X0.8 10 SST PSVT
14	1		G1314-87012	Label Standard RFID Bio
15	1		0515-4780	Screw for Clip, M2.2, 4.5 mm long
16	1		0960-2971	RF Transponder
17	1		G1314-44010	Clip for RFI ID tag
	1		0470-0960	Sealant Methacrylate Ester Liquid

Bio Micro Flow Cell

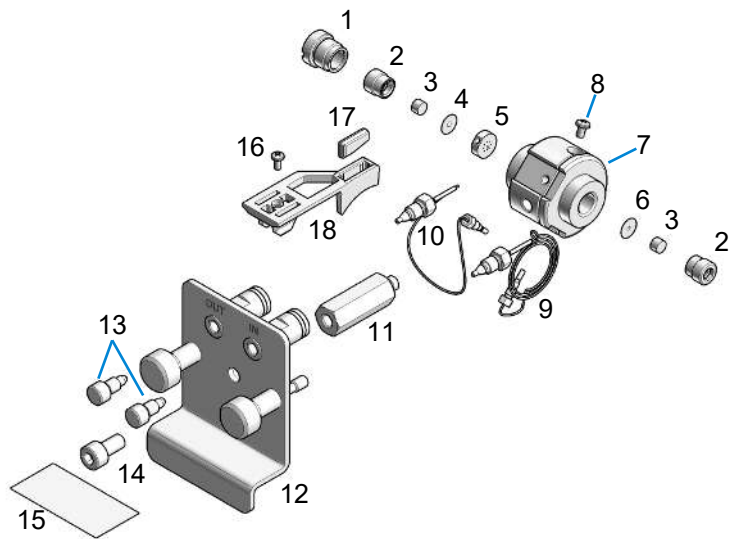











Figure 45: Bio Micro Flow Cell

#	Qty.	p/n	Description
1	1	79883-21701	Screw-Bushing 6
2	2	G1315-60021	Cell screw assembly (comprises window screw, spring washers, compression washer, window holder and quartz window)
3	2	5190-0921	Sapphire window
4	1	79883-07101	Gasket
5	1	G1314-27201	VWD Cell Insert 3 mm MP35N
6	1	G1315-07101	Gasket Front 1.3
7	1	G1314-27700	VWD Cell Body 3 mm
8	1	0515-1056	Screw M 2.5, 4 mm lg
9	1	G1314-67003	Cap In VWD 3 mm MP35N
10	1	G1314-67004	Cap Out VWD 3 mm MP35N

Parts and Materials for Maintenance

Bio Micro Flow Cell

#	Qty.	p/n	Description
11	1	 G1314-20011	Cell shaft
12	1	 G1314-60058	Cell-Plate Assembly Bio
13	2	 0100-1259	Plug-Screw 1032- Fitting
14	1	 0515-1096	SCR-SKT-HD HEX M5X0.8 10 SST PSVT
15	1	 G1314-87013	Label micro RFID Bio
16	1	 0515-4780	Screw for Clip, M2.2, 4.5 mm long
17	1	 0960-2971	RF Transponder
18	1	 G1314-44010	Clip for RFI ID tag
	1	 0470-0960	Sealant Methacrylate Ester Liquid

Cuvette Holder (Parts)

For information the use of the cuvette holder, refer to [Using the Cuvette Holder](#) on page 165.

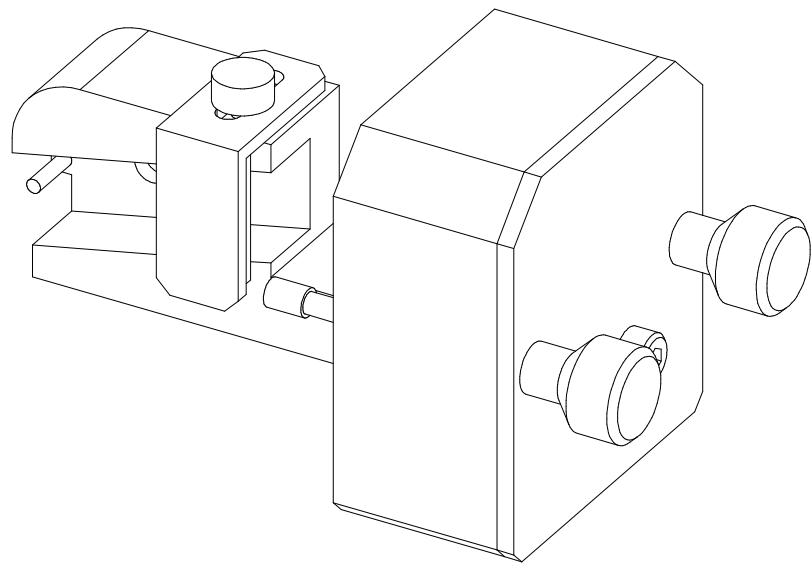









Figure 46: Cuvette Holder

p/n	Description
 G1314-60200	Cuvette Holder

Accessory Kit

Accessory kit (G7114-68755) contains some accessories and tools needed for installation and repair of the module.

p/n	Description
 5062-8535	Waste accessory kit (Flow Cell to waste)
 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm (see Item 4 in Figure 39 on page 174)
 5181-1516	CAN cable, Agilent module to module, 0.5 m
 5500-1155	Tube Connector, 90 degree, ID 6.4
 5043-1013	Tubing Clip (see item 7 in Figure 39 on page 174)
 0100-1516	Finger-tight fitting PEEK, 2/pk

10

Identifying Cables

This chapter provides information on cables used with the modules.

Cable Overview 190

Analog Cables 192

Remote Cables 194

BCD Cables 198

CAN/LAN Cables 200

RS-232 Cables 201

USB 202

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 35900A A/D converter
01046-60105	Analog cable (BNC to general purpose, spade lugs)

Remote cables

p/n	Description
5188-8029	ERI to general purpose
5188-8044	Remote Cable ERI – ERI
5188-8045	Remote Cable APG – ERI
5188-8059	ERI-Extension-Cable 1.2 m
5061-3378	Remote Cable to 35900 A/D converter
01046-60201	Agilent module to general purpose
5188-8057	Fraction Collection ERI remote Y-cable

CAN cables

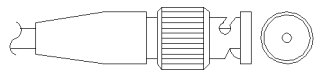
p/n	Description
5181-1516	CAN cable, Agilent module to module, 0.5 m
5181-1519	CAN cable, Agilent module to module, 1 m

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)

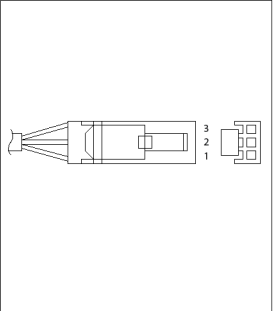
RS-232 cables	p/n	Description
	RS232-61601	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It is also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3, 3-2, 4-6, 5-5, 6-4, 7-8, 8-7, 9-9.
	5181-1561	RS-232 cable, 8 m
USB cables	p/n	Description
	5188-8050	USB A M-USB Mini B 3 m (PC-Module)
	5188-8049	USB A F-USB Mini B M OTG (Module to Flash Drive)

Analog Cables

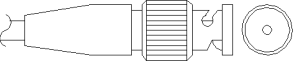


One end of these cables provides a BNC connector to be connected to Agilent modules. The other end depends on the instrument to which connection is being made.

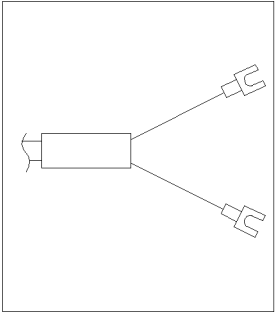
Agilent Module to 35900 A/D converters

p/n 35900-60750	35900	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 +

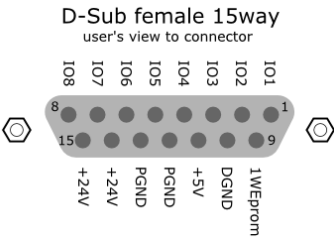
Agilent Module to General Purpose

p/n 01046-60105	Pin	Pin Agilent module	Signal Name
	1		Not connected
	2	Black	Analog -
	3	Red	Analog +

Remote Cables

ERI (Enhanced Remote Interface)

- 5188-8029 ERI to general purpose (D-Sub 15 pin male - open end)
- 5188-8044 ERI to ERI (D_Sub 15 pin male - male)
- 5188-8059 ERI-Extension-Cable 1.2 m (D-Sub15 pin male / female)

p/n 5188-8029	pin	Color code	Enhanced Remote	Classic Remote	Active (TTL)
	1	white	IO1	START REQUEST	Low
	2	brown	IO2	STOP	Low
	3	green	IO3	READY	High
	4	yellow	IO4	PEAK DETECT	Low
	5	grey	IO5	POWER ON	High
	6	pink	IO6	SHUT DOWN	Low
	7	blue	IO7	START	Low
	8	red	IO8	PREPARE	Low
	9	black	1wire DATA		
	10	violet	DGND		
	11	grey-pink	+5V ERI out		
	12	red-blue	PGND		
	13	white-green	PGND		
	14	brown-green	+24V ERI out		
	15	white-yellow	+24V ERI out		
	NC	yellow-brown			


NOTE

Configuration is different with old firmware revisions.
The configuration for IO4 and IO5 is swapped for modules with firmware lower than D.07.10.

NOTE


Peak Detection is used for LCMS systems connected with the Fraction Collection Remote Y-Cable (5188-8057).

- 5188-8045 ERI to APG (Connector D_Subminiature 15 pin (ERI), Connector D_Subminiature 9 pin (APG))

p/n 5188-8045		Pin (ERI)	Signal	Pin (APG)	Active (TTL)
		10	GND	1	
		1	Start Request	9	Low
		2	Stop	8	Low
		3	Ready	7	High
		5	Power on	6	High
		4	Future	5	
		6	Shut Down	4	Low
		7	Start	3	Low
		8	Prepare	2	Low
		Ground	Cable Shielding	NC	

- 5188-8057 ERI to APG and RJ45 (Connector D_Subminiature 15 pin (ERI), Connector D_Subminiature 9 pin (APG), Connector plug Cat5e (RJ45))

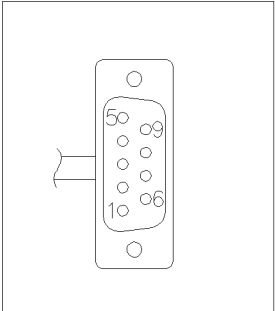
Table 19: 5188-8057 ERI to APG and RJ45

p/n 5188-8057	Pin (ERI)	Signal	Pin (APG)	Active (TTL)	Pin (RJ45)
	10	GND	1		5
	1	Start Request	9	High	
	2	Stop	8	High	
	3	Ready	7	High	
	4	Fraction Trigger	5	High	4
	5	Power on	6	High	
	6	Shut Down	4	High	
	7	Start	3	High	
	8	Prepare	2	High	
	Ground	Cable Shielding	NC		

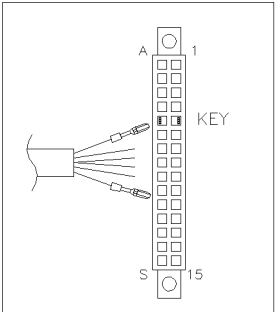


One end of these cables provides an Agilent Technologies APG (Analytical Products Group) remote connector to be connected to Agilent modules. The other end depends on the instrument to be connected to.

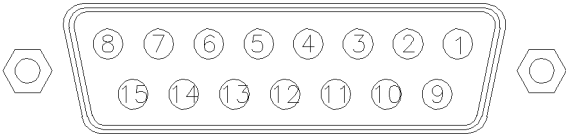
Agilent Module to Agilent 35900 A/D Converters

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

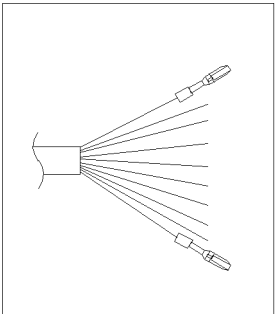
p/n 01046-60201	Wire Color	Pin Agilent module	Signal Name	Active (TTL)
	White	1	Digital ground	
	Brown	2	Prepare run	Low
	Gray	3	Start	Low
	Blue	4	Shut down	Low
	Pink	5	Not connected	
	Yellow	6	Power on	High
	Red	7	Ready	High
	Green	8	Stop	Low
	Black	9	Start request	Low

BCD Cables

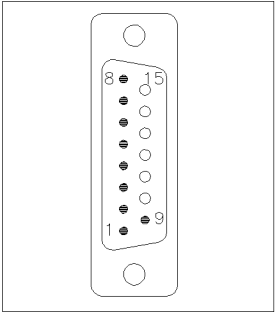


One end of these cables provides a 15-pin BCD connector to be connected to the Agilent modules. The other end depends on the instrument to be connected to

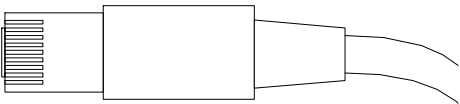
Agilent Module to General Purpose

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

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)

RS-232 Cables

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 is 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

USB

To connect a USB Flash Drive use a USB OTG cable with Mini-B plug and A socket.

p/n	Description
5188-8050	USB A M-USB Mini B 3 m (PC-Module)
5188-8049	USB A F-USB Mini B M OTG (Module to Flash Drive)

This chapter describes the module in more detail on hardware and electronics.

General Hardware Information 204

Firmware Description 204

Electrical Connections 206

Interfaces 208

Instrument Layout 215

Early Maintenance Feedback (EMF) 216

Module-Specific Hardware Information 218

Setting the 6-bit Configuration Switch 218

Early Maintenance Feedback (EMF) 220

General Hardware Information

This section provides detailed hardware information on firmware that is valid for this module.

Firmware Description

The firmware of the instrument consists of two independent sections:

- a non-instrument specific section, called *resident system*
- an instrument specific section, called *main system*

Resident System

This resident section of the firmware is identical for all Agilent 1100/1200/1220/1260/1290 series modules. Its properties are:

- the complete communication capabilities (CAN, LAN, USB and RS- 232)
- memory management
- ability to update the firmware of the 'main system'

Main System

Its properties are:

- the complete communication capabilities (CAN, LAN, USB and RS- 232)
- memory management
- ability to update the firmware of the 'resident system'

In addition the main system comprises the instrument functions that are divided into common functions like

- run synchronization through APG/ERI remote,
- error handling,
- diagnostic functions,

- or module specific functions like
 - internal events such as lamp control, filter movements,
 - raw data collection and conversion to absorbance.

Firmware Updates

Firmware updates can be done with the Agilent Lab Advisor software with files on the hard disk (latest version should be used).

Required tools, firmware and documentation are available from the Agilent web: <https://www.agilent.com/en-us/firmwareDownload?whid=69761>

The file naming conventions are:

PPPP_RVVV_XXX.dlb, where

- PPPP is the product number, for example, 1315B for the G1315B DAD,
- R the firmware revision, for example, A for G1315B or B for the G1315C DAD,
- VVV is the revision number, for example 650 is revision 6.50,
- XXX is the build number of the firmware.

For instructions on firmware updates refer to section *Replacing Firmware* in chapter *Maintenance* or use the documentation provided with the *Firmware Update Tools*.

NOTE

Update of main system can be done in the resident system only. Update of the resident system can be done in the main system only.
Main and resident firmware must be from the same set.

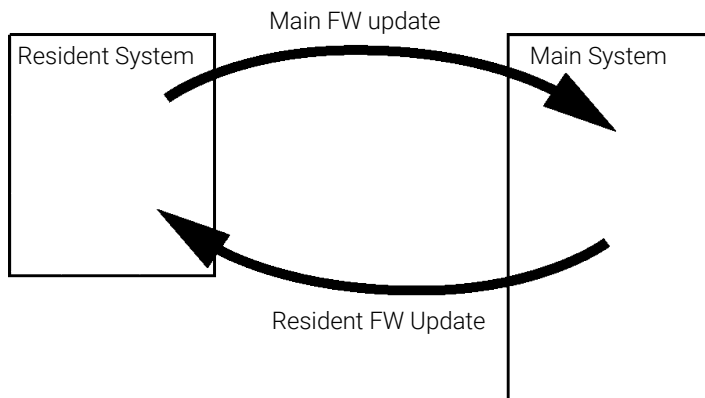


Figure 47: Firmware update mechanism

NOTE

Some modules are limited in downgrading due to their mainboard version or their initial firmware revision. For example, a G1315C DAD SL cannot be downgraded below firmware revision B.01.02 or to a A.xx.xx.

Some modules can be re-branded (e.g. G1314C to G1314B) to allow operation in specific control software environments. In this case, the feature set of the target type is used and the feature set of the original one is lost. After re-branding (e.g. from G1314B to G1314C), the original feature set is available again.

All this specific information is described in the documentation provided with the firmware update tools.

The firmware update tools, firmware and documentation are available from the Agilent web.

- <https://www.agilent.com/en-us/firmwareDownload?whid=69761>

Electrical Connections

- The CAN bus is a serial bus with high-speed data transfer. The two connectors for the CAN bus are used for internal module data transfer and synchronization.
- One analog output provides signals for integrators or data handling systems.
- The ERI connector may be used in combination with other analytical instruments from Agilent Technologies if you want to use features such as start, stop, common shut down, prepare, and so on.

- With the appropriate software, the LAN connector may be used to control the module from a computer through a LAN connection. This connector is activated and can be configured with the configuration switch.
- With the appropriate software, the USB connector may be used to control the module from a computer through a USB connection.
- The power input socket accepts a line voltage of 100 – 240 VAC \pm 10 % with a line frequency of 50 or 60 Hz. Maximum power consumption varies by module. There is no voltage selector on your module because the power supply has wide-ranging capability. There are no externally accessible fuses because automatic electronic fuses are implemented in the power supply.

WARNING

Electric shock due to insufficient insulation of connected instruments

Personal injury or damage to the instrument

- **Any other instruments connected to this instrument shall be approved to a suitable safety standard and must include reinforced insulation from the mains.**

NOTE

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

Rear View of the Module

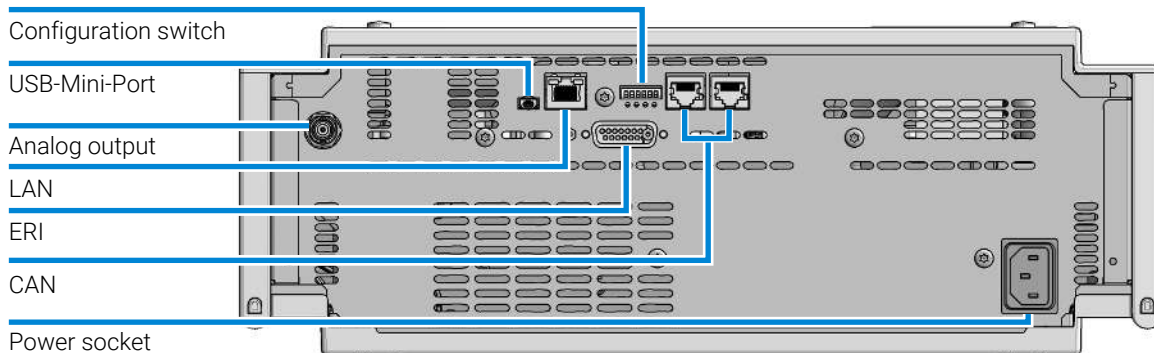


Figure 48: Rear view of detector (example shows a G7114A/B VWD) – electrical connections and label

Serial Number Information

The serial number information on the instrument labels provide the following information:

CCXZZ00000	Format
CC	Country of manufacturing <ul style="list-style-type: none">• DE = Germany• JP = Japan• CN = China
X	Alphabetic character A-Z (used by manufacturing)
ZZ	Alpha-numeric code 0-9, A-Z, where each combination unambiguously denotes a module (there can be more than one code for the same module)
00000	Serial number

Interfaces

The Agilent InfinityLab LC Series modules provide the following interfaces:

Table 20: Agilent InfinityLab LC Series interfaces

Module	CAN	USB	LAN (on-board)	RS-232	Analog	APG (A) / ERI (E)	Special
Pumps							
G7104A/C	2	No	Yes	Yes	1	A	
G7110B	2	Yes	Yes	No	No	E	
G7111A/B, G5654A	2	Yes	Yes	No	No	E	
G7112B	2	Yes	Yes	No	No	E	
G7120A, G7132A	2	No	Yes	Yes	1	A	
G7161A/B	2	Yes	Yes	No	No	E	
Samplers							
G7129A/B/C	2	Yes	Yes	No	No	E	
G7167A/B/C, G7137A, G5668A, G3167A	2	Yes	Yes	No	No	E	

Module	CAN	USB	LAN (on-board)	RS-232	Analog	APG (A) / ERI (E)	Special
G7157A	2	Yes	Yes	No	No	E	
Detectors							
G7114A/B	2	Yes	Yes	No	1	E	
G7115A	2	Yes	Yes	No	1	E	
G7117A/B/C	2	Yes	Yes	No	1	E	
G7121A/B	2	Yes	Yes	No	1	E	
G7162A/B	2	Yes	Yes	No	1	E	
G7165A	2	Yes	Yes	No	1	E	
Fraction Collectors							
G7158B	2	Yes	Yes	No	No	E	
G7159B	2	Yes	Yes	No	No	E	
G7166A	2	No	No	No	No	No	Requires a host module with on-board LAN with minimum FW B.06.40 or C.06.40, or with additional G1369C LAN Card
G1364E/F, G5664B	2	Yes	Yes	No	No	E	THERMOSTAT for G1330B
Others							
G1170A	2	No	No	No	No	No	Requires a host module with on-board LAN or with additional G1369C LAN Card.
G7116A/B	2	No	No	No	No	No	Requires a host module with on-board LAN or with additional G1369C LAN Card.
G7122A	No	No	No	Yes	No	A	
G7170B	2	No	No	No	No	No	Requires a host module with on-board LAN with minimum FW B.06.40 or C.06.40, or with additional G1369C LAN Card

NOTE

LAN connection is made between at least one of the Agilent modules and the Control PC.

- If an Assist Hub is installed, connect the LAN to the Lab LAN port of this module.
 - If an Assist Hub is NOT installed and a detector (DAD/MWD/FLD/VWD/RID) is installed, connect the LAN to this module.
 - If an Assist Hub is NOT installed and there are multiple detectors with spectral capabilities, consider using additional LAN connections for each detector.
 - If an Assist Hub is installed, connect additional LAN connections from the detectors and pumps to the Assist Hub.
-
- CAN connectors as interface to other modules
 - LAN connector as interface to the control software
 - RS-232C as interface to a computer
 - USB (Universal Series Bus) as interface to a computer
 - REMOTE connector as interface to other Agilent products
 - Analog output connector for signal output

Overview Interfaces

CAN

The CAN is inter-module communication interface. It is a 2-wire serial bus system supporting high speed data communication and real-time requirement.

LAN

The modules have either an interface slot for a LAN card (e.g. Agilent G1369B/C LAN Interface) or they have an on-board LAN interface (e.g. detectors G1315C/D DAD and G1365C/D MWD). This interface allows the control of the module/system via a PC with the appropriate control software. Some modules have neither on-board LAN nor an interface slot for a LAN card (e.g. G1170A Valve Drive or G4227A Flexible Cube). These are hosted modules and require a Host module with firmware B.06.40 or later or with additional G1369C LAN Card.

NOTE

LAN connection is made between at least one of the Agilent modules and the Control PC.

- If an Assist Hub is installed, connect the LAN to the Lab LAN port of this module.
- If an Assist Hub is NOT installed and a detector (DAD/MWD/FLD/VWD/RID) is installed, connect the LAN to this module.
- If an Assist Hub is NOT installed and there are multiple detectors with spectral capabilities, consider using additional LAN connections for each detector.
- If an Assist Hub is installed, connect additional LAN connections from the detectors and pumps to the Assist Hub.

USB

The USB interface replaces the RS-232 Serial interface in new generation modules. For details on USB refer to **USB (Universal Serial Bus)** on page 215.

Analog Signal Output

The analog signal output can be distributed to a recording device. For details refer to the description of the module's mainboard.

Remote (ERI)

The ERI (Enhanced Remote Interface) connector may be used in combination with other analytical instruments from Agilent Technologies if you want to use features as common shut down, prepare, and so on.

It allows easy connection between single instruments or systems to ensure coordinated analysis with simple coupling requirements.

The subminiature D connector is used. The module provides one remote connector which is inputs/outputs (wired- or technique).

To provide maximum safety within a distributed analysis system, one line is dedicated to **SHUT DOWN** the system's critical parts in case any module detects a serious problem. To detect whether all participating modules are switched on or properly powered, one line is defined to summarize the **POWER ON** state of all connected modules. Control of analysis is maintained by signal readiness **READY**

for next analysis, followed by **START** of run and optional **STOP** of run triggered on the respective lines. In addition **PREPARE** and **START REQUEST** may be issued. The signal levels are defined as:

- standard TTL levels (0 V is logic true, + 5.0 V is false),
- fan-out is 10,
- input load is 2.2 kOhm against + 5.0 V, and
- output are open collector type, inputs/outputs (wired- or technique).

NOTE

All common TTL circuits operate with a 5 V power supply. A TTL signal is defined as "low" or L when between 0 V and 0.8 V and "high" or H when between 2.0 V and 5.0 V (with respect to the ground terminal).

Table 21: ERI signal distribution

Pin	Signal	Description
1	START REQUEST	(L) Request to start injection cycle (for example, by start key on any module). Receiver is the autosampler.
2	STOP	(L) Request to reach system ready state as soon as possible (for example, stop run, abort or finish and stop injection). Receiver is any module performing run-time controlled activities.
3	READY	(H) System is ready for next analysis. Receiver is any sequence controller.
4	POWER ON	(H) All modules connected to system are switched on. Receiver is any module relying on operation of others.
5		Not used
6	SHUT DOWN	(L) System has serious problem (for example, leak: stops pump). Receiver is any module capable to reduce safety risk.
7	START	(L) Request to start run / timetable. Receiver is any module performing run-time controlled activities.
8	PREPARE	(L) Request to prepare for analysis (for example, calibration, detector lamp on). Receiver is any module performing pre-analysis activities.

Special Interfaces

There is no special interface for this module.

ERI (Enhanced Remote Interface)

ERI replaces the AGP Remote Interface that is used in the HP 1090/1040/1050/1100 HPLC systems and Agilent 1100/1200/1200 Infinity HPLC modules. All new InfinityLab LC Series products using the communication board core electronics use ERI. This interface is already used in the Agilent Universal Interface Box 2 (UIB2)

ERI Description

The ERI interface contains eight individual programmable input/output pins. In addition, it provides 24 V power and 5 V power and a serial data line to detect and recognize further add-ons that could be connected to this interface. This way the interface can support various additional devices like sensors, triggers (in and out) and small controllers, etc.

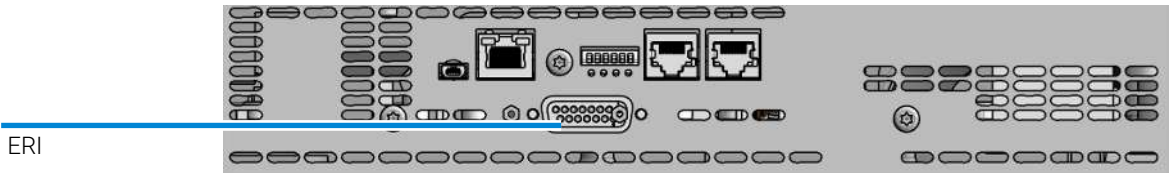


Figure 49: Location of the ERI interface

	Pin	Enhanced Remote
<p>D-Sub female 15way user's view to connector</p> <p>101 102 103 104 105 106 107 108 1 2 3 4 5 6 7 8 9 10 11 12 13 14 15 1Wireprom DGND +5V PGND PGND +24V +24V</p>	1	IO 1 (START REQUEST)
	2	IO 2 (STOP)
	3	IO 3 (READY)
	4	IO 4 (POWER ON)
	5	IO 5 (NOT USED)
	6	IO 6 (SHUT DOWN)
	7	IO 7 (START)
	8	IO 8 (PREPARE)
	9	1 wire DATA
	10	DGND
	11	+5 V ERI out
	12	PGND

	Pin	Enhanced Remote
	13	PGND
	14	+24 V ERI out
	15	+24 V ERI out

IO (Input/Output) Lines

- Eight generic bi-directional channels (input or output).
- Same as the APG Remote.
- Devices like valves, relays, ADCs, DACs, controllers can be supported/ controlled.

1-Wire Data (Future Use)

This serial line can be used to read out an EPROM or write into an EPROM of a connected ERI-device. The firmware can detect the connected type of device automatically and update information in the device (if required).

5V Distribution (Future Use)

- Available directly after turning on the hosting module (assures that the firmware can detect certain basic functionality of the device).
- For digital circuits or similar.
- Provides 500 mA maximum.
- Short-circuit proof with automatic switch off (by firmware).

24V Distribution (Future Use)

- Available by firmware command (defined turn on/off).
- For devices that need higher power
 - Class 0: 0.5 A maximum (12 W)
 - Class 1: 1.0 A maximum (24 W)
 - Class 2: 2.0 A maximum (48 W)
- Class depends on hosting module’s internal power overhead.

- If a connected device requires more power the firmware detects this (overcurrent detection) and provides the information to the user interface.
- Fuse used for safety protection (on board).
- Short circuit will be detected through hardware.

USB (Universal Serial Bus)

USB (Universal Serial Bus) - replaces RS232, supports:

- a PC with control software (for example Agilent Lab Advisor)
- USB Flash Disk

Instrument Layout

The industrial design of the module incorporates several innovative features. It uses Agilent's E-PAC concept for the packaging of electronics and mechanical assemblies. This concept is based upon the use of expanded polypropylene (EPP) layers of foam plastic spacers in which the mechanical and electronic boards components of the module are placed. This pack is then housed in a metal inner cabinet which is enclosed by a plastic external cabinet. The advantages of this packaging technology are:

- virtual elimination of fixing screws, bolts or ties, reducing the number of components and increasing the speed of assembly/disassembly,
- the plastic layers have air channels molded into them so that cooling air can be guided exactly to the required locations,
- the plastic layers help cushion the electronic and mechanical parts from physical shock, and
- the metal inner cabinet shields the internal electronics from electromagnetic interference and also helps to reduce or eliminate radio frequency emissions from the instrument itself.

Early Maintenance Feedback (EMF)

Maintenance requires the exchange of components that are subject to wear or stress. Ideally, the frequency at which components are exchanged should be based on the intensity of use of the module and the analytical conditions, and not on a predefined time interval. The early maintenance feedback (EMF) feature monitors the use of specific components in the instrument, and provides feedback when the user-selectable limits have been exceeded. The visual feedback in the user interface provides an indication that maintenance procedures should be scheduled.

EMF Counters

EMF counters increment with use and can be assigned a maximum limit which provides visual feedback in the user interface when the limit is exceeded. Some counters can be reset to zero after the required maintenance procedure.

Lamp Type	Counter Reset	Comment
Lamp with RFID tag	NO	
Lamp without RFID tag	YES	Via LabAdvisor or Instant Pilot

The detector provides the following EMF counters:

- Deuterium Lamp On-Time
- Number of UV lamp ignitions

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 needs 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.

Module-Specific Hardware Information

Setting the 6-bit Configuration Switch

The 6-bit configuration switch is located at the rear of the module with communication board electronics. Switch settings provide configuration parameters for LAN and instrument specific initialization procedures.

All modules with communication board electronics:

- Default is ALL switches DOWN (best settings).
 - Default IP address for LAN 192.168.254.11
- For specific LAN modes switches 4-5 must be set as required.
- For boot resident/cold start modes switches 1+2 or 6 must be UP.

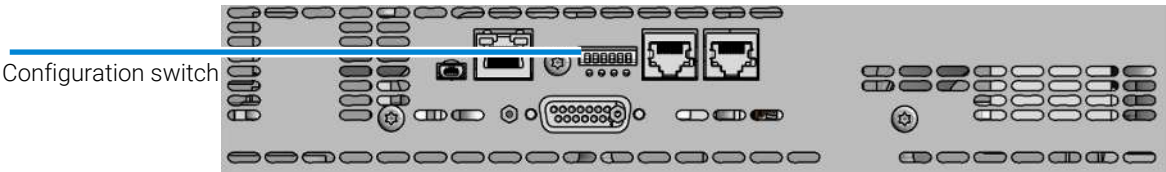


Figure 50: Location of configuration switch

Table 22: 6-bit configuration switch

SW1	SW2	SW3	SW4	SW5	SW6	Mode	Init Mode
0	0	0	0	0	0	COM	Use Default IP Address (192.168.254.11, Subnet mask: 255.255.255.0)
0	0	0	0	1	0	COM	Use Stored IP Address
0	0	0	1	0	0	COM	USE DHCP to request IP Address (Host name will be the MAC address)
1	0	0	0	0	0	Test	Boot Main System/Keep Data
1	1	0	0	0	0	Test	Boot Resident System/Keep Data

SW1	SW2	SW3	SW4	SW5	SW6	Mode	Init Mode
1	0	0	0	0	1	Test	Boot Main System/Revert to Default Data
1	1	0	0	0	1	Test	Boot Resident System/Revert to Default Data

Legend:

0 (switch down), 1 (switch up), SW (switch)

Special Settings

Boot-Resident/Main

Firmware update procedures may require this mode in case of firmware loading errors (main/resident firmware part).

If you use the following switch settings and power the instrument up again, the instrument firmware stays in the resident/main mode. In resident mode, it is not operable as a module. It only uses basic functions of the operating system for example, for communication. In this mode the main firmware can be loaded (using update utilities).

Forced Cold Start

A forced cold start can be used to bring the module into a defined mode with default parameter settings.

- Boot Main System / Revert to Default Data
The instrument will boot to main mode and changes to the module’s default parameter. May be also required to load resident firmware into the module.
- Boot Resident System / Revert to Default Data
The instrument will boot to resident mode and changes to the module’s default parameter. May be also required to load main firmware into the module.

CAUTION

Loss of data
Forced cold start erases all methods and data stored in the non-volatile memory. Exceptions are calibration settings, diagnosis and repair log books which will not be erased.

— Save your methods and data before executing a forced cold start.

Early Maintenance Feedback (EMF)

Maintenance requires the exchange of components which are subject to wear or stress. Ideally, the frequency at which components are exchanged should be based on the intensity of usage of the module and the analytical conditions, and not on a predefined time interval. The early maintenance feedback (EMF) feature monitors the usage of specific components in the instrument, and provides feedback when the user-selectable limits have been exceeded. The visual feedback in the user interface provides an indication that maintenance procedures should be scheduled.

EMF Counters

EMF counters increment with use and can be assigned a maximum limit which provides visual feedback in the user interface when the limit is exceeded. Some counters can be reset to zero after the required maintenance procedure.

Lamp Type	Counter Reset	Comment
lamp with RFID tag	NO	
lamp without RFID tag	YES	via Lab Advisor or Instant Pilot

The detector provides the following EMF counters:

- Deuterium Lamp On-Time
- Number of UV lamp ignitions

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 needs 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.

NOTE

This function is only available via Agilent Lab Advisor or Instant Pilot.

12

LAN Configuration

This chapter provides information on connecting the module to the control software.

What You Have to Do First 223

TCP/IP Parameter Configuration 224

Configuration Switch 225

Initialization Mode Selection 226

Dynamic Host Configuration Protocol (DHCP) 228

General Information (DHCP) 228

Setup (DHCP) 228

Manual Configuration 231

With Telnet 231

With the Instant Pilot (G4208A) 234

PC and User Interface Software Setup 236

PC Setup for Local Configuration 236

What You Have to Do First

The module has an on-board LAN communication interface.

NOTE

This chapter is generic and may show figures that differ from your module. The functionality is the same.

- 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 of the module underneath the configuration switch (see [Figure 52](#) on page 223).



Part number of the detector mainboard
Revision Code, Vendor, Year and Week of assembly
MAC address
Country of Origin

Figure 51: 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.

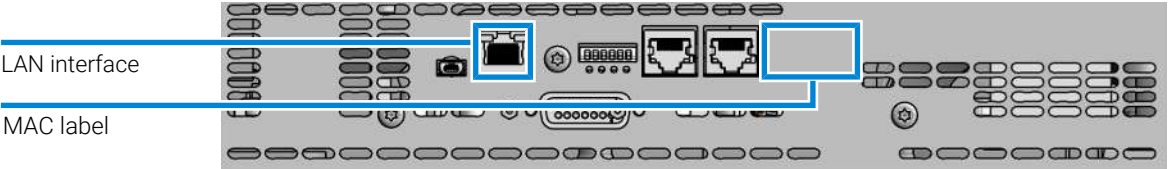


Figure 52: Location of LAN interfaces and MAC label

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 228
- by manually setting the parameters using Telnet
- by manually setting the parameters using the Local Controller

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 non-volatile memory or initialized with known default values. The initialization mode is selected by the configuration switch, see [Table 23](#) on page 226.

Configuration Switch

The configuration switch can be accessed at the rear of the module.

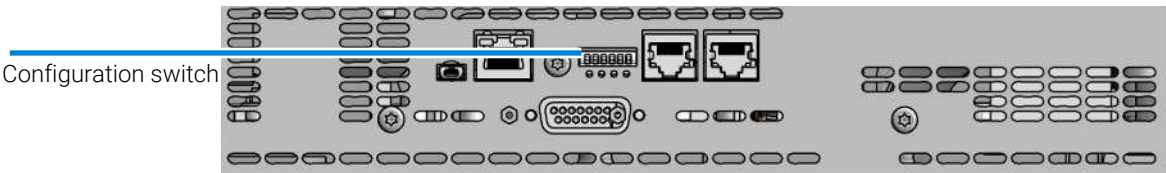


Figure 53: Location of configuration switch

The module is shipped with all switches set to OFF, as shown above.


NOTE

To perform any LAN configuration, SW1 and SW2 must be set to OFF.

Initialization Mode Selection

The following initialization (init) modes are selectable:

Table 23: Initialization mode switches

	SW1	SW2	SW3	SW4	SW5	SW6	Init Mode
	0	0	0	0	0	0	Use Default IP Address
	0	0	0	0	1	0	Use Stored IP Address
	0	0	0	1	0	0	Use DHCP

Note: The setting '0' (down) is essential.

Legend:

0 (switch down), 1 (switch up), SW (switch)
Default IP address for LAN is 192.168.254.11.
DHCP address is the module’s LAN MAC address.

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.

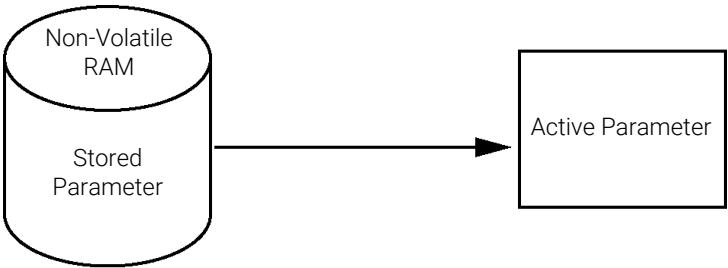


Figure 54: Using Stored (principle)

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 24](#) on page 227.

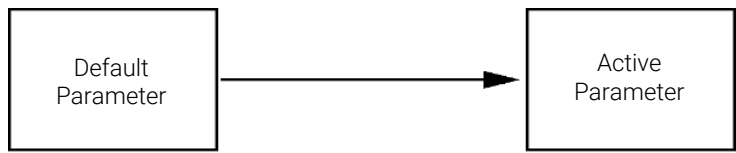


Figure 55: Using Default (principle)

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 24: 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**.

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).

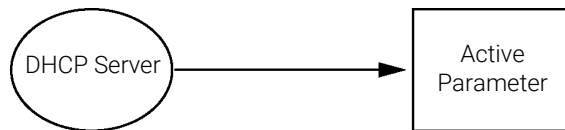


Figure 56: DHCP (principle)

NOTE

- It may take some time until the DHCP server has updated the DNS server with the hostname information.
- It may be necessary to fully qualify the hostname with the DNS suffix, e.g. 0030d3177321.country.company.com.
- 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 mainboard). This MAC address is on a label on the card or at the rear of the mainboard, for example, 0030d3177321.

On the Local Controller the MAC address can be found under **Details** in the LAN section.

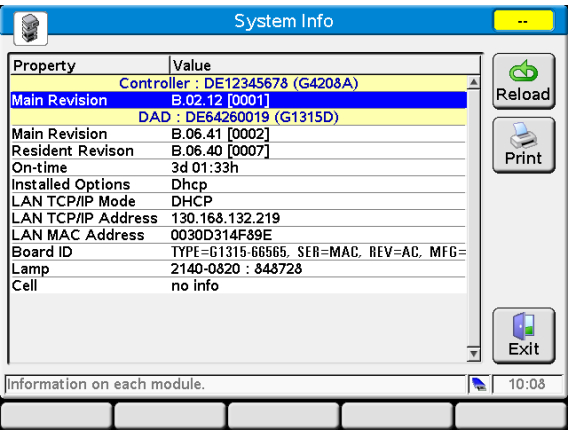


Figure 57: LAN setting on Instant Pilot

- 2 Set the configuration switch to DHCP either on the G1369C LAN Interface Card or the mainboard of above mentioned modules.

Table 25: 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 26: 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 228).

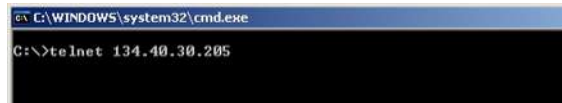
Manual Configuration

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.

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>



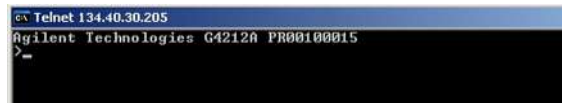
```

C:\WINDOWS\system32\cmd.exe
C:\>telnet 134.40.30.205
  
```

Figure 58: 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 Switch](#) on page 225).

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

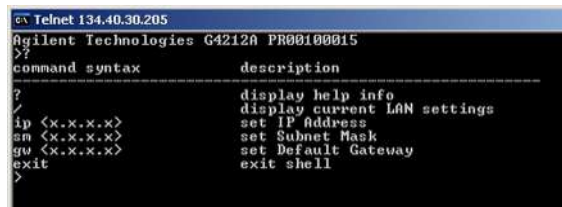


```

c\ Telnet 134.40.30.205
Agilent Technologies G4212A PR00100015
>_
  
```

Figure 59: A connection to the module is made

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



```


c\ Telnet 134.40.30.205
Agilent Technologies G4212A PR00100015
>?
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 60: Telnet commands

Table 27: 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 134.40.28.56`
- 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.



information about the LAN interface
MAC address, initialization mode
Initialization mode is Using Stored
active TCP/IP settings

Telnet - Current settings in "Using Stored" mode

TCP/IP status - here ready
connected to PC with controller software
(e.g. Agilent ChemStation), here not
connected

- 6 Change the IP address (in this example 192.168.254.12) and type “/” to list current settings.

```

c:\ Telnet 134.40.30.205
>ip 192.168.254.12
>
LAN Status Page
-----
MAC Address   : 0030D317521C
Init Mode    : Using Stored
TCP/IP Properties
- active -
IP Address   : 134.40.30.205
Subnet Mask  : 255.255.248.0
Def. Gateway : 134.40.24.1
- stored -
IP Address   : 192.168.254.12
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

Telnet - Change IP settings

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

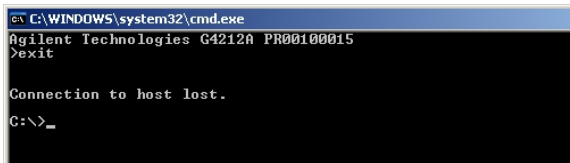


Figure 61: Closing the Telnet session

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 192.168.254.12.

With the Instant Pilot (G4208A)

To configure the TCP/IP parameters before connecting the module to the network, the Instant Pilot (G4208A) can be used.

- 1 From the Welcome screen press the **More** button.
- 2 Select **Configure**.
- 3 Press the module button of the module that hosts the LAN interface (usually the detector).
- 4 Scroll down to the LAN settings.

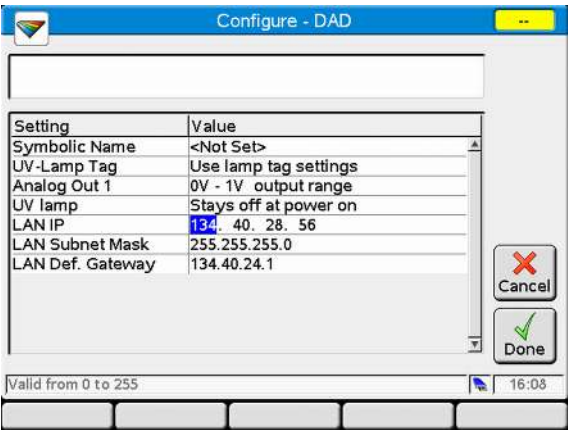


Figure 62: Instant Pilot - LAN configuration (edit mode)

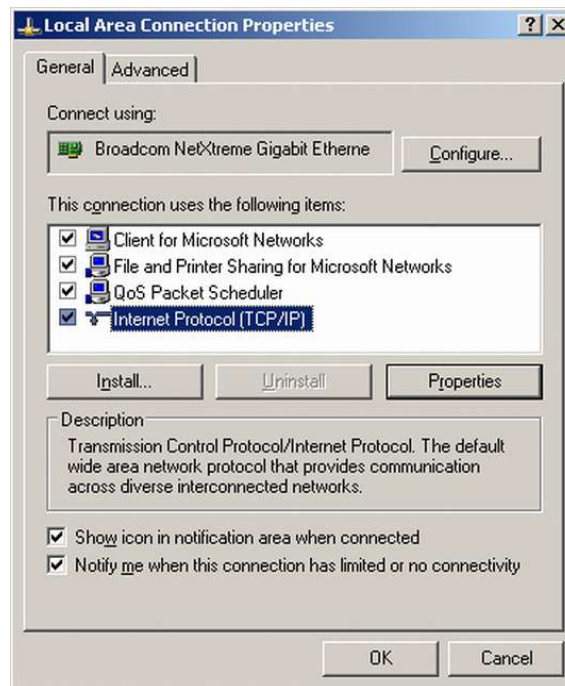
- 5 Press the **Edit** button (only visible if not in Edit mode), perform the required changes and press the **Done** button.
- 6 Leave the screen by clicking **Exit**.

PC and User Interface Software Setup

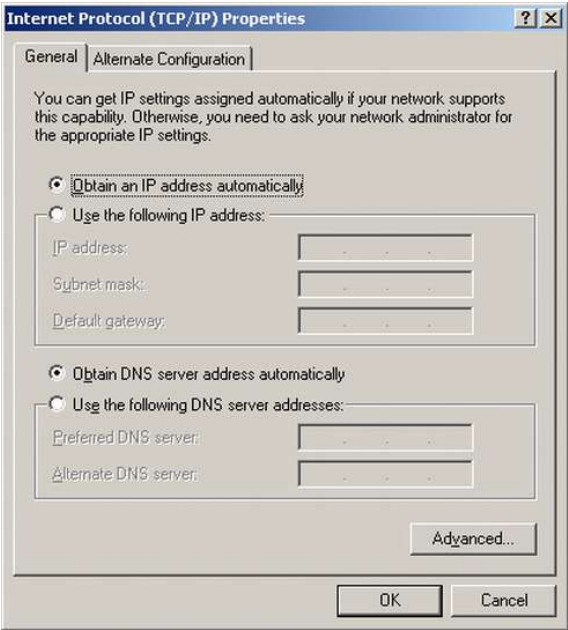
PC Setup for Local Configuration

This procedure describes the change of the TCP/IP settings on your PC to match the module's default parameters in a local configuration (see [Table 24](#) on page 227).

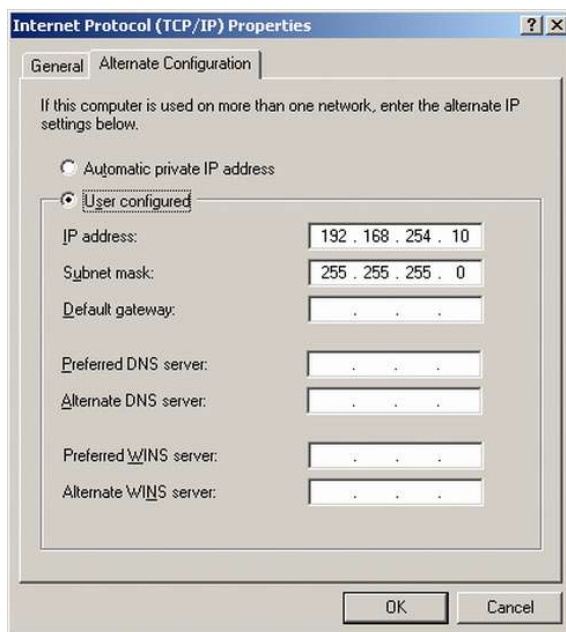
- 1 Open the Local Area Connection Properties and select **Internet Protocol (TCP/IP)**. Then click on **Properties**.



- 2 You may enter here the fixed IP address of the module or use the Alternative Configuration.



- 3 We will use the direct LAN access via Cross-over LAN cable with the module's IP address.



- 4 Click on **OK** to save the configuration.

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

General Safety Information 240

Safety Standards 240

General 240

Before Applying Power 241

Ground the Instrument 241

Do Not Operate in an Explosive Atmosphere 242

Do Not Remove the Instrument Cover 242

Do Not Modify the Instrument 242

In Case of Damage 242

Solvent Information 243

Magnets 245

Safety Symbols 245

Material Information 248

General Information About Solvent/Material Compatibility 248

At-a-Glance Details About Agilent Capillaries 254**Waste Electrical and Electronic Equipment (WEEE) Directive 258****Radio Interference 259****RFID Statement 260****Sound Emission 262****Declaration of Conformity for HOX2 Filter 263****Agilent Technologies on Internet 265**

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.

General

Do not use this product in any manner not specified by the manufacturer. The protective features of this product may be impaired if it is used in a manner not specified in the operation instructions.

Before Applying Power

WARNING

Wrong voltage range, frequency or cabling

Personal injury or damage to the instrument

- Verify that the voltage range and frequency of your power distribution matches to the power specification of the individual instrument.
- Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.
- Make all connections to the unit before applying power.

WARNING

Use of unsupplied cables

Using cables not 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.

NOTE

Note the instrument's external markings described under [Safety Symbols](#) on page 245.

Ground the Instrument

WARNING

Missing electrical ground

Electrical shock

- If your product is provided with a grounding type power plug, the instrument chassis and cover must be connected to an electrical ground to minimize shock hazard.
- The ground pin must be firmly connected to an electrical ground (safety ground) terminal at the power outlet. Any interruption of the protective (grounding) conductor or disconnection of the protective earth terminal will cause a potential shock hazard that could result in personal injury.

Do Not Operate in an Explosive Atmosphere

WARNING

Presence of flammable gases or fumes

Explosion hazard

- Do not operate the instrument in the presence of flammable gases or fumes.
-

Do Not Remove the Instrument Cover

WARNING

Instrument covers removed

Electrical shock

- Do Not Remove the Instrument Cover
 - Only Agilent authorized personnel are allowed to remove instrument covers. Always disconnect the power cables and any external circuits before removing the instrument cover.
-

Do Not Modify the Instrument

Do not install substitute parts or perform any unauthorized modification to the product. Return the product to an Agilent Sales and Service Office for service and repair to ensure that safety features are maintained.

In Case of Damage

WARNING

Damage to the module

Personal injury (for example electrical shock, intoxication)

- Instruments that appear damaged or defective should be made inoperative and secured against unintended operation until they can be repaired by qualified service personnel.
-

Solvent Information

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.
- Do not use solvents with an auto-ignition temperature below 200 °C (392 °F). Do not use solvents with a boiling point below 56 °C (133 °F).
- Avoid high vapor concentrations. Keep the solvent temperature at least 40 °C (72 °F) below the boiling point of the solvent used. This includes the solvent temperature in the sample compartment. For the solvents methanol and ethanol keep the solvent temperature at least 25 °C (45 °F) below the boiling point.
- Do not operate the instrument in an explosive atmosphere.
- Do not use solvents of ignition Class IIC according IEC 60079-20-1 (for example, carbon disulfide).
- Reduce the volume of substances to the minimum required for the analysis.
- Never exceed the maximum permissible volume of solvents (8 L) in the solvent cabinet. Do not use bottles that exceed the maximum permissible volume as specified in the usage guideline for solvent cabinet.
- Ground the waste container.
- Regularly check the filling level of the waste container. The residual free volume in the waste container must be large enough to collect the waste liquid.
- To achieve maximal safety, regularly check the tubing for correct installation.

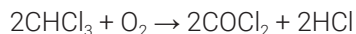
NOTE

For details, see the usage guideline for the solvent cabinet. A printed copy of the guideline has been shipped with the solvent cabinet, electronic copies are available in the Agilent Information Center or via the Internet.

Recommendations on the Use of Solvents

Observe the following recommendations on the use of solvents.

- Brown glass ware can avoid growth of algae.
- Follow the recommendations for avoiding the growth of algae, see the pump manuals.
- Small particles can permanently block capillaries and valves. Therefore, always filter solvents through 0.22 µm filters.
- Avoid or minimize the use of solvents that may corrode parts in the flow path. Consider specifications for the pH range given for different materials such as flow cells, valve materials etc. and recommendations in subsequent sections.
- 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 sulfuric acid and nitric acid, especially at higher temperatures (if your chromatography method allows, replace 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,

- chromatographic grade ethers, which can contain peroxides (for example, THF, dioxane, diisopropyl ether) should be filtered through dry aluminium oxide which adsorbs the peroxides,
- solvents containing strong complexing agents (e.g. EDTA),
- mixtures of carbon tetrachloride with 2-propanol or THF.
- Avoid the use of dimethyl formamide (DMF). Polyvinylidene fluoride (PVDF), which is used in leak sensors, is not resistant to DMF.

Flow cell

To protect optimal functionality of your flow-cell:

- Avoid the use of alkaline solutions (pH > 9.5) which can attack quartz and thus impair the optical properties of the flow cell.

Magnets

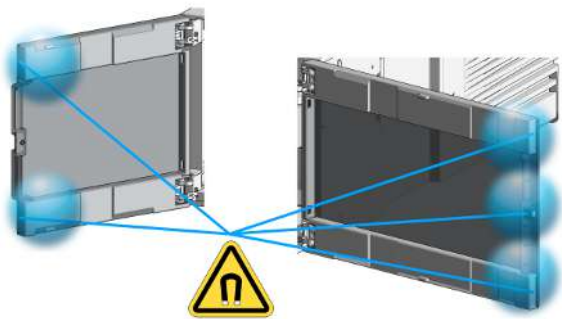


Figure 63: Magnets in doors of pumps, autosamplers, detectors, and fraction collectors

Safety Symbols

Table 28: Symbols

	The apparatus is marked with this symbol when the user shall 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.

Appendix

General Safety Information



The apparatus is marked with this symbol when hot surfaces are available and the user should not touch it when heated up.



Indicates flammable material used. Consult the Agilent Information Center / User Manual before attempting to install or service this equipment. Follow all safety precautions.



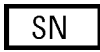
Confirms that a manufactured product complies with all applicable European Community directives. The European Declaration of Conformity is available at: <http://regulations.corporate.agilent.com/DoC/search.htm>



Manufacturing date.



Product Number



Serial Number



Power symbol indicates On/Off.

The apparatus is not completely disconnected from the mains supply when the on/off switch is in the Off position



Pacemaker

Magnets could affect the functioning of pacemakers and implanted heart defibrillators. A pacemaker could switch into test mode and cause illness. A heart defibrillator may stop working. If you wear these devices keep at least 55 mm distance to magnets. Warn others who wear these devices from getting too close to magnets.



Magnetic field

Magnets produce a far-reaching, strong magnetic field. They could damage TVs and laptops, computer hard drives, credit and ATM cards, data storage media, mechanical watches, hearing aids and speakers. Keep magnets at least 25 mm away from devices and objects that could be damaged by strong magnetic fields.



Indicates a pinching or crushing hazard



Indicates a piercing or cutting hazard.

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.
-

Material Information

This section provides detailed information about materials used in the HPLC system and general information about solvent/material compatibility.

General Information About Solvent/Material Compatibility

Materials in the flow path are carefully selected based on Agilent's experiences in developing highest-quality instruments for HPLC analysis over several decades. These materials exhibit excellent robustness under typical HPLC conditions. For any special condition, please consult the material information section or contact Agilent.

Disclaimer

Subsequent data was collected from external resources and is meant as a reference. Agilent cannot guarantee the correctness and completeness of such information. Data is based on compatibility libraries, which are not specific for estimating the long-term life time under specific but highly variable conditions of UHPLC systems, solvents, solvent mixtures, and samples. Information also cannot be generalized due to catalytic effects of impurities like metal ions, complexing agents, oxygen etc. Apart from pure chemical corrosion, other effects like electro corrosion, electrostatic charging (especially for nonconductive organic solvents), swelling of polymer parts etc. need to be considered. Most data available refers to room temperature (typically 20 – 25 °C, 68 – 77 °F). If corrosion is possible, it usually accelerates at higher temperatures. If in doubt, please consult technical literature on chemical compatibility of materials.

MP35N

MP35N is a nonmagnetic, nickel-cobalt-chromium-molybdenum alloy demonstrating excellent corrosion resistance (for example, against nitric and sulfuric acids, sodium hydroxide, and seawater) over a wide range of concentrations and temperatures. In addition, this alloy shows exceptional

resistance to high-temperature oxidation. Due to excellent chemical resistance and toughness, the alloy is used in diverse applications: dental products, medical devices, nonmagnetic electrical components, chemical and food processing equipment, marine equipment. Treatment of MP35N alloy samples with 10 % NaCl in HCl (pH 2.0) does not reveal any detectable corrosion. MP35N also demonstrates excellent corrosion resistance in a humid environment. Although the influence of a broad variety of solvents and conditions has been tested, users should keep in mind that multiple factors can affect corrosion rates, such as temperature, concentration, pH, impurities, stress, surface finish, and dissimilar metal contacts.

Polyphenylene Sulfide (PPS)

Polyphenylene sulfide has outstanding stability even at elevated temperatures. It is resistant to dilute solutions of most inorganic acids, but it can be attacked by some organic compounds and oxidizing reagents. Nonoxidizing inorganic acids, such as sulfuric acid and phosphoric acid, have little effect on polyphenylene sulfide, but at high concentrations and temperatures, they can still cause material damage. Nonoxidizing organic chemicals generally have little effect on polyphenylene sulfide stability, but amines, aromatic compounds, and halogenated compounds may cause some swelling and softening over extended periods of time at elevated temperatures. Strong oxidizing acids, such as nitric acid (> 0.1 %), hydrogen halides (> 0.1 %), peroxy acids (> 1 %), or chlorosulfuric acid degrade polyphenylene sulfide. It is not recommended to use polyphenylene sulfide with oxidizing material, such as sodium hypochlorite and hydrogen peroxide. However, under mild environmental conditions, at low concentrations and for short exposure times, polyphenylene sulfide can withstand these chemicals, for example, as ingredients of common disinfectant solutions.

PEEK

PEEK (Polyether-Ether Ketones) combines excellent properties regarding biocompatibility, chemical resistance, mechanical and thermal stability. PEEK is therefore the material of choice for UHPLC and biochemical instrumentation.

It is stable in the specified pH range (for the Bio-Inert LC system: pH 1 – 13 , see bio-inert module manuals for details), and inert to many common solvents.

There are still some known incompatibilities with chemicals such as chloroform, methylene chloride, THF, DMSO, strong acids (nitric acid > 10 %, sulfuric acid > 10 %, sulfonic acids, trichloroacetic acid), halogens or aqueous halogen solutions, phenol and derivatives (cresols, salicylic acid, and so on).

When used above room temperature, PEEK is sensitive to bases and various organic solvents, which can cause it to swell. Under such conditions, normal PEEK capillaries are sensitive to high pressure. Therefore, Agilent uses stainless steel clad PEEK capillaries in bio-inert systems. The use of stainless steel clad PEEK capillaries keeps the flow path free of steel and ensures pressure stability up to 600 bar. If in doubt, consult the available literature about the chemical compatibility of PEEK.

Polyimide

Agilent uses semi-crystalline polyimide for rotor seals in valves and needle seats in autosamplers. One supplier of polyimide is DuPont, which brands polyimide as Vespel, which is also used by Agilent.

Polyimide is stable in a pH range between 1 and 10 and in most organic solvents. It is incompatible with concentrated mineral acids (e.g. sulphuric acid), glacial acetic acid, DMSO and THF. It is also degraded by nucleophilic substances like ammonia (e.g. ammonium salts in basic conditions) or acetates.

Polyethylene (PE)

Agilent uses UHMW (ultra-high molecular weight)-PE/PTFE blends for yellow piston and wash seals, which are used in 1290 Infinity pumps, 1290 Infinity II/III pumps, the G7104C and for normal phase applications in 1260 Infinity pumps.

Polyethylene has a good stability for most common inorganic solvents including acids and bases in a pH range of 1 to 12.5. It is compatible with many organic solvents used in chromatographic systems like methanol, acetonitrile and isopropanol. It has limited stability with aliphatic, aromatic and halogenated hydrocarbons, THF, phenol and derivatives, concentrated acids and bases. For normal phase applications, the maximum pressure should be limited to 200 bar.

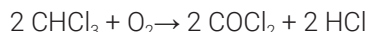
Tantalum (Ta)

Tantalum is inert to most common HPLC solvents and almost all acids except fluoric acid and acids with free sulfur trioxide. It can be corroded by strong bases (e.g. hydroxide solutions > 10 %, diethylamine). It is not recommended for the use with fluoric acid and fluorides.

Stainless Steel (SST)

Stainless steel is inert against many common solvents. It is stable in the presence of acids and bases in a pH range of 1 to 12.5. It can be corroded by acids below pH 2.3. It can also corrode in following solvents:

- Solutions of alkali halides, their respective acids (for example, lithium iodide, potassium chloride) and aqueous solutions of halogens.
- High concentrations of inorganic acids like nitric acid, sulfuric acid, and organic solvents 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.

- Chromatographic grade ethers, which can contain peroxides (for example, THF, dioxane, diisopropyl ether). Such ethers should be filtered through dry aluminum 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, ethylenediaminetetraacetic acid).
- Mixtures of carbon tetrachloride with isopropanol or THF.

Titanium (Ti)

Titanium is highly resistant to oxidizing acids (for example, nitric, perchloric and hypochlorous acid) over a wide range of concentrations and temperatures. This is due to a thin oxide layer on the surface, which is stabilized by oxidizing compounds. Non-oxidizing acids (for example, hydrochloric, sulfuric and phosphoric acid) can cause slight corrosion, which increases with acid concentration and temperature. For example, the corrosion rate with 3 % HCl (about pH 0.1) at room temperature is about 13 $\mu\text{m}/\text{year}$. At room temperature, titanium is resistant to concentrations of about 5 % sulfuric acid (about pH 0.3). Addition of nitric acid to hydrochloric or sulfuric acids significantly reduces corrosion rates. Titanium is sensitive to acidic metal chlorides like FeCl_3 or CuCl_2 .

Titanium is subject to corrosion in anhydrous methanol, which can be avoided by adding a small amount of water (about 3 %). Slight corrosion is possible with ammonia > 10 %.

Diamond-Like Carbon (DLC)

Diamond-Like Carbon is inert to almost all common acids, bases, and solvents. There are no documented incompatibilities for HPLC applications.

Fused Silica and Quartz (SiO₂)

Fused silica is used in Max Light Cartridges. Quartz is used for classical flow cell windows. It is inert against all common solvents and acids except hydrofluoric acid and acidic solvents containing fluorides. It is corroded by strong bases and should not be used above pH 12 at room temperature. The corrosion of flow cell windows can negatively affect measurement results. For a pH greater than 12, the use of flow cells with sapphire windows is recommended.

Gold

Gold is inert to all common HPLC solvents, acids, and bases within the specified pH range. It can be corroded by complexing cyanides and concentrated acids like aqua regia.

Zirconium Oxide (ZrO₂)

Zirconium Oxide is inert to almost all common acids, bases, and solvents. There are no documented incompatibilities for HPLC applications.

Platinum/Iridium

Platinum/Iridium is inert to almost all common acids, bases, and solvents. There are no documented incompatibilities for HPLC applications.

Fluorinated Polymers (PTFE, PFA, FEP, FFKM, PVDF)

Fluorinated polymers like PTFE (polytetrafluorethylene), PFA (perfluoroalkoxy), and FEP (fluorinated ethylene propylene) are inert to almost all common acids, bases, and solvents. FFKM is perfluorinated rubber, which is also resistant to most chemicals. As an elastomer, it may swell in some organic solvents like halogenated hydrocarbons.

TFE/PDD copolymer tubings, which are used in all Agilent degassers except G1322A/G7122A, are not compatible with fluorinated solvents like Freon, Fluorinert, or Vertrel. They have limited life time in the presence of hexafluoroisopropanol (HFIP). To ensure the longest possible life with HFIP, it is best to dedicate a particular chamber to this solvent, not to switch solvents, and not to let dry out the chamber. For optimizing the life of the pressure sensor, do not leave HFIP in the chamber when the unit is off.

The tubing of the leak sensor is made of PVDF (polyvinylidene fluoride), which is incompatible with the solvent DMF (dimethylformamide).

Sapphire, Ruby, and Al₂O₃-Based Ceramics

Sapphire, ruby, and ceramics based on aluminum oxide Al₂O₃ are inert to almost all common acids, bases, and solvents. There are no documented incompatibilities for HPLC applications.


At-a-Glance Details About Agilent Capillaries

The following section provides useful information about Agilent capillaries and its characteristics.

Syntax for capillary description

Type - Material - Capillary dimensions - Fitting Left/Fitting right

Table 29: Example for a capillary description



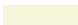










Code provided with the part	Meaing of the code
Color code: 	Material of the product is MP35N, the inner diameter is 0.20 or 0.25 mm
Capillary	The part is a connection capillary
MP35N	Material of the part is MP35N
0.25 x 80 mm	The part has an inner diameter of 0.25 mm and a length of 80 mm
SI/SI	Left fitting: Swagelok + 1.6 mm Port id, Intermediate Right fitting: Swagelok + 1.6 mm Port id, Intermediate

To get an overview of the code in use, see

- Color: [Table 30](#) on page 255
- Type: [Table 31](#) on page 255
- Material: [Table 32](#) on page 256
- Dimension: [Table 33](#) on page 256
- Fittings: [Table 34](#) on page 257

Color Coding Guide

Table 30: Color-coding key for Agilent capillary tubing

Internal diameter in mm		Color code	
0.015			Orange
0.025			Yellow
0.05			Beige
0.075			Black
0.075	MP35N		Black with orange stripe
0.1			Purple
0.12			Red
0.12	MP35N		Red with orange stripe
0.17			Green
0.17	MP35N		Green with orange stripe
0.20 /0.25			Blue
0.20 /0.25	MP35N		Blue with orange stripe
0.3			Grey
0.50			Bone White

NOTE

As you move to smaller-volume, high efficiency columns, you'll want to use narrow id tubing, as opposed to the wider id tubing used for conventional HPLC instruments.

Abbreviation Guide for Type

Table 31: Type (gives some indication on the primary function, like a loop or a connection capillary)

Key	Description
Capillary	Connection capillaries
Loop	Loop capillaries
Seat	Autosampler needle seats

Key	Description
Tube	Tubing
Heat exchanger	Heat exchanger

Abbreviation Guide for Material

Table 32: Material (indicates which raw material is used for the capillary)

Key	Description
ST	Stainless steel
Ti	Titanium
PK	PEEK
FS/PK	PEEK-coated fused silica ²
PK/ST	Stainless steel-coated PEEK ³
PFFE	PTFE
FS	Fused silica
MP35N	Nickel-cobalt-chromium-molybdenum alloy

Abbreviation Guide for Capillary Dimensions

Table 33: Capillary dimensions (indicates inner diameter (id), length, and volume of the capillary)

Description
id (mm) x Length (mm)
Volume (μL)

2 Fused silica in contact with solvent
3 Stainless steel-coated PEEK

Abbreviation Guide for Fitting Left/Fitting Right

Table 34: Fitting left/fitting right (indicates which fitting is used on both ends of the capillary)

Key	Description
W	Swagelok + 0.8 mm Port id
S	Swagelok + 1.6 mm Port id
M	Metric M4 + 0.8 mm Port id
E	Metric M3 + 1.6 mm Port id
U	Swagelok union
L	Long
X	Extra long
H	Long head
G	Small head SW 4
N	Small head SW 5
F	Finger-tight
V	1200 bar
B	Bio
P	PEEK
I	Intermediate

Waste Electrical and Electronic Equipment (WEEE) Directive

This product complies with the European WEEE Directive marking requirements. The affixed label indicates that you must not discard this electrical/electronic product in domestic household waste.

**NOTE**

Do not dispose of in domestic household waste

To return unwanted products, contact your local Agilent office, or see <https://www.agilent.com> for more information.

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.

RFID Statement

Brasil

Este equipamento não tem direito à proteção contra interferência prejudicial e não pode causar interferência em sistemas devidamente autorizados. Para mais informações, consulte o site da Anatel: <https://www.gov.br/anatel/pt-br>.

Este produto não é apropriado para uso em ambientes domésticos, pois poderá causar interferências eletromagnéticas que obrigam o usuário a tomar medidas necessárias para minimizar estas interferências.

Canada

Statement according to RSS GEN Issue 5:

This device contains licence-exempt transmitter(s)/receiver(s) that comply with Innovation, Science and Economic Development Canada's licence-exempt RSS(s). Operation is subject to the following two conditions:

1. This device may not cause interference
2. This device must accept any interference, including interference that may cause undesired operation of the device.

Cet appareil contient des émetteurs / récepteurs exemptés de licence conformes aux RSS (RSS) d'Innovation, Sciences et Développement économique Canada. Le fonctionnement est soumis aux deux conditions suivantes:

1. Cet appareil ne doit pas causer d'interférences
2. Cet appareil doit accepter toutes les interférences, y compris celles susceptibles de provoquer un fonctionnement indésirable de l'appareil.

Mexico

La operación de este equipo está sujeta a las siguientes dos condiciones:

1. es posible que este equipo o dispositivo no cause interferencia perjudicial y
2. este equipo o dispositivo debe aceptar cualquier interferencia, incluyendo la que pueda causar su operación no deseada.

Thailand

เครื่องโทรคมนาคมและอุปกรณ์นี้มีความสอดคล้องตามมาตรฐานหรือข้อกำหนดทางเทคนิคของ กสทช.
This telecommunication equipment conforms to NTC/NBTC technical requirement.

USA

- 1. User Information according to FCC 15.21:Changes or modifications not expressly approved by the party responsible for compliance could void the user’s authority to operate the equipment.
- 2. Part 15 Statement according to FCC 15.19:

This device complies with Part 15 of the FCC Rules. Operation is subject to the following two conditions:
 - This device may not cause interference.
 - This device must accept any interference, including interference that may cause undesired operation.

CAUTION

Do not change or modify the equipment.
Changes or modifications not expressly approved by Agilent could void your authority to operate the equipment.

NOTE

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules.
These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

Table 35: Operating frequencies and maximum power levels

Technology	Operating Frequencies/ Bands	Maximum Transmit Power Level
RFID	125 kHz	26.8 dBm

Sound Emission

Sound Pressure

Sound pressure $L_p < 70 \text{ dB(A)}$ according to DIN EN ISO 7779

Schalldruckpegel

Schalldruckpegel $L_p < 70 \text{ dB(A)}$ nach DIN EN ISO 7779

Declaration of Conformity for HOX2 Filter

Declaration of Conformity

We herewith inform you that the

Holmium Oxide Glass Filter

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.

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.

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.

Test wavelengths:

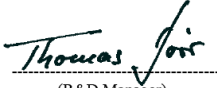
Where "x" can be any alphanumeric character

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			


*) The variation in Measured Wavelength depends on the different Optical Bandwidth.

28-Oct-2014


(Date)




(R&D Manager)



(Quality Manager)

P/N 89550-90501


Revision: G
Effective by: 28-Oct-2014



Agilent Technologies on Internet

For the latest information on products and services visit our worldwide web site on the Internet at:

<https://www.agilent.com>

In This Book

This manual contains technical reference information about the Agilent 1290 Infinity III Variable Wavelength Detector (G7114B) and the Agilent 1260 Infinity III Variable Wavelength Detector (G7114A).

The manual describes the following:

- introduction and specifications,
- installation,
- using and optimizing,
- troubleshooting and diagnose,
- maintenance and repair,
- parts identification,
- hardware information,
- safety and related information.

www.agilent.com

© Agilent Technologies Inc. 2014-2024
Edition: 10/2024

Document No: SD-29000240 Rev. E

