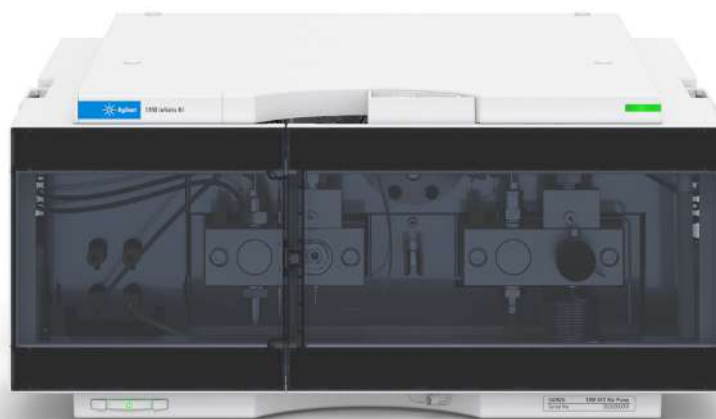




Agilent InfinityLab LC Series

1260 Infinity III SFC Binary Pump

User Manual



Notices

Document Information

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

Document No: SD-29000180 Rev. B
Edition: 10/2024

Copyright

© Agilent Technologies, Inc. 2017 - 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

In This Book 7

1 Introduction 8

- Introduction to the Pump 9
- Product Description of the 1260 Infinity III SFC Binary Pump (G4782A) 10
- Features of the 1260 Infinity III SFC Binary Pump (G4782A) 11
- Operating Principle 12

2 Site Requirements and Specifications 18

- Site Requirements 19
- Specifications of the 1260 Infinity III SFC Binary Pump (G4782A) 22

3 Installation 25

- Installing Capillaries 26
- Handling Leak and Waste 30
- Connecting Modules and Control Software 40

4 Using the Module 41

- General Information 42
- Best Practices 45
- Preparation of the System 51
- Preparation of the Module 60

5 Optimizing the Performance of the Module 64

- When to Use a Vacuum Degasser 65
- When to Use the Active Seal Wash Option 66
- How to Optimize the Compressibility Compensation Setting 67

6 Diagnostics and Troubleshooting 70

- Diagnostic Features 71
- Maintenance and Troubleshooting Tools of the Module 72
- Agilent Lab Advisor Software 74

7 Error Information 75

What Are Error Messages 77

General Error Messages 78

Pump Error Messages 90

8 Maintenance 110

Safety Information Related to Maintenance 112

Introduction to Maintenance 114

Overview of Maintenance and Simple Repairs 115

Maintenance Procedures 116

Cleaning the Module 117

Remove and Install Doors 118

Exchanging the Purge Valve Frit or the Purge Valve 121

Replace the O-Ring on the Purge Valve 124

Remove the Pump Head Assembly 127

Maintenance of a Pump Head without Seal Wash 129

Maintenance of a Pump Head with Seal Wash 134

Reinstall the Pump Head Assembly 140

Seal Wear-in Procedure 142

Exchange the Seal Wash Cartridge 143

Replace Leak Handling System Parts 145

Exchange the Inlet Valve 146

Exchange the Outlet Valve 148

Exchange the Solvent Selection Valve 150

Replace the Module Firmware 154

9 Parts and Materials for Maintenance 155

Accessory Kit G7111-68755 156

Pump Head Assembly Channel B 157

Pump Head Assembly Channel A 159

Outlet Valve 161

Purge Valve Assembly 162

Active Seal Wash Option 163

- HPLC System Tool Kit 164
- Solvent Cabinet 165
- Bottle Head Assembly 166
- Hydraulic Path with Solvent Selection Valve 167
- Cover Parts 169

10 Identifying Cables 170

- Cable Overview 171
- Analog Cables 173
- Remote Cables 175
- BCD Cables 179
- CAN/LAN Cables 181
- RS-232 Cables 182
- USB 183

11 Hardware Information 184

- General Hardware Information 185
- Module-Specific Hardware Information 199

12 LAN Configuration 202

- What You Have to Do First 203
- TCP/IP Parameter Configuration 204
- Configuration Switch 205
- Initialization Mode Selection 206
- Dynamic Host Configuration Protocol (DHCP) 208
- Manual Configuration 211
- PC and User Interface Software Setup 215

13 Appendix 218

- General Safety Information 219
- Material Information 227
- At-a-Glance Details About Agilent Capillaries 234
- Waste Electrical and Electronic Equipment (WEEE) Directive 238
- Radio Interference 239

Sound Emission 240

Agilent Technologies on Internet 241



In This Book

This manual covers the Agilent InfinityLab LC Series 1260 Infinity III SFC Binary Pump (G4782A).

1

Introduction

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

Introduction to the Pump 9

Product Description of the 1260 Infinity III SFC Binary Pump (G4782A) 10

Features of the 1260 Infinity III SFC Binary Pump (G4782A) 11

Operating Principle 12

Principle of Operation 12

Overview of the Hydraulic Path 13

Introduction to the Pump

The binary pump comprises two identical pumps integrated into one housing. Binary gradients are created by high-pressure mixing. A built-in degasser is available for applications that require best flow stability, especially at low flow rates, for maximum detector sensitivity. The pump is capable of delivering flow in the range of 0.1 – 5 mL/min against up to 600 bar. A solvent selection valve allows to form binary mixtures (isocratic or gradient). Active seal wash is available for up to three modifier solvents on channel B, while channel A is always the CO₂ delivering channel.

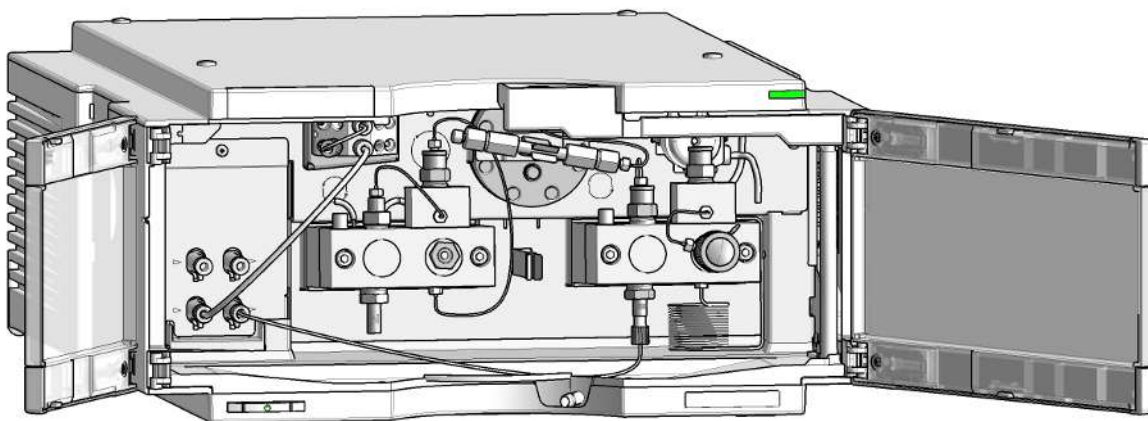


Figure 1: Overview of the pump

Product Description of the 1260 Infinity III SFC Binary Pump (G4782A)

The 1260 Infinity III SFC Binary Pump is equipped with passive inlet valves and with special seals and valves to allow for CO₂ pumping in channel A while channel B adds organic modifier for either isocratic or gradient performance. Pump head B is also equipped with a purge valve to allow for quick changeover of the organic modifier. In addition, it has an integrated solvent selection 2-channel degasser and built-in active seal-wash on channel A for increased uptime.

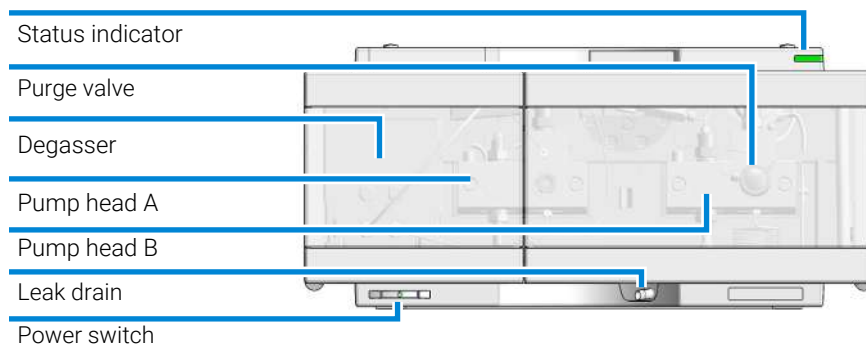


Figure 2: Overview of the pump

Features of the 1260 Infinity III SFC Binary Pump (G4782A)

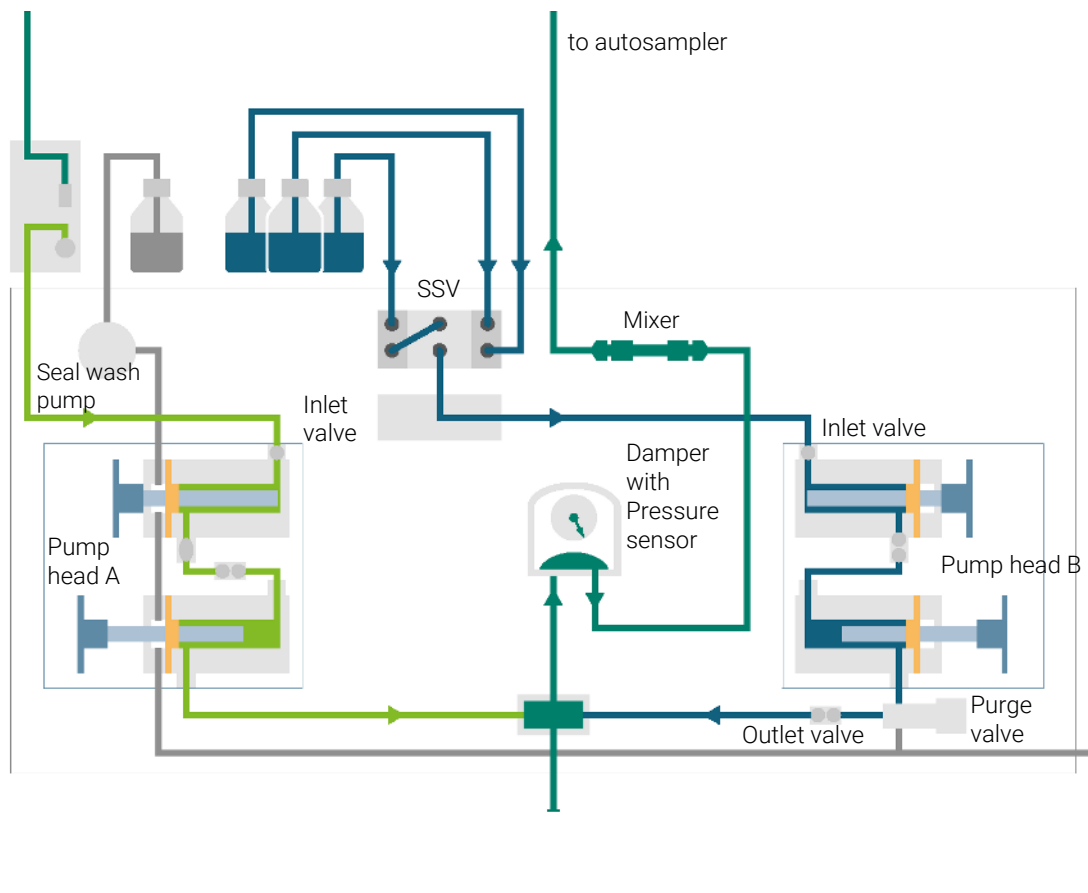
- Fast and precise gradients - the pump is the perfect choice for fast and precise gradients using UV-only, as well as SFC/MS systems.
- Power range combining high pressure up to 600 bar and flow rates up to 5 mL/min for maximum SFC performance.
- Integrated 2-channel degasser change-over of solvents for purging and priming the pump.
- Built-in active seal-wash for increased uptime.
- Integrated Solvent Selection Valve.
- InfinityLab Level Sensing adds weight controlled level sensing and solvent prediction of the equipped HPLC solvents to avoid downtime by running out solvent.

Operating Principle

Principle of Operation

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

Each channel comprises a pump assembly including pump drive, pump head, and outlet valve. The two channels are fed into cross connector which is connected to a damping unit and a mixer. A pressure sensor monitors the pump pressure. A purge valve with integrated PTFE frit is fitted to the Pump Head B outlet for convenient priming of the modifier channel.



For pump specifications, see [Specifications of the 1260 Infinity III SFC Binary Pump \(G4782A\)](#) on page 22.

Overview of the Hydraulic Path

The solvent from the bottle in the solvent cabinet enters the pump through an active inlet valve. Each side of the binary pump comprises two substantially identical pump units. Both pump units comprise a ball-screw drive and a pump head with two sapphire pistons for reciprocating movement.

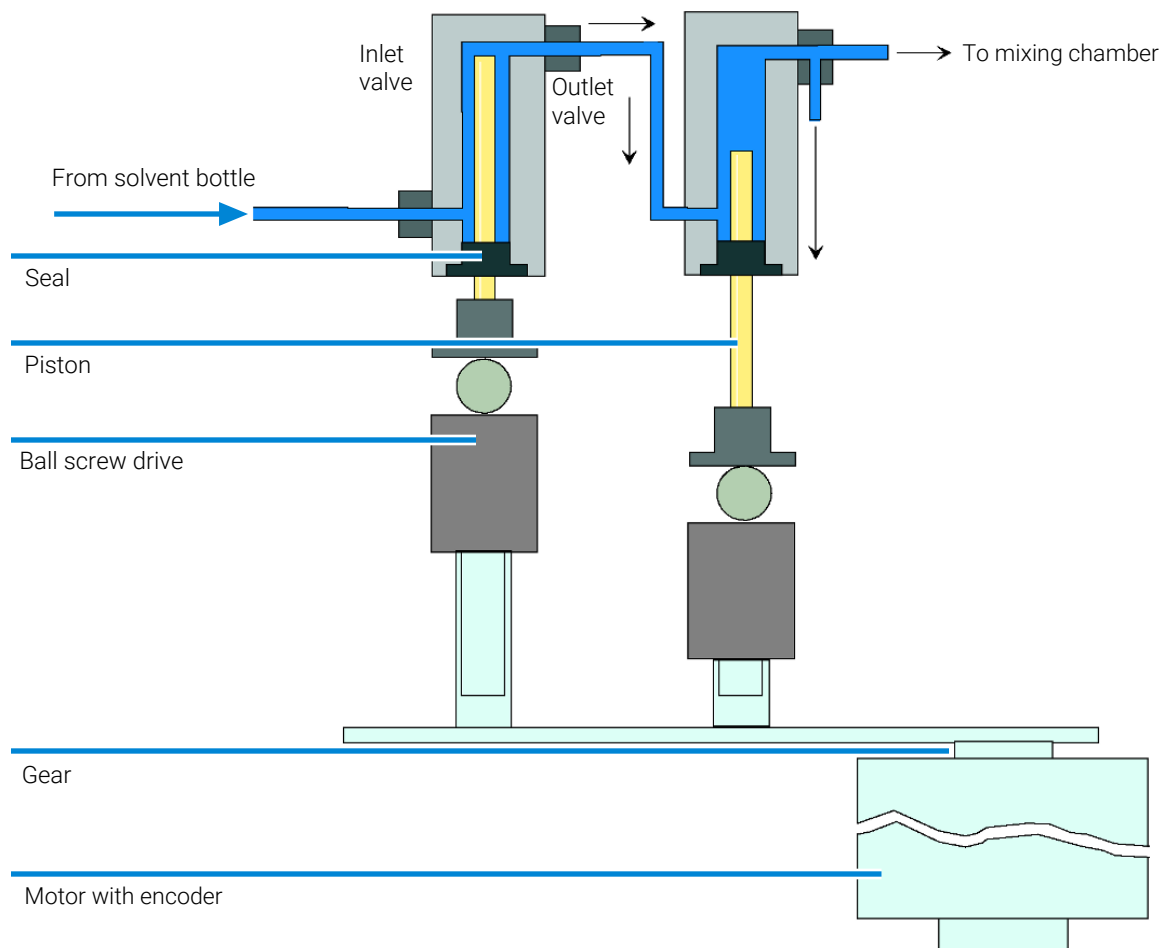


Figure 3: Pump head

A servo-controlled variable reluctance motor drives the two ball-screw drives in opposite directions. The gears for the ball-screw drives have different circumferences (ratio 2:1) allowing the first piston to move at double the speed of the second piston. The solvent enters the pump heads close to the bottom limit and leaves it at its top. The outer diameter of the piston is smaller than the inner diameter of the pump-head chamber allowing the solvent to fill the gap in between. The first piston has a stroke volume in the range of 20 μL to 100 μL depending on the flow rate. The microprocessor controls all flow rates in a range

Introduction

Operating Principle

of 1 $\mu\text{L}/\text{min}$ to 5 mL/min . The inlet of the first pumping unit is connected to the active inlet valve which is processor-controlled opened or closed allowing solvent to be drawn into the first pump unit.

When turned on, the pump runs through an initialization procedure to determine the upper dead center of the first piston of both pump channels. The first piston moves slowly upwards to the mechanical stop of the pump head and from there it moves back a predetermined path length. The controller stores this piston position in memory. After this initialization the pump starts operation with the set parameters for the two pump channels.

The inlet valve is opened and the down moving piston draws solvent into the first pump head. At the same time the second piston is moving upwards delivering into the system. After a controller defined stroke length (depending on the flow rate) the drive motors are stopped and the inlet valve is closed. The motor direction is reversed and moves the first piston up until it reaches the stored upper limit and at the same time moving the second piston downwards.

Then the sequence starts again moving the pistons up and down between the two limits. During the delivery stroke of the first piston the solvent in the pump head is pressed through the outlet valve into the second pumping unit. The second piston draws in half of the volume displaced by the first piston and the remaining half volume is directly delivered into the system. During the drawing stroke of the first piston, the second piston delivers the drawn volume into the system.

What is Solvent Compressibility Compensation?

Although the compressibility of liquids is orders of magnitude lower than the compressibility of gases, without correction a noticeable volume error would be seen if typical chromatographic solvents are compressed to operating pressures as high as 600 bar. In addition, the compressibility depends on pressure, temperature and the amount of dissolved gas. In order to minimize the influence of the latter, the use of a vacuum degasser is mandatory for a high flow and composition precision. Unfortunately, the influence of the temperature on compressibility is non-linear and cannot be calculated.

The Agilent 1260 Infinity III Binary Pump features a multi point compressibility calibration. The compressibility of a solvent is determined at different pressures from 0 – 600 bar and stored in an XML file. This file can be distributed to other pumps because the solvent compressibility is independent from the pump.

Introduction

Operating Principle

The binary pump and ChemStation come with predetermined solvent compressibility data for the most common HPLC solvents like water, acetonitrile, methanol, etc. Users can calibrate their own solvent mixtures with the help of an easy to use calibration procedure in the Agilent Lab Advisor software.

Let us use the practical example from the last section once again to understand how compressibility compensation works:

Piston 1 draws solvent at ambient pressure. The movement direction is reversed and piston 1 now compresses the solvent until the operating pressure of the HPLC system is reached. The outlet valve opens, and solvent is pumped by piston 1 into pump chamber 2.

Without any compensation, the delivered volume at operating pressure would be too low. In addition, it would take a noticeable amount of time to recompress the solvent to operating pressure. During this time frame, no solvent would be delivered into the system and as a result a high pressure fluctuation (known as *pressure ripple*) would be observed.

When both solvent compressibility at the current operating pressure and pump elasticity are known, the pump can automatically correct for the missing volume by drawing the appropriate larger solvent volume at ambient pressure and speed up the piston during the recompression phase in the first pump chamber. As a result, the pump delivers the accurate volume with any (calibrated) solvent at any pressure at a greatly reduced pressure ripple. For applications that require lowest transition volume of the pump, damper and mixer can be bypassed.

How Does Variable Stroke Volume Work?

The smaller the solvent volume in the pump chamber is, the faster it can be recompressed to operating pressure. The binary pump allows to manually or automatically adjust the pump stroke volume of the first piston in the range of 20 – 100 μL . Due to the compression of the solvent volume in the first pump chamber, each piston stroke of the pump will generate a small pressure pulsation, influencing the flow ripple of the pump. The amplitude of the pressure pulsation mainly depends on the stroke volume and the compressibility compensation for the solvent in use. Small stroke volumes generate less pressure pulsation than larger stroke volumes at the same flow rate. In addition, the frequency of the pressure pulsation will be higher. This will decrease the influence of flow pulsations on retention times.

In gradient mode, a smaller stroke volume results in less flow ripple and reduces the composition ripple.

Introduction

Operating Principle

The binary pump uses a processor-controlled ball screw system for driving its pistons. The normal stroke volume is optimized for the selected flow rate. Small flow rates use a small stroke volume while higher flow rates use a higher stroke volume.

The stroke volume for the pump is by default set to AUTO mode. This means that the stroke is optimized for the flow rate in use. A change to larger stroke volumes is possible but not recommended.



2

Site Requirements and Specifications

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

Site Requirements 19

Power Considerations 19

Power Cords 20

Bench Space 21

Condensation 21

Specifications of the 1260 Infinity III SFC Binary Pump (G4782A) 22

Site Requirements

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

Power Considerations

The module power supply has wide ranging capability. It accepts any line voltage in the range described in [Table 1](#) on page 22. Consequently there is no voltage selector in the rear of the module. There are also no externally accessible fuses, because automatic electronic fuses are implemented in the power supply.

WARNING

Incorrect line voltage at the module

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

- Connect your module to the specified line voltage.

WARNING

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

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

- Make sure that it is always possible to access the power plug.
- Remove the power cable from the instrument before opening the cover.
- Do not connect the power cable to the Instrument while the covers are removed.

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.
-

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 module dimensions and weight (see [Table 1](#) on page 22) allow you to place the module on almost any desk or laboratory bench. It needs an additional 2.5 cm (1.0 inches) of space on either side and approximately 8 cm (3.1 inches) in the rear for air circulation and electric connections.

If the bench shall carry a complete HPLC system, make sure that the bench is designed to bear the weight of all modules.

The module should be operated in a horizontal position.

NOTE

Agilent recommends that you install the HPLC instrument in the InfinityLab Flex Bench rack. This option helps to save bench space as all modules can be placed into one single stack. It also allows to easily relocate the instrument to another lab.

Condensation

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 SFC Binary Pump (G4782A)

Table 1: Physical Specifications of the 1260 Infinity III SFC Binary Pump (G4782A)

Type	Specification	Comments
Weight	17.8 kg (39.2 lbs)	
Dimensions (height x width x depth)	180 x 396 x 436 mm (7.1 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	90 VA / 74 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	Installation category II, Pollution degree 2	For indoor use only
ISM Classification	ISM Group 1 Class B	According to CISPR 11

Site Requirements and Specifications

Specifications of the 1260 Infinity III SFC Binary Pump (G4782A)

Table 2: Performance Specifications of the 1260 Infinity III SFC Binary Pump (G4782A)

Type	Specification	Comments
Hydraulic system	Two dual piston in series pumps with servo-controlled variable stroke drive, power transmission by gears and ball screws, floating pistons	
Designed for use with Agilent InfinityLab Assist Technology	Intuitive User Interface, Automated Workflows, Predictive Maintenance & Assisted Troubleshooting	
Flow range	Settable: 0.001 – 5 mL/min	Recommended: 1.0 – 5.0 mL/min (in 0.001 mL/min increments)
Flow precision	$\leq 0.07\%$ RSD or < 0.02 min SD, whichever is greater	Based on retention time at constant temperature
Pressure operating range	Up to 60 MPa (600 bar, 8702 psi) up to 5 mL/min	
Compressibility compensation	Pre-defined, based on mobile phase compressibility	
Recommended pH range	1.0 – 12.5	Solvents with pH < 2.3 should not contain acids that attack stainless steel
Gradient formation	High-pressure binary mixing	
Delay volume	<i>Standard delay volume configuration:</i> 600 – 900 μ L, (includes 400 μ L mixer), dependent on back pressure	
Composition range	settable: 0 – 100 %	
Composition precision	$< 0.15\%$ RSD or < 0.04 min SD, whichever is greater	Based on retention time at constant temperature
Integrated degassing unit	Number of channels: 2 Internal volume per channel: 1.5 mL	
Instrument Control	LC & CE Drivers A.02.16 or above Instrument Control Framework (ICF) A.02.04 or above Instant Pilot (G4208A) with firmware B.02.21 or above InfinityLab Assist (G7180A) with firmware D.07.40 or above Lab Advisor software B.02.09 or above	For details about supported software versions refer to the compatibility matrix of your version of the LC and CE Drivers
Communication	Controller-area network (CAN), Extended Remote Interface (ERI), Local Area Network (LAN)	

Site Requirements and Specifications

Specifications of the 1260 Infinity III SFC Binary Pump (G4782A)

Type	Specification	Comments
Safety features and maintenance	Leak detection, safe leak handling, leak output signal for shutdown of the pumping system. Low voltage in major maintenance areas. Extensive diagnostics, error detection and display with Agilent Lab Advisor software.	
GLP features	Early maintenance feedback (EMF) for continuous tracking of instrument usage in terms of seal wear and volume of pumped mobile phase with pre-defined and user settable limits and feedback messages. Electronic records of maintenance and errors.	
Housing	All materials are recyclable	



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 26

Install Capillaries 26

Handling Leak and Waste 30

Drain Connectors Installation 33

SFC Waste Concept 37

Waste Guidance 39

Leak Sensor 39

Connecting Modules and Control Software 40

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 234.

Table 3: 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

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

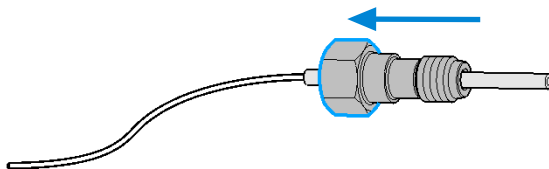
Installation

Installing Capillaries

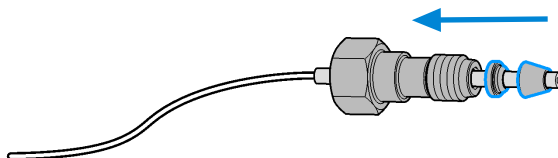
- 1 Select a nut that is long enough for the fitting you'll be using.



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



- 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.



Installation

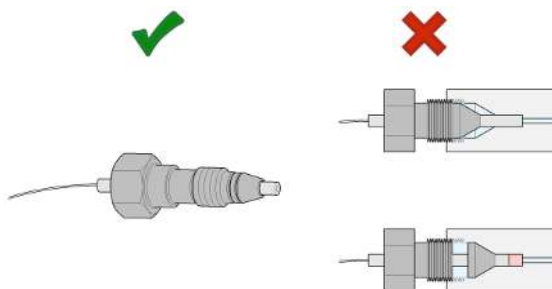
Installing Capillaries

- 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

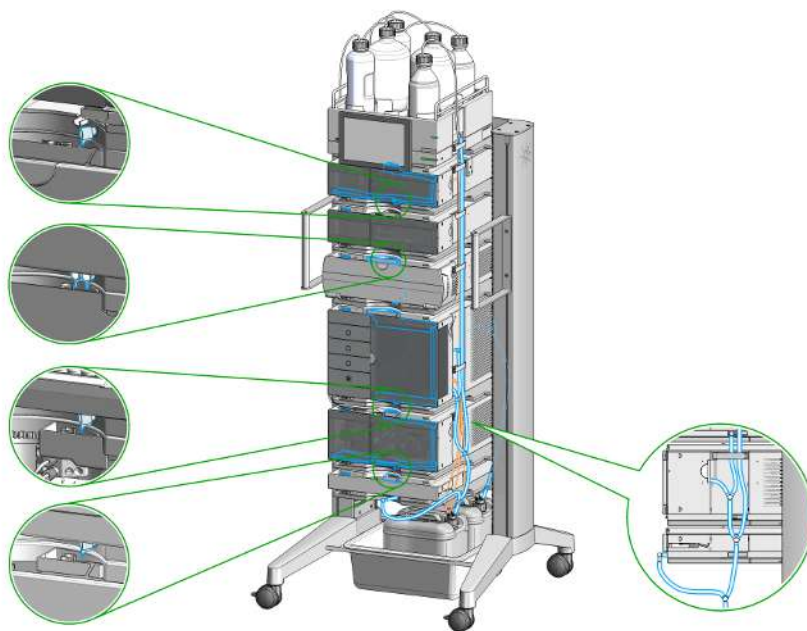


Figure 4: Infinity III Leak Waste Concept (Flex Bench installation)

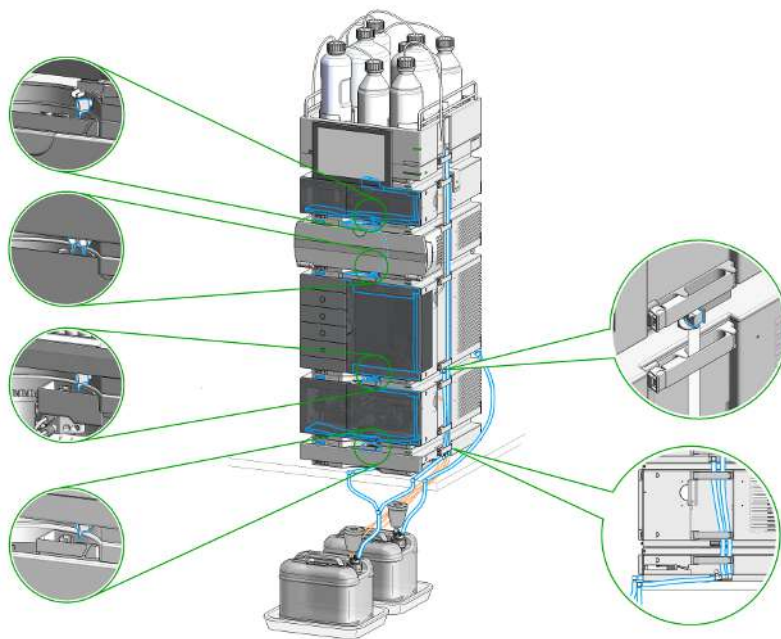


Figure 5: Infinity III Single Stack Leak Waste Concept (bench installation)

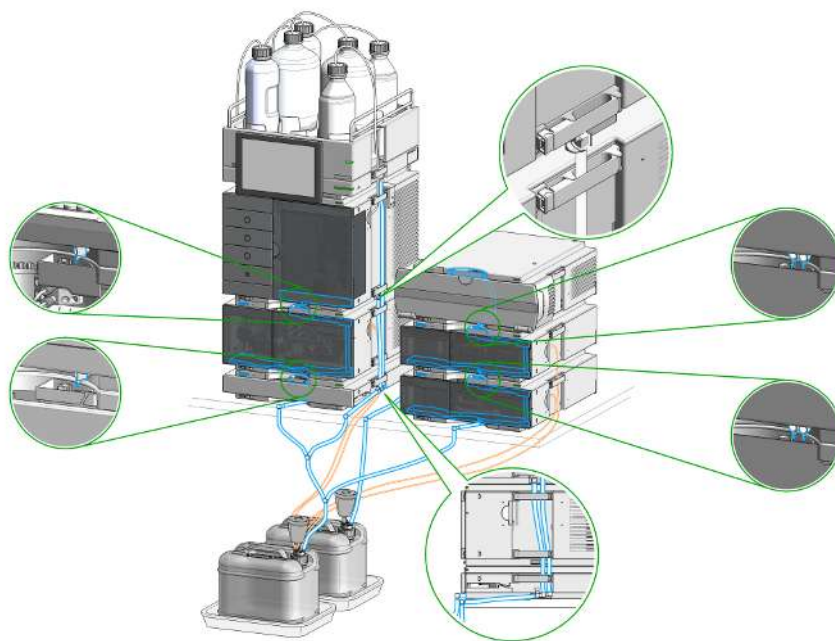


Figure 6: Infinity III Two Stack Leak Waste Concept (bench installation)

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 30.

Parts required

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

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

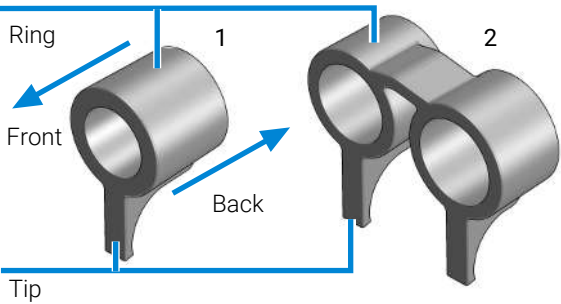


Figure 7: Overview of Drain Connectors: Single (left) and Double (right)



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 4: 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

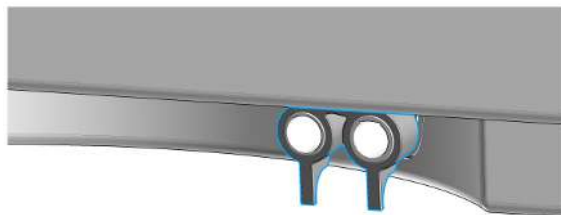
In case of incorrect installation, drain connectors cannot fully perform the intended function.

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)**

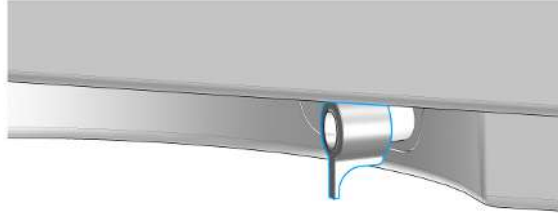
- 1 Align the rings with the leak drain outlets of the module, press slightly with the fingers, and slide the connector along the leak drain outlets until it is aligned with the front of the leak drain.

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

Installation

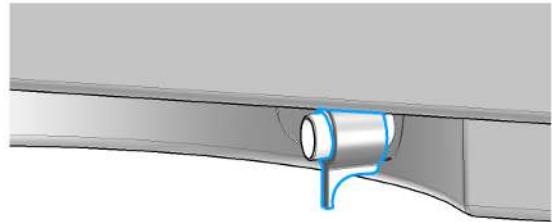
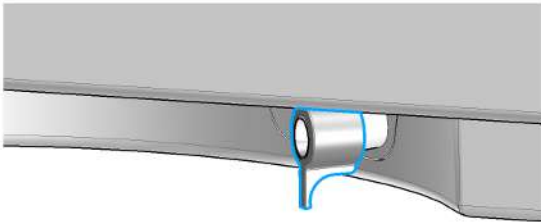
Handling Leak and Waste

- 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.



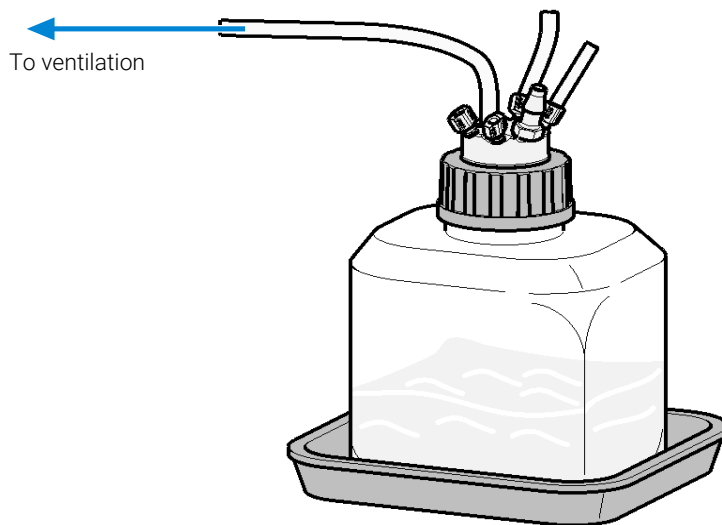
Make sure that the following requirements are covered:

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

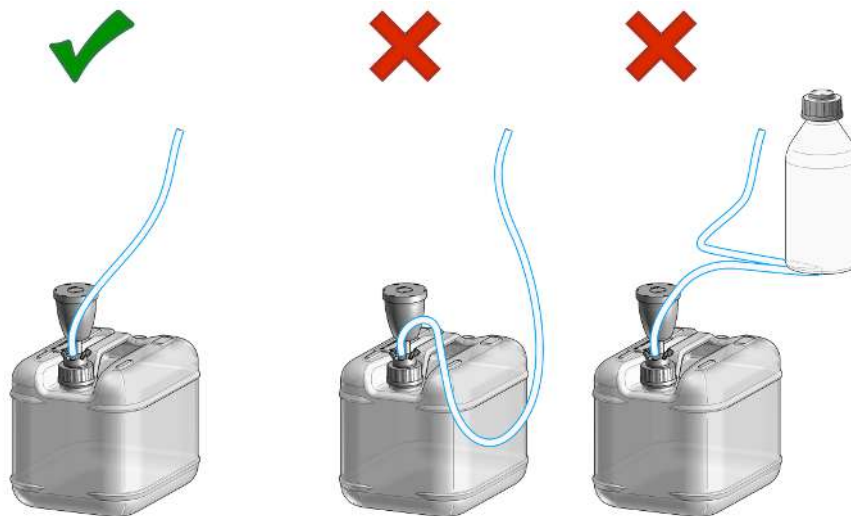


SFC Waste Concept

- 1 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.
-



4 Using the Module

This chapter provides information on how to use the module.

General Information 42

Turn on/off 42

Status Indicators 44

Best Practices 45

Daily / Weekly Tasks 45

Power-Up/Shut-Down the Pump 45

Prepare the Pump 46

How to Deal With Solvents 46

Hints for Successful Use of the Binary Pump 47

Prevent Blocking of Solvent Filters 47

Normal Phase Applications 49

Preparation of the System 51

Prepare a Run 51

Prime and Purge the System 58

Preparation of the Module 60

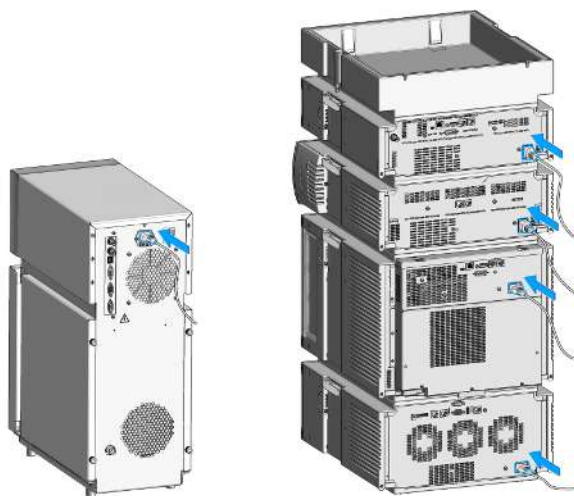
Setting up the Pump with the G4208A Instant Pilot 60

Setting up the Pump with the Instrument Control Interface 60

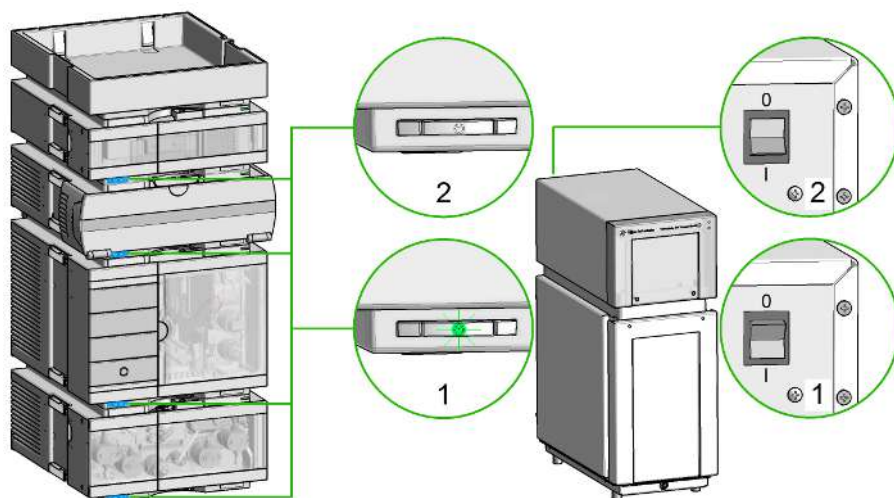
General Information

Turn on/off

- 1 Plug in power cord



2

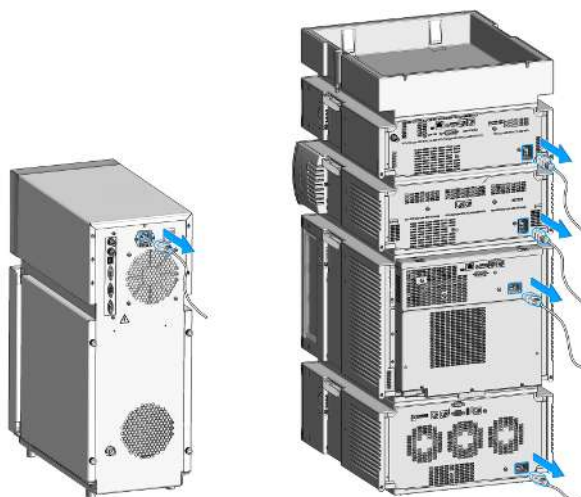


Power switch

(1) On

(2) Off

3 Plug off the module.



Status Indicators

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

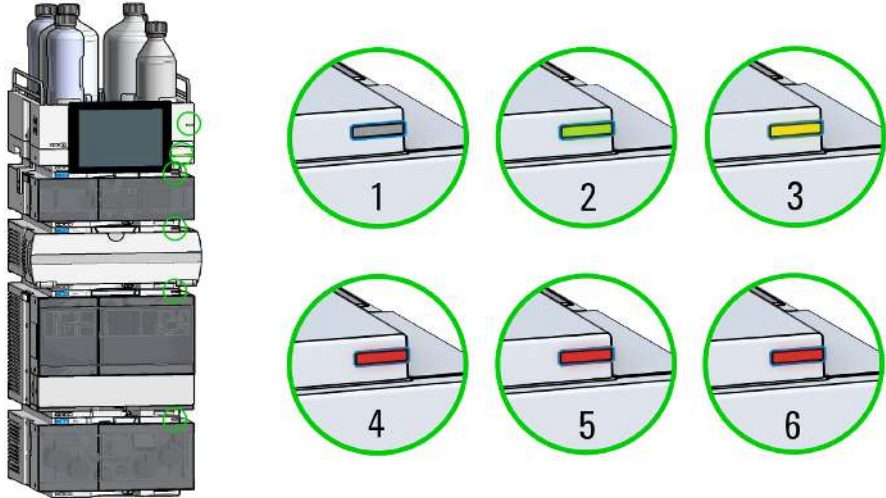


Figure 8: Arbitrary LC stack configuration (example)

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**.

Best Practices

Daily / Weekly Tasks

Daily Tasks

- Replace mobile phase based on water/buffer.
- Replace organic mobile phase latest every second day.
- Check seal wash solvent.

Weekly Tasks

- Change seal wash solvent (10 % isopropanol in water) and bottle.
- If applications with salts were used, flush all channels with water and remove possible salt deposits manually.
- Inspect solvent filters for dirt or blockages. Exchange if no flow is coming out of the solvent line when removed from the degasser inlet.

Power-Up/Shut-Down the Pump

Power Up the Pump

- Use new or different mobile phase (as required).
- Purge pump heads with 2.5 – 3 mL/min for 5 min.
- Stabilize the system by running for 10 – 20 min.

Long-Term Shut-Down of the System

- Flush system with water to remove buffer.
- Remove all samples from the sampler and store according to good laboratory practice.
- Use recommended solvents to store the system.

- Power off the system.

Prepare the Pump

Purge

Use the Purge function to:

- fill the pump,
- exchange a solvent,
- remove air bubbles in tubes and pump heads.

Seal Wash

Seal Wash guarantees a maximum seal life time. Use Seal Wash:

- When using buffers with elevated salt concentrations
- When using volatile solvents with non-volatile additives

CAUTION

Contaminated seal wash solvent

- Do not recycle seal wash solvent to avoid contamination.
- Weekly exchange seal wash solvent.

How to Deal With Solvents

- Use clean bottles only.
- Exchange water-based solvents daily.
- Select solvent volume to be used up within 1 – 2 days.
- Use only HPLC-grade solvents and water filtered through 0.2 µm filters.
- Label bottles correctly with bottle content, and filling date / expiry date.
- Use solvent inlet filters.
- Reduce risk of algae growth: use brown bottles for aqueous solvents, avoid direct sunlight.

Hints for Successful Use of the Binary Pump

- Place solvent cabinet with the solvent bottles always on top (or at a higher level) of the pump.
- Flush the degasser with at least 5 mL per channel before operating the pump, especially when the pumping system had been turned off for a certain length of time (for example, overnight) and volatile solvent mixtures are used in the channels.
- Prevent blocking of solvent inlet filters (never use the pump without solvent inlet filters). Growth of algae should be avoided (see [Prevent Blocking of Solvent Filters](#) on page 47).
- Check purge valve frit and column frit in regular time intervals. A blocked purge valve frit can be identified by black, yellow or greenish layers on its surface or by a pressure greater than 10 bar in low delay volume configuration and 20 bar in standard configuration when pumping distilled water at a rate of 5 mL/min with an open purge valve.
- Whenever possible use a minimum flow rate of 5 μ L/min per solvent channel to avoid crossflow of solvent into the unused pump channel.
- Whenever exchanging the pump seals, the purge valve frit should be exchanged, too.
- Check the pump pistons for scratches, grooves and dents when changing the piston seals. Damaged pistons cause micro leaks and will decrease the lifetime of the seals.
- After changing the piston seals, apply the seal wear-in procedure (see [Seal Wear-in Procedure](#) on page 142).

Prevent Blocking of Solvent Filters

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

- Use a sterile, if possible amber, solvent bottle to slow down algae growth.
- Filter solvents through filters or membranes that remove algae.
- Exchange solvents every two days or refilter.

- If the application permits add 0.0001 – 0.001 M sodium azide to the solvent.
- Place a layer of argon on top of your solvent.
- Avoid exposure of the solvent bottle to direct sunlight.

NOTE

Never use the system without solvent filter installed.

Checking the Solvent Filters

The solvent filters are located on the low-pressure side of the binary pump. A blocked filter therefore does not necessarily affect the high pressure readings of the pump. The pressure readings cannot be used to check whether the filters are blocked or not. If the solvent cabinet is placed on top of the binary pump, the filter condition can be checked in the following way:

Remove the solvent inlet tube from the inlet port of the solvent selection valve or the degasser. If the filter is in good condition, the solvent will freely drip out of the solvent tube (due to hydrostatic pressure). If the solvent filter is partly blocked only very little solvent will drip out of the solvent tube.

WARNING

When opening capillary or tube fittings, solvents may leak out.

The handling of toxic and hazardous solvents and reagents can carry health risks.

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

Cleaning the Solvent Filters

- Remove the blocked solvent filter from the bottle-head assembly and place it in a beaker with concentrated nitric acid (35%) for one hour.
- Thoroughly flush the filter with HPLC-grade water (remove all nitric acid, some capillary columns can be damaged by nitric acid).
- Replace the filter.

NOTE

Never use the system without solvent filter installed.

Normal Phase Applications

Current passive inlet valves and outlet ball valves used with 1260 and 1290 Infinity pumps do not work well with applications using non-polar solvents as for normal phase applications (e.g. hexane, heptane and CO₂). With such applications, pressure drops could be observed. They are a result of particles electrostatically charging up in insulating solvents and sticking to the balls inside the valves, such that the valves do not close properly any more after some time of use (can be hours).

For normal phase applications, a second type of valves is available, which has a design based on the existing one for 1260 and 1290 Infinity valves. These valves use a new material for valve balls, which is a conductive ceramic and replaces non-conductive ruby balls. The balls do not charge up electrostatically and show good performance in normal phase.

The valves are marked with N for non-polar or normal phase.

Agilent recommends using these valves for (and only for) normal phase applications.

CAUTION

Corrosion of valves

Normal phase balls/valves corrode quickly in aqueous solutions and acids (at or below pH 7).

- Do not use normal phase valves in applications running with aqueous solutions.

The N-Valves have been tested successfully in using hexane at pressures below 100 bar; heptane can be used as a substitute for neurotoxic hexane.

Seals for Normal Phase Applications

For running normal phase applications on 1200 Infinity Series pumps, yellow PE seals are required, which exist as piston seals and wash seals. Seal wash is very uncommon for normal phase applications (no buffers needed), but wash seals are needed for seal wash pump heads.

1290 Infinity pumps use PE seals by default. In combination with ceramic pistons, PE seals are used for both reversed phase (1200 bar) and normal phase applications.

1260 Infinity pumps use sapphire pistons and black PTFE piston and wash seals by default (600 bar). Such PTFE seals create small wear particles in normal phase applications, which can clog valves and other parts in the flow path.

PE seals have a limited life time when used with normal phase solvents and sapphire pistons. Agilent recommends a maximum pressure of 200 bar for this combination, which shall also be applied for pressure tests.

Choice of Normal Phase Valves and Seals

Table 5: Recommended valves and seals for normal phase applications

	1260 Infinity	1290 Infinity
Inlet valves	G1312-60166 (1260 Inlet Valve Type N)	G4220-60122 (1290 Inlet Valve Type N) G4204-60122 (1290 Quat Inlet Valve Type N)
Outlet valves	G1312-60167 (Outlet Valve Type N/ SFC)	G4220-60128 (1290 Outlet Valve Type N)
Seals	0905-1420 (PE seal (pack of 2)) 0905-1718 (Wash Seal PE)	

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

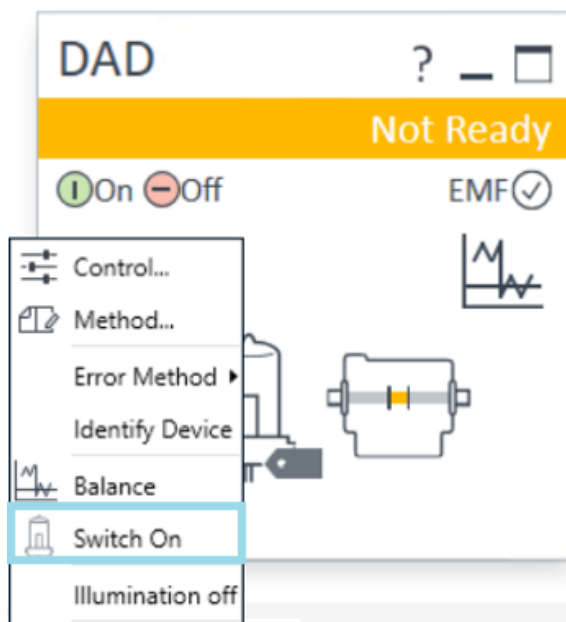
The handling of solvents, samples and reagents can hold health and safety risks.

- 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

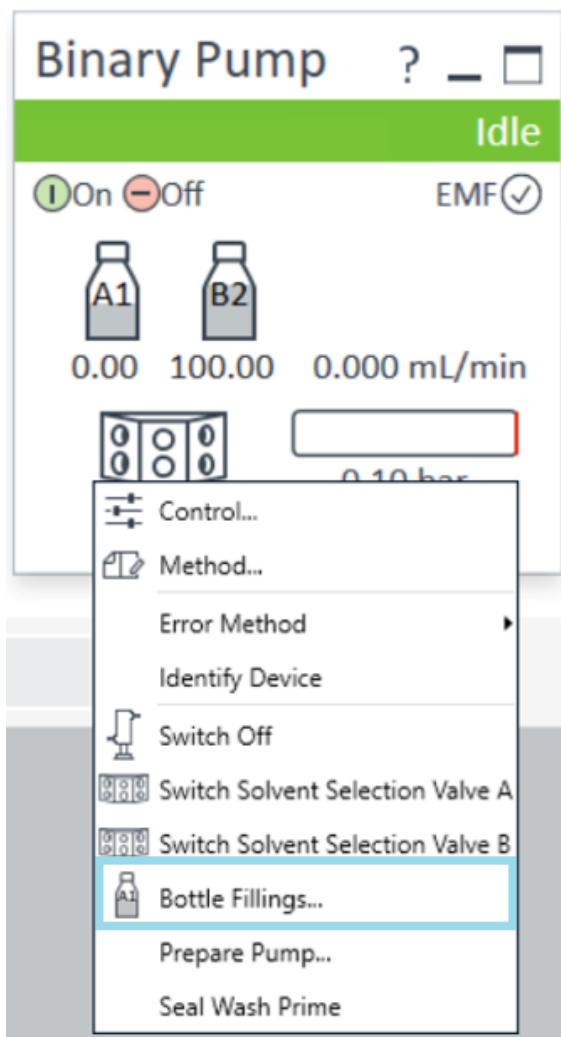
Preparation of the System

- 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 0.00 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 0.00 liter

☐ Turn pump off if waste volume has reached maximum limit

Ok Cancel Help

6 Purge the pump.

NOTE

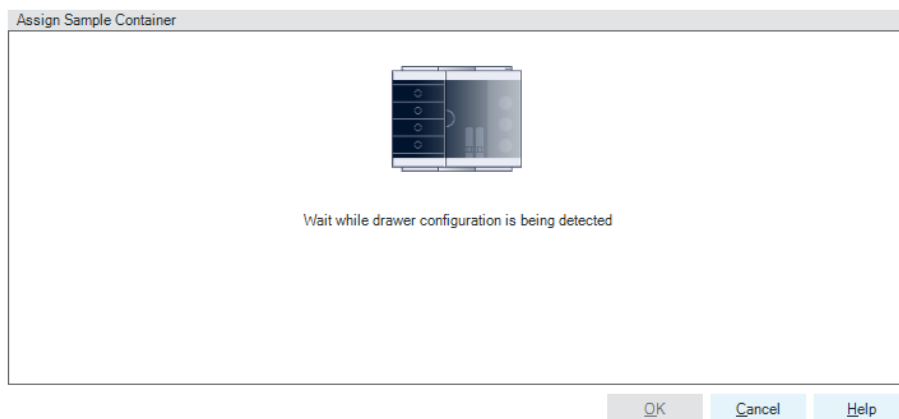
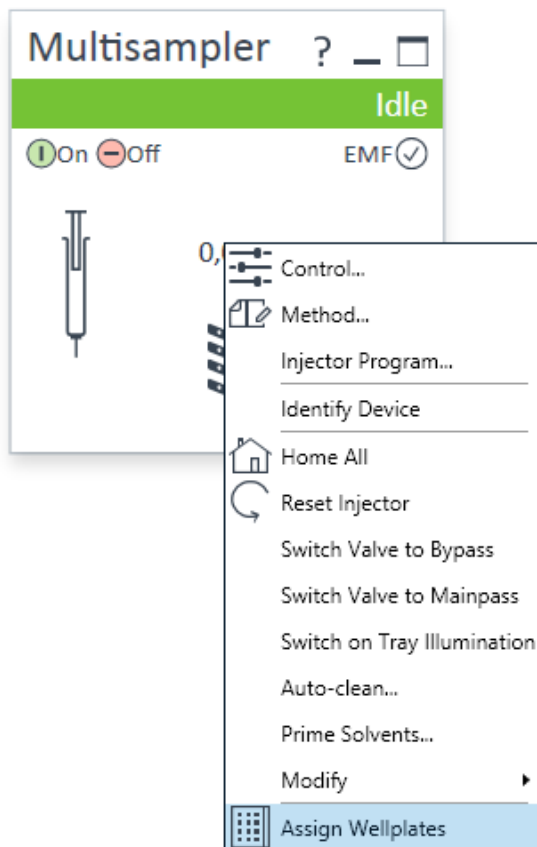
For details on priming and purging, refer to the technical note *Best Practices for Using an Agilent LC System Technical Note (InfinityLab-BestPractice-en-SD-29000194.pdf, SD-29000194)*.

7 Change solvent type if necessary.

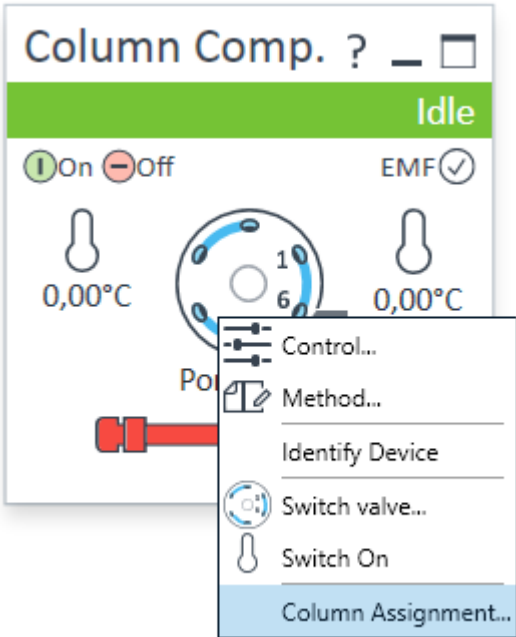
Using the Module

Preparation of the System

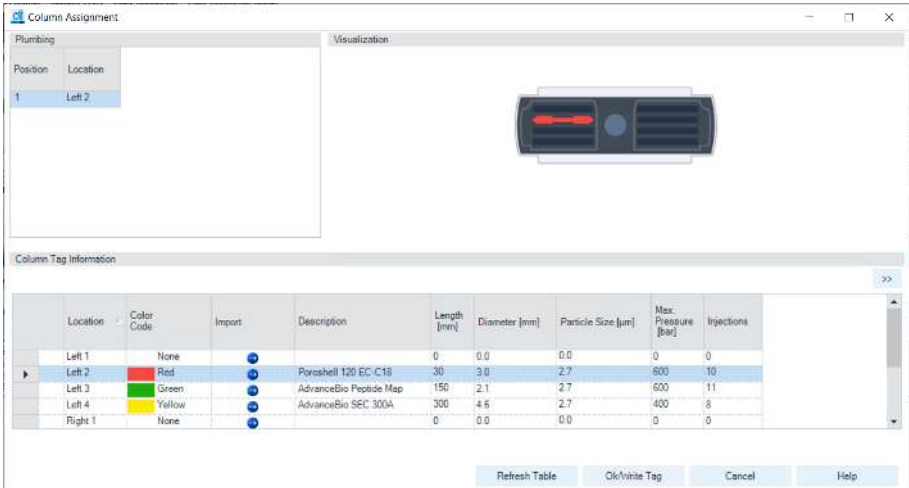
- 8 Choose the tray format of the sampler.



9 Add a new column.



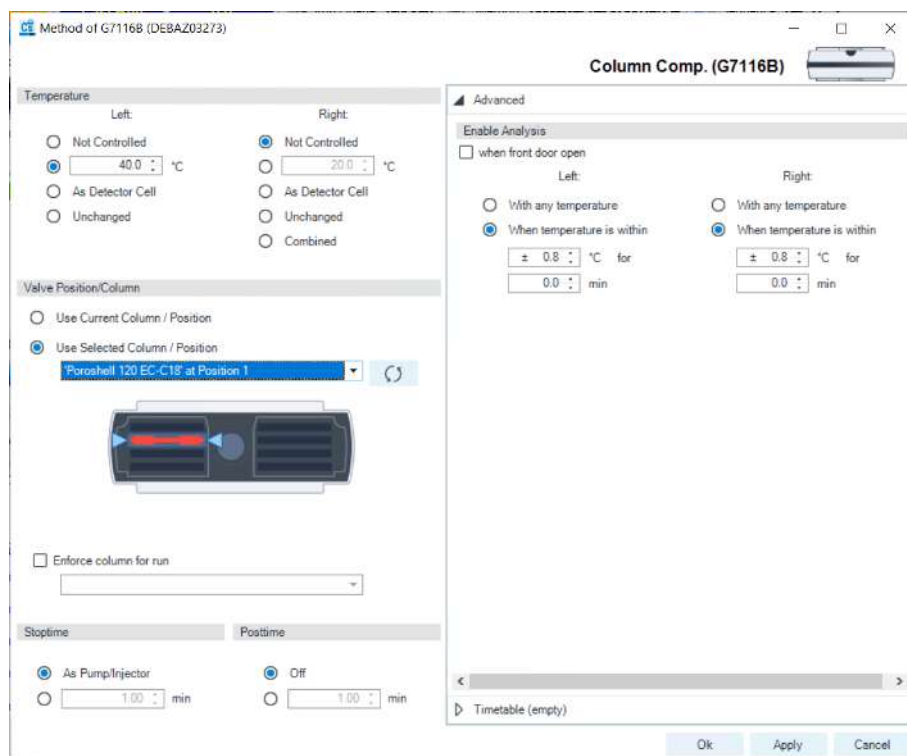
10 Enter the column information.



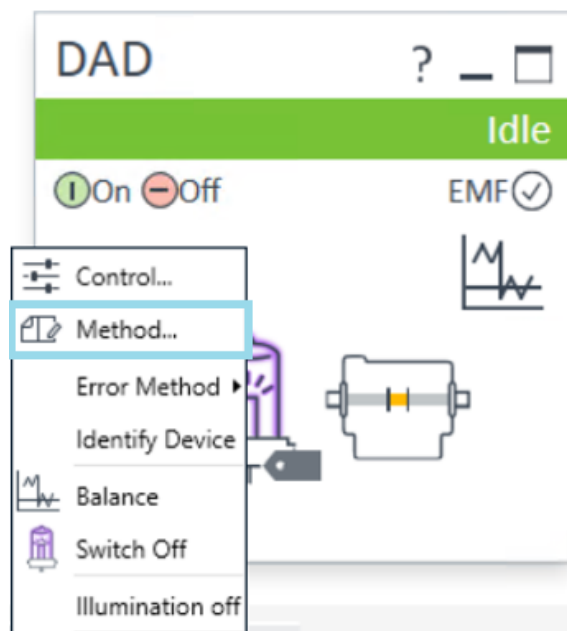
Using the Module

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 6: 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.

Preparation of the Module

Setting up the Pump with the G4208A Instant Pilot

Generic operation of the G4208A Instant Pilot is covered in the G4208-90006 (Agilent Instant Pilot G4208A User's Guide) . Details about setting up module specific parameters can be found in the Instant Pilot online help.

The pump parameters are described in detail in [Overview](#) on page 60.

Setting up the Pump with the Instrument Control Interface

Overview

The instrument control interface offers the parameters described in the following sections, and can usually be accessed through Agilent instrument control software. For details, please see manuals and online help of respective user interfaces.

Setup of Basic Pump Parameters

The most important parameters of the pump are listed in the following table.

Table 7: Basic pump parameters

Parameter	Limits	Description
• Flow	0.001 – 5 mL/min	Total flow rate of the pump.
• Stop Time	0.01 min - no limit	The stop time of the pump usually controls the run time of the whole LC system. Use no limit to stop the run manually (useful for method development).
• Post Time	off - 99999 min	Time between the end of a run and the start of the next. Used for column equilibration after a gradient.

Parameter	Limits	Description
• Pressure Limits	Max: 0 – 600 bar Min: 0 – 600 bar	Max must be bigger than Min ! Set max pressure to the maximum operating pressure of your column. A min pressure setting of e.g. 10 bar will turn off your pump automatically when running out of solvent. A smarter way, however, is to use the bottle fillings function (see Bottle Filling on page 63).
• Solvent A	0 – 100 %	Although channel A can be set to 0 %, it cannot be turned off. This channel should be used for the aqueous phase (water).
• Solvent B	off - 100 %	The percentage of channel B is automatically complemented by channel A to give 100 %.
• Solvent type	H ₂ O, ACN, MeOH, IPA CO ₂ pre-compressed	Select the solvent you are using in the respective solvent channel from the drop-down list. In case your solvent is not listed, perform a solvent compressibility calibration.
• Solvent Comment		Free text field for a description of the solvent. This description will show up in method printouts, etc.
• Timetable	max. number of lines depends on free space in pump memory	Use the timetable to build solvent gradients, flow gradients, or combinations of both. Gradients are always linear. Use multiple timetable entries to mimic exponential or parabolic gradients.
• Display		There are three ways to display the timetable: <ul style="list-style-type: none"> • in tabular form • as flow/pressure graph • as solvent percentage plot Values can only be changed in tabular view.

Pump Control

The pump can be switched between following states: **On**, **Off** or to **Standby**. In **Standby**, the pump motor is still controlled. When the pump is switched on from standby, it does not re-initialize.

CAUTION

Upon initialization, the pump ignores the Maximum Flow Gradient value. This can result in a rapid and uncontrolled pressure increase.

- To prevent harm to the column, open the purge valve until the initialization is finished.

The optional seal wash pump can be controlled by either switching it off, using it for a single time or specifying frequency and duration of periodic wash intervals.

Auxiliary Pump Parameters

The auxiliary pump parameters are pre-set to fit most applications. Adjustments should only be made when required. Auxiliary pump parameters shows the available auxiliary parameters with their default values.

CAUTION

Upon initialization, the pump ignores the Maximum Flow Gradient value.

This can result in a rapid and uncontrolled pressure increase.

- To prevent harm to the column, open the purge valve until the initialization is finished.

Table 8: Auxiliary pump parameters

Parameter	Limits	Description
• Maximum Flow Gradient	0.1 – 100 mL/min ² default: 100 mL/min ²	With this parameter flow rate changes can be ramped up and down slowly to avoid pressure shocks to the column. The default value is 100 mL/min ² which in fact turns the function off.
• Minimum Stroke	20 – 100 µL default: Auto	The volume one pump piston delivers per stroke. In general, a smaller stroke volume results in lower pump ripple. The Auto setting adjusts the strokes dynamically to the lowest possible value. The strokes can be set individually for pump heads A and B.
• Compressibility	0 - 150·10 ⁻⁶ /bar or enhanced compressibility calibration default: use enhanced comp. calibration	For best performance, check option Use enhanced compressibility calibration . With this option, the pump will use solvent data libraries provided by Agilent or data generated by using solvent compressibility calibrations. For backward compatibility to 400 bar pumps, the solvent compressibility can still be set manually for each channel when the box is unticked.

Data Curves

The binary pump provides the possibility to store the following operational data in the data file of the Agilent data system:

- Solvent percentage for each channel,
- pump flow,
- pressure

NOTE

The pressure data curve is *generated* from the pressure sensor readings, while %A, %B and flow are *calculated* from the method settings of the pump.

For details, please refer to the online help or manual of your instrument control software.

Bottle Filling

The pump offers a powerful feature to monitor the liquid level in the solvent bottles. With total bottle volume and initial filling volume set correctly, the pump subtracts the consumed volume continuously from the initial value and stops the pump and method/sequence execution before the system runs dry or an analysis is corrupted.

CAUTION

The bottle filling feature fails if multiple solvent inlets are put into one solvent bottle!

- In that case implement a minimum pressure limit to avoid that the pump runs dry when solvents are empty.

[Table 9](#) on page 63 lists the available bottle filling parameters.

Table 9: Bottle Filling Parameters

Parameter	Limits	Description
• Total Volume	0 – 1000 L default: 0 L	This is the capacity (maximum possible volume) in liter of the solvent bottle. In combination with the actual volume, this parameter is used for calculating and displaying the relative liquid level.
• Actual Volume	0 – 1000 L default: 0 L	After filling the solvent bottles, enter the actual volumes into these boxes. The Actual Volume must not be larger than the Total Volume of the bottle.
• Prevent analysis...default: unchecked		If this option is checked, the pump won't start a new run if the solvent level in one or more bottles is below the minimum volume. Enter a minimum volume in liter, which considers the position of the solvent inlet and size/shape of the solvent bottle such that no air is drawn if the actual volume gets close to this limit.
• Turn pump off... default: unchecked		If this option is checked, the pump will turn off before air is aspirated. However, the residual solvent volume has been calculated for 1 L solvent bottles and may be too small for large bottles or other vessels.

5

Optimizing the Performance of the Module

This chapter provides information on how to optimize the module.

When to Use a Vacuum Degasser 65

Operational Hints for the Vacuum Degasser 65

When to Use the Active Seal Wash Option 66

How to Optimize the Compressibility Compensation Setting 67

Optimization of Legacy Compressibility Settings 67

When to Use a Vacuum Degasser

A degasser removes air, which is dissolved in any solvent. When solvents are heated or mixed with other solvents, air can leave the solvent and form small bubbles. Over time, these bubbles accumulate and can cause pressure fluctuations which may finally result in retention time shifts.

All Agilent 1200 Infinity II/III Series Pumps have a built-in degasser. While a degasser is needed for low pressure mixing pumps like Agilent quaternary pumps, high pressure mixing pumps like Agilent binary pumps are more robust with respect to bubble formation. However, a degasser is recommended for best performance.

Additionally, a degasser is highly recommended for the following applications:

- Your detector is used with maximum sensitivity in the low UV wavelength range.
- Your application requires highest injection precision.
- Your application requires highest retention-time reproducibility (flow rates below 0.5 mL/min).
- The binary pump is used with bypassed damper and mixer.

The external G7122A Degasser is recommended for use with applications using highly volatile solvents like Hexane or DCM, solvents with special characteristics like THF, or applications using refractive index detection.

Operational Hints for the Vacuum Degasser

If you are using the vacuum degasser for the first time, if the vacuum degasser was switched off for any length of time (for example, overnight), or if the vacuum degasser chambers are empty, you have to prime the vacuum degasser before running an analysis. Priming is usually done by pumping at a high flow rate (3 – 5 mL/min). Alternatively, a syringe can be used to draw the solvent through the (empty) degasser if the pump does not aspirate the solvent by itself.

When to Use the Active Seal Wash Option

Concentrated buffer solutions will reduce the lifetime of the seals and pistons in your binary pump. The active seal wash option allows to maintain the seal lifetime by flushing the low pressure side of the seals with a wash solvent.

The seal wash option is strongly recommended if buffer concentrations of 0.1 M or higher are used regularly with the pump.

The seal wash option comprises a peristaltic pump, secondary seals, gaskets, seal holders and tubing for both pump heads. A bottle of premixed water/isopropanol (90 /10 vol%) is placed in the solvent cabinet and connected to the peristaltic pump.

Always use a mixture of HPLC-grade water (90 %) and isopropanol (10 %) as wash solvent. This mixture prevents bacteria growth in the wash bottle and reduces the surface tension of the water.

NOTE

In order to avoid accumulation of buffer salts or impurities, regularly replace the washing solution using fresh solvents.

The operation of the peristaltic pump can be controlled from the data system or the Instant Pilot.

For adding a seal-wash option, please contact your local Agilent Technologies service representative.

How to Optimize the Compressibility Compensation Setting

When a solvent is metered at ambient pressure and compressed to a higher pressure, the volume decreases depending on its compressibility. Solvent compressibility is a non-linear function of pressure and temperature. It is specific for each solvent.

In order to deliver the desired flow accurately at all pressures, Agilent pumps use a compressibility compensation. For standard LC applications, e.g. using a 400 bar binary pump, an average compressibility value for the solvent is sufficient.

For the 600 bar 1260 Infinity III Binary Pump, the pressure-dependency of a solvent compressibility needs to be considered. It is determined at different pressures between 0 – 600 bar. The pump uses the obtained non-linear function to select the correct compressibility value for the actual pump pressure. Compressibility data for the most common solvents is readily available in the pump firmware.

The compensation algorithm is so powerful that the damper and mixer can be removed from the pump flow path at low flow rate while the pressure ripple and composition ripple remain at low levels.

For method compatibility reasons, the legacy compressibility compensation is still available.

Optimization of Legacy Compressibility Settings

The compressibility compensation default settings are 50×10^{-6} /bar (best for most aqueous solutions) for pump head A and 115×10^{-6} /bar (to suit organic solvents) for pump head B. The settings represent average values for aqueous solvents (A side) and organic solvents (B side). Therefore it is always recommended to use the aqueous solvent on the A side of the pump and the organic solvent on the B side. Under normal conditions, the default settings reduce the pressure pulsation to below 2 % of system pressure, which is sufficient for most applications. If the compressibility values for the solvents used differ from the default settings, it is recommended to change the

compressibility values accordingly. Compressibility settings can be optimized by using the values for various solvents described in [Table 10](#) on page 68. If the solvent in use is not listed in the compressibility table, when using premixed solvents and if the default settings are not sufficient for your application, the following procedure can be used to optimize the compressibility settings:

1. Start channel A of the binary pump with the required flow rate.
2. Before starting the optimization procedure, the flow must be stable. Use degassed solvent only.
Check the tightness of the system with the pressure test.
3. Your pump must be connected to an Agilent data system or Instant Pilot, the pressure- and %-ripple can be monitored with one of these instruments.
4. Start the recording device in plot mode.
5. Starting with a compressibility setting of 40×10^{-6} /bar, increase the value in steps of 10. The compressibility compensation setting that generates the smallest pressure ripple is the optimum value for your solvent composition.
6. Repeat step 1 through step 5 for the B channel of your binary pump.

Table 10: Solvent compressibility

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

¹ Please check section "Solvent Information" for compatibility to your specific LC system.

Optimizing the Performance of the Module

How to Optimize the Compressibility Compensation Setting

Solvent (pure)	Compressibility ($10^{-6}/\text{bar}$)
1-Propanol	100
Toluene	87
Water	46



6 **Diagnostics and Troubleshooting**

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

Diagnostic Features 71

User Interfaces 71

Troubleshooting With HPLC Advisor 71

Maintenance and Troubleshooting Tools of the Module 72

Overview of the Module's Indicators and Test Functions 72

Pump Leak Rate Test 72

Agilent Lab Advisor Software 74

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 74.
- 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.

Maintenance and Troubleshooting Tools of the Module

Overview of the Module's Indicators and Test Functions

Status Indicators

The module is provided with two status indicators which indicate the operational state of the module. The status indicators provide a quick visual check of the operation of the module.

Error Messages

In the event of an electronic, mechanical or hydraulic failure, the module generates an error message in the user interface. For each message, a short description of the failure, a list of probable causes of the problem, and a list of suggested actions to fix the problem are provided (see chapter Error Information).

Test Functions

A series of test functions are available for troubleshooting and operational verification after exchanging internal components (see Tests and Calibrations).

Pump Leak Rate Test

Introduction

The **Pump Leak Rate Test** is used for verifying the internal tightness of the pump and helps identifying parts which may have caused a leak.

Minimum firmware revisions:

- D.07.01

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.

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

What Are Error Messages 77

General Error Messages 78

- Timeout 78
- Shutdown 78
- Remote Timeout 79
- Lost CAN Partner 80
- Leak 81
- Leak Sensor Open 82
- Leak Sensor Short 83
- Compensation Sensor Open 84
- Compensation Sensor Short 85
- Fan Failed 86
- Open Cover 87
- ERI Messages 88

Pump Error Messages 90

- Solvent Zero Counter 90
- Pressure Above Upper Limit 90
- Pressure Below Lower Limit 91
- Pressure Signal Missing 92
- Valve Failed 93
- Missing Pressure Reading 94
- Wrong Pump Configuration 95
- Electronic Fuse of SSV Open 96
- AIV Fuse 97
- Temperature Out of Range 98
- Temperature Limit Exceeded 99
- Motor-Drive Power 100

Encoder Missing 101
Servo Restart Failed 102
Pump Head Missing 103
Index Limit 104
Index Adjustment 105
Index Missing 106
Stroke Length 107
Initialization Failed 108

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.	<ul style="list-style-type: none">• Ensure all fittings are tight.
2	Broken capillary.	<ul style="list-style-type: none">• Exchange defective capillaries.
3	Loose or leaking purge valve, inlet valve, or outlet valve.	<ul style="list-style-type: none">• Ensure pump components are seated correctly. If there are still signs of a leak, exchange the appropriate seal (purge valve, inlet valve, outlet valve).
4	Defective pump seals.	<ul style="list-style-type: none">• Exchange the pump seals.

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 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.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
2	Foam not activating the sensor.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
3	Defective sensor or main board.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
4	Rear of the module is exposed to strong direct sunlight.	<ul style="list-style-type: none">• Ensure that the rear of module is not directly exposed to strong sunlight.

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.

Pump Error Messages

Solvent Zero Counter

Error ID: 2055

The error message is triggered if the remaining volume in a solvent bottle falls below the set limit.

Probable cause		Suggested actions
1	Volume in bottle below specified volume.	<ul style="list-style-type: none">• Refill bottles and reset solvent counters.
2	Incorrect setting.	<ul style="list-style-type: none">• Make sure the set solvent volume matches the actual bottle filling• Make sure the set solvent volume matches the actual bottle filling and set the shutoff limit to a reasonable value (e.g. 100 mL for 1 L bottles).

Pressure Above Upper Limit

Error ID: 2014, 2500

Probable cause		Suggested actions
1	Upper pressure limit set too low.	<ul style="list-style-type: none"> Ensure the upper pressure limit is set to a value suitable for the analysis.
2	Blockage in the flowpath (after the pressure sensor).	<ul style="list-style-type: none"> Check for blockage in the flow path. The following components are particularly subject to blockage: inline filter frit, needle (autosampler), seat capillary (autosampler), sample loop (autosampler), column frits and capillaries with small internal diameters (e.g. 50 ID). Check for blockage in the flow path. The following components are particularly subject to blockage: inline filter frit, needle (autosampler), seat capillary (autosampler), sample loop (autosampler), column frits and capillaries with small internal diameters (e.g. 50 µm ID).
3	Defective pressure sensor.	<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.

Pressure Below Lower Limit

Error ID: 2015, 2501

The system pressure has fallen below the lower pressure limit.

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

Pressure Signal Missing

Error ID: 2016

The pressure signal is missing.

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

Probable cause		Suggested actions
1	Pressure sensor disconnected.	• Please contact your Agilent service representative.
2	Defective pressure sensor.	• Please contact your Agilent service representative.

Valve Failed

Error ID: 2040

Valve 0 Failed: valve A1

Valve 1 Failed: valve A2

Valve 2 Failed: valve B2

Valve 3 Failed: valve B1

One of the solvent selection valves in the module failed to switch correctly.

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

Probable cause		Suggested actions
1	Solvent selection valve disconnected.	• Please contact your Agilent service representative.
2	Connection cable (inside instrument) not connected.	• Please contact your Agilent service representative.
3	Connection cable (inside instrument) defective.	• Please contact your Agilent service representative.
4	Solvent selection valve defective.	• Exchange the solvent selection valve.

Missing Pressure Reading

Error ID: 2054

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

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

Probable cause		Suggested actions
1	Pressure sensor disconnected.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
2	Defective pressure sensor.	<ul style="list-style-type: none">• Please contact your Agilent service representative.
3	Defective mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Wrong Pump Configuration

Error ID: 2060

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

The binary pump is assigned its configuration at the factory. If the active inlet valve and pump encoder of channel B are disconnected, and the binary pump is rebooted, the error message is generated.

Probable cause		Suggested actions
1	Active-inlet valve and pump encoder of channel B disconnected.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Electronic Fuse of SSV Open

Error ID: 2049

Valve Fuse 0: Channels A1 and A2

Valve Fuse 1: Channels B1 and B2

One of the solvent-selection valves in the module has drawn excessive current causing the selection-valve electronic fuse to open.

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

AIV Fuse

Error ID: 2044

Inlet-Valve Fuse 0: Pump channel A

Inlet-Valve Fuse 1: Pump channel B

One of the active-inlet valves in the module has drawn excessive current causing the inlet-valve electronic fuse to open.

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

Temperature Out of Range

Error ID: 2517

Temperature Out of Range 0: Pump channel A

Temperature Out of Range 1: Pump channel B

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

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

Probable cause		Suggested actions
1	Defective mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Temperature Limit Exceeded

Error ID: 2517

Temperature Limit Exceeded 0: Pump channel A

Temperature Limit Exceeded 1: Pump channel B

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

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

Probable cause		Suggested actions
1	High friction (partial mechanical blockage) in the pump drive assembly.	
2	Partial blockage of the flowpath in front of the pressure sensor.	<ul style="list-style-type: none">• Ensure the outlet valve is not blocked.
3	Defective pump drive assembly.	<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.

Motor-Drive Power

Error ID: 2041, 2042

Motor-Drive Power: Pump channel A

B: Motor-Drive Power: Pump channel B

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

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

Probable cause		Suggested actions
1	Flow path blockage in front of the pressure sensor.	• Ensure the capillaries and frits between the pump head and pressure sensor inlet are free from blockage.
2	Blocked (passive or active) inlet valve.	• Exchange the (passive or active) inlet valve.
3	Blocked outlet valve.	• Exchange the outlet valve.
4	High friction (partial mechanical blockage) in the pump drive assembly.	
5	Defective pump drive assembly.	• Please contact your Agilent service representative.
6	Defective mainboard.	• Please contact your Agilent service representative.
7	Restriction capillary blocked at pre-mixing union.	• Exchange restriction capillary.

Encoder Missing

Error ID: 2046, 2050, 2510

Encoder Missing: Pump channel A

B: Encoder Missing: Pump channel B

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

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

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

Servo Restart Failed

Error ID: 2201, 2211

Servo Restart Failed: Pump channel A

B: Servo Restart Failed: Pump channel B

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

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

Probable cause		Suggested actions
1	Mechanical blockage of the module.	
2	Disconnected or defective cable.	• Please contact your Agilent service representative.
3	Blocked (passive or active) inlet valve.	• Exchange the (passive or active) inlet valve.
4	Defective pump drive assembly.	• Please contact your Agilent service representative.
5	Defective mainboard.	• Please contact your Agilent service representative.

Pump Head Missing

Error ID: 2202, 2212

Pump Head Missing: Pump channel A

B: Pump Head Missing: Pump channel B

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

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

Probable cause		Suggested actions
1	Pump head not installed correctly (screws not secured, or pump head not seated correctly).	<ul style="list-style-type: none">• Install the pump head correctly. Ensure nothing (e.g. capillary) is trapped between the pump head and body.
2	Broken piston.	<ul style="list-style-type: none">• Exchange the piston.

Index Limit

Error ID: 2203, 2213

Index Limit: Pump channel A

B: Index Limit: Pump channel B

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

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

Probable cause		Suggested actions
1	Irregular or sticking drive movement.	<ul style="list-style-type: none">Remove the pump head, and examine the seals, pistons, and internal components for signs of wear, contamination or damage. Exchange components as required.
2	Defective pump drive assembly.	<ul style="list-style-type: none">Please contact your Agilent service representative.

Index Adjustment

Error ID: 2204, 2214

Index Adjustment: Pump channel A

B: Index Adjustment: Pump channel B

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

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

Probable cause		Suggested actions
1	Irregular or sticking drive movement.	<ul style="list-style-type: none">Remove the pump head, and examine the seals, pistons, and internal components for signs of wear, contamination or damage. Exchange components as required.
2	Defective pump drive assembly.	<ul style="list-style-type: none">Please contact your Agilent service representative.

Index Missing

Error ID: 2205, 2215, 2505

Index Missing: Pump channel A

B: Index Missing: Pump channel B

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

During initialization, the first piston is moved to the mechanical stop. After reaching the mechanical stop, the piston reverses direction until the encoder index position is reached. If the index position is not recognized within a defined time, the error message is generated.

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

Stroke Length

Error ID: 2206, 2216

Stroke Length: Pump channel A

B: Stroke Length: Pump channel B

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

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

Probable cause		Suggested actions
1	Defective pump drive assembly.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Initialization Failed

Error ID: 2207, 2217

Initialization Failed: Pump channel A

B: Initialization Failed: Pump channel B

The module failed to initialize successfully within the maximum time window.

A maximum time is assigned for the complete pump-initialization cycle. If the time is exceeded before initialization is complete, the error message is generated.

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

8 Maintenance

It is necessary to perform periodic inspection of the instrument to ensure its safe use. It is possible to have these periodic inspections performed by Agilent service representatives on a contractual basis. For information regarding the maintenance inspection contract, contact your Agilent representative.

Safety Information Related to Maintenance 112

Introduction to Maintenance 114

Overview of Maintenance and Simple Repairs 115

Maintenance Procedures 116

Cleaning the Module 117

Remove and Install Doors 118

Exchanging the Purge Valve Frit or the Purge Valve 121

Replace the O-Ring on the Purge Valve 124

Remove the Pump Head Assembly 127

Maintenance of a Pump Head without Seal Wash 129

Maintenance of a Pump Head with Seal Wash 134

Reinstall the Pump Head Assembly 140

Seal Wear-in Procedure 142

Exchange the Seal Wash Cartridge 143

Replace Leak Handling System Parts 145

Exchange the Inlet Valve 146

Exchange the Outlet Valve 148

Exchange the Solvent Selection Valve 150

Replace the Module Firmware 154

Safety Information Related to Maintenance

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.

Introduction to Maintenance

The pump is designed for easy maintenance. The most frequent maintenance procedures such as piston seal replacement and purge valve frit exchange can be done from the front side without removing the pump from the system stack.

Overview of Maintenance and Simple Repairs

Figure 9 on page 115 shows the main user accessible assemblies of the binary pump. The pump heads and its parts require normal maintenance (for example, seal exchange) and can be accessed from the front (simple repairs). Replacement of valve or filters don't require to remove the pump from the system stack.

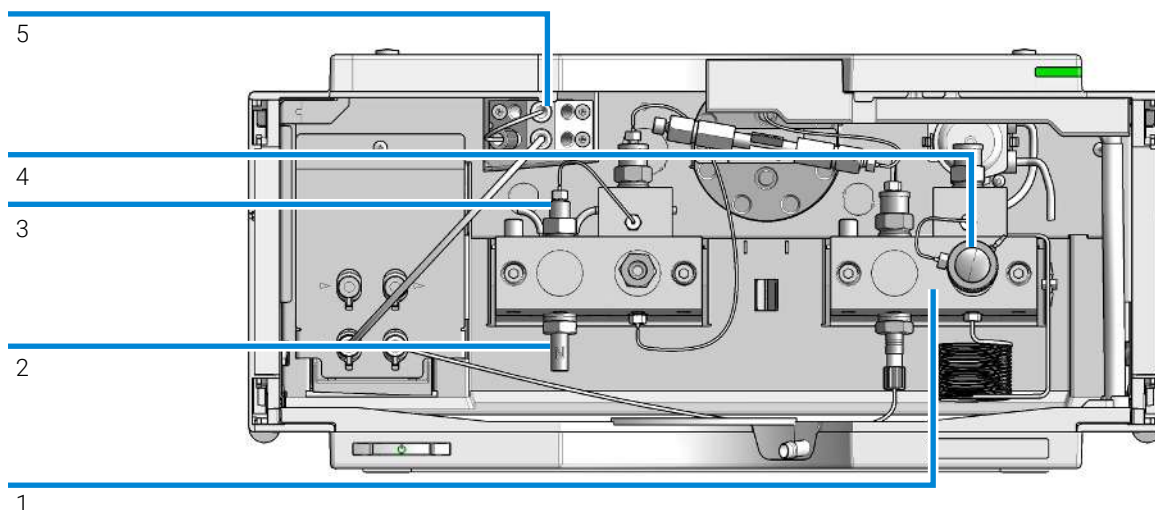


Figure 9: Overview of Maintenance and Simple Repairs

1	Pump head, see Remove the Pump Head Assembly on page 127
2	Inlet valve
3	Outlet valve, see Exchange the Outlet Valve on page 148
4	Purge valve, see Exchanging the Purge Valve Frit or the Purge Valve on page 121
5	Solvent selection valve, see Exchange the Solvent Selection Valve on page 150

Maintenance Procedures

The procedures described in this section can be done with the binary pump in place in the system stack.

Table 11: Maintenance procedures

Procedure	Typical Frequency	Notes
Exchanging the Purge Valve Frit or the Purge Valve on page 121	Yearly, or if the frit shows indication of contamination or blockage If internally leaking	A pressure drop of > 10 bar in low delay volume configuration and > 20 bar in standard configuration across the frit (5 mL/min H ₂ O with purge valve open) indicates blockage Solvent dripping out of waste outlet when valve is closed
Remove the Pump Head Assembly on page 127	During yearly maintenance	Necessary to get access to pump seals and pistons
Maintenance of a Pump Head without Seal Wash on page 129	Yearly, or if pump performance indicates seal wear	Leaks at lower pump head side, unstable retention times, pressure ripple unstable — run Valve Test for verification Seal life time shorter than normally expected — check pistons while changing the seals
Maintenance of a Pump Head with Seal Wash on page 134	Yearly, or if pump performance indicates seal wear	Only necessary when Seal Wash Option is installed. Leaks at lower pump head side, loss of wash solvent
Exchange the Outlet Valve on page 148	If internally leaking	Pressure ripple unstable, run Valve Test for verification
Exchange the Solvent Selection Valve on page 150	If internally leaking If solenoid is defective	Cross port flow Error message "Valve Failed"

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.

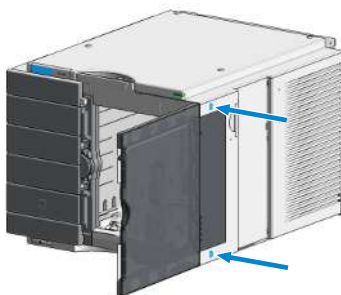
Remove and Install Doors

Parts required	Qty.	p/n	Description
	1	 5360-0017	Door 180mm left Infinity III
	1	 5360-0018	Door 180mm right Infinity III

NOTE

The figures shown in this procedure exemplarily show the Infinity III Multisampler 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 move the door in until the release buttons click into their final position.







Exchanging the Purge Valve Frit or the Purge Valve

When

- Frit – when piston seals are exchanged or when contaminated or blocked (pressure drop of > 10 bar in low delay volume configuration and > 20 bar in standard configuration across the frit at a flow rate of 5 mL/min of water with purge valve opened)
- Purge valve – if internally leaking

Tools required

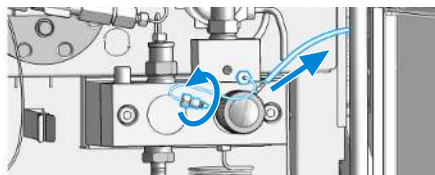
Qty.	p/n	Description
1	 8710-0510	Open-end wrench 1/4-5/16 inch
1	 8710-1924	Open-end wrench 14 mm
1		Pair of tweezers, OR
1		Toothpick

Parts required

Qty.	p/n	Description
1		Frits
1		Purge Valve
1	 5067-4728	Seal cap assembly

Preparations

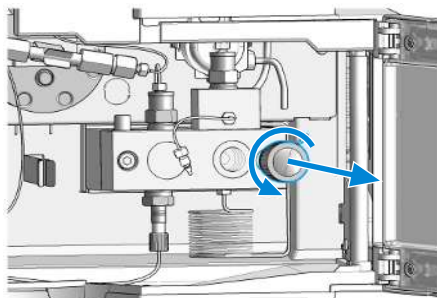
- Switch off pump at the main power switch
 - Open the doors
 - Use an optional solvent shutoff valve or lift up solvent filters in solvent reservoirs for avoiding leakages
- 1 Using a 1/4 inch wrench disconnect the pump outlet capillary from the purge valve. Disconnect the waste tube. Beware of leaking solvents due to hydrostatic pressure.



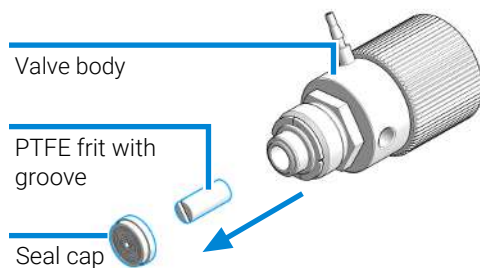
Maintenance

Exchanging the Purge Valve Frit or the Purge Valve

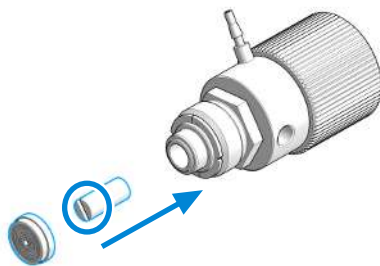
- 2 Using the 14 mm wrench, unscrew the purge valve and remove it from the purge valve holder.



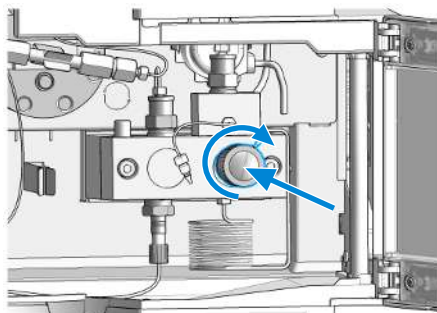
- 3 Remove the seal cap from the purge valve. Using a pair of tweezers or a toothpick, remove the frit.



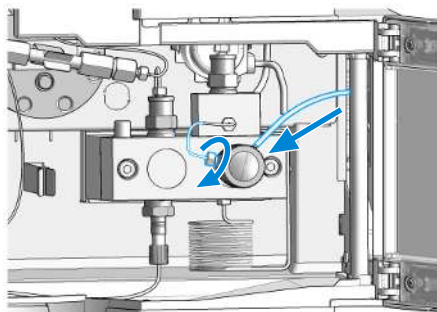
- 4 Place a new frit into the purge valve with the orientation of the frit as shown below (slit in frit points to the front). Reinstall the seal cap including the gold seal.



- 5 Insert the purge valve into the purge valve holder.



- 6 Tighten the purge valve and reconnect outlet capillary and waste tubing.





Replace the O-Ring on the Purge Valve




When

- If the original o-ring is damaged and needs to be replaced

Tools required

Qty.	p/n	Description
1	 8710-0510	Open-end wrench 1/4-5/16 inch
1	 8710-1924	Open-end wrench 14 mm
1		Pair of tweezers , OR
1		Toothpick

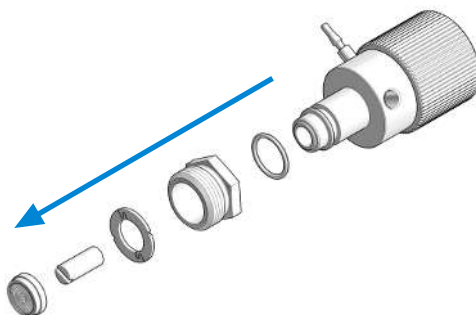
Parts required

Qty.	p/n	Description
1	 5067-6595	1260 PV O-ring FKM 5/pack
1	 01018-22707	PTFE Frit (5/Pk) (optional)
1	 5067-4728	Seal cap assembly (optional)

Preparations

- Switch off pump at the main power switch.
- Open the doors of the module.
- Use an optional solvent shutoff valve or lift up solvent filters in solvent reservoirs for avoiding leakages.
- Remove the purge valve from the pump head.

- 1 Disassemble the purge valve.



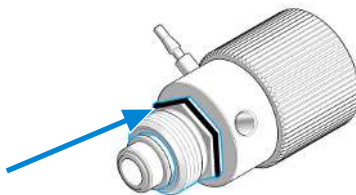
- 2 Remove the old o-ring from the purge valve.
- 3 Clean the purge valve parts.

Maintenance**Replace the O-Ring on the Purge Valve**

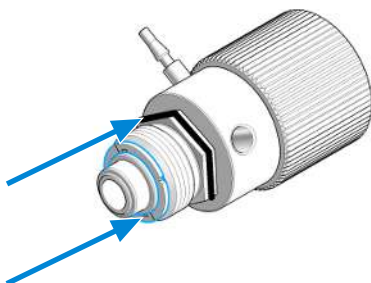
- 4 Place the new o-ring on the Screw Purge Valve.



- 5 Place the screw with o-ring on the Purge Valve Body.

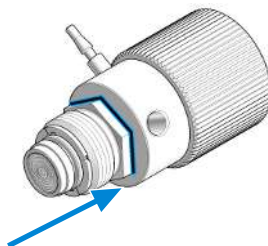


- 6 Place the mounting ring on the screw and push down the o-ring.

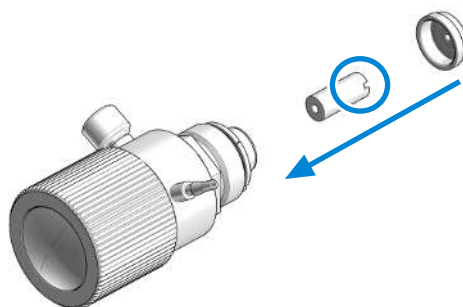


Replace the O-Ring on the Purge Valve

- 7 Push the screw up and guide the o-ring into the gap.



- 8 Place a new frit into the purge valve with the orientation of the frit as shown below (slit in frit points to the front). Reinstall the seal cap including the gold seal.

**NOTE**

Before reinstallation always check the gold seal in the seal cap. A deformed seal cap should be exchanged.

- 9 Install the purge valve to the pump. Make sure not to turn the purge valve body when the screw is fixed to the pump. The o-ring will take damage.
- 10 Run the System Pressure Test (see System Pressure Test).

Remove the Pump Head Assembly

- When**
- Exchanging pump seals
 - Exchanging pistons
 - Exchanging seals of the seal wash option

Tools required	Qty.	p/n	Description
	1	 8710-0510	Open-end wrench 1/4-5/16 inch
	1		Hexagonal key, 4 mm
	1	 5023-0240	Hex driver, ¼", slitted

- Preparations**
- Switch off the pump at the main power switch.

CAUTION

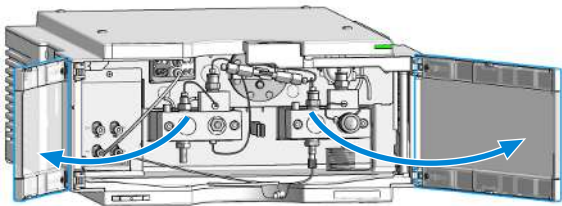
Damage of the pump drive
Starting the pump when the pump head is removed may damage the pump drive.

— Never start the pump when the pump head is removed.

NOTE

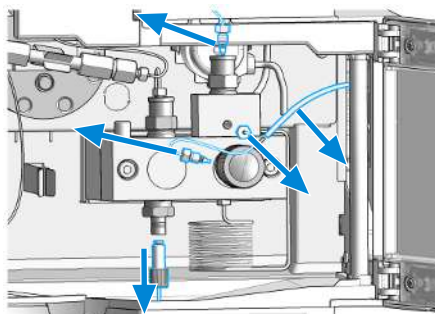
Both pump head assemblies use the same internal components. In addition, pump head A is fitted with the purge valve. The following procedure describes the removal and disassembly of pump head A (left). For pump head B (right) proceed in the same way and skip steps that deal with the purge valve.

- 1 Open the doors.

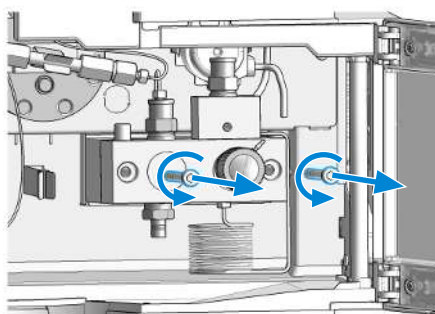


Remove the Pump Head Assembly

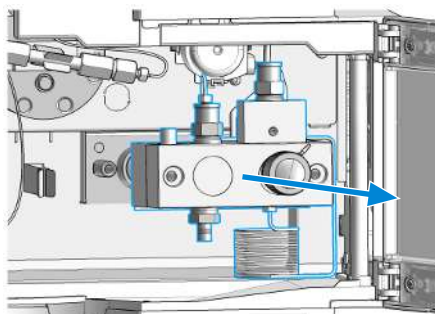
- 2 Disconnect the capillaries. Beware of leaking solvents.



- 3 Using a 4 mm hexagonal key stepwise loosen and remove the two pump head screws.



- 4 Remove the pump head assembly from the module.





Maintenance of a Pump Head without Seal Wash




When

- In case of maintenance or pump head internal leaks

Tools required

Qty.	p/n	Description
1	 8710-0510	Open-end wrench 1/4-5/16 inch
1		Hexagonal key, 4 mm
1	 01018-23702	Insert tool

Parts required

Qty.	p/n	Description
1	 5063-6589	Metering seal (pack of 2) for 100 μ L analytical head, OR
1	 0905-1420	PE seal (pack of 2)
1	 5063-6586	Sapphire piston, 100 μ L

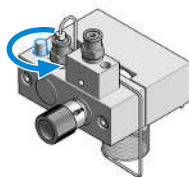
Preparations

- Switch off the pump at the main power switch.
- Open the doors of the module.
- Use an optional solvent shutoff valve or lift up solvent filters for avoiding leakages.
- Remove the Pump Head Assembly.

NOTE

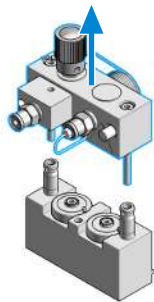
Both pump head assemblies use the same internal components. In addition, pump head A is fitted with the purge valve. The following procedure describes the removal and disassembly of pump head A (left). For pump head B (right) proceed in the same way and skip steps that deal with the purge valve.

- 1 Place the pump head on a flat surface. Loosen the lock screw (two turns).

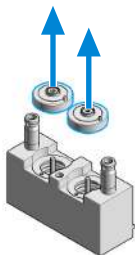


Maintenance**Maintenance of a Pump Head without Seal Wash**

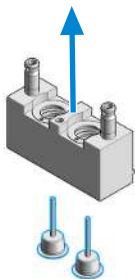
- 2 While holding the lower half of the assembly (piston housing), carefully pull the pump housing away from the piston housing.



- 3 Remove the support rings from the piston housing.



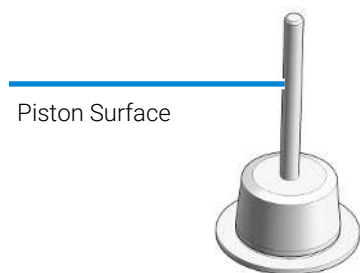
- 4 Lift the housing away from the pistons.



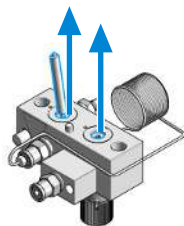
Maintenance

Maintenance of a Pump Head without Seal Wash

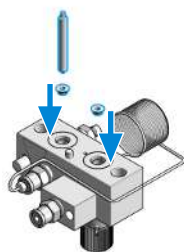
- 5 Check the piston surface and remove any deposits or layers: clean the piston surface with abrasive paper and rinse with 2-propanol. Replace piston if scratched.



- 6 Using the steel side of the insert tool, carefully remove the seal from the pump housing.

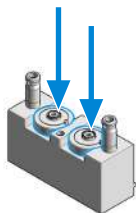


- 7 Using the plastic side of the insert tool, insert new seals into the pump head.

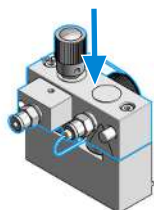


Maintenance of a Pump Head without Seal Wash

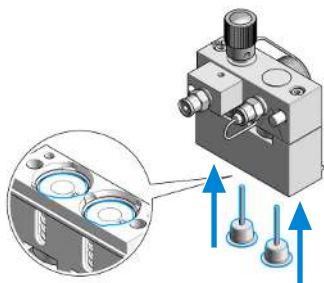
- 8 Place the support rings on the piston housing. Note the correct position of the pins on the support ring.



- 9 Place the pump housing onto the piston housing.

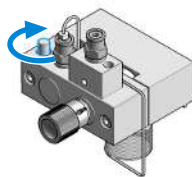


- 10 Insert the pistons and carefully press them into the seals.



Maintenance**Maintenance of a Pump Head without Seal Wash**

11 Tighten the lock screw.




12 Install the pump head, see [Reinstall the Pump Head Assembly](#) on page 140.

Maintenance of a Pump Head with Seal Wash

When

- In case of maintenance or pump head internal leaks

Tools required

Qty.	p/n	Description
1		Hexagonal key, 4 mm
1	 01018-23702	Insert tool
1		Screwdriver, small flat head

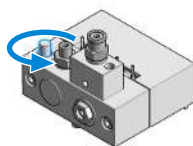
Parts required

Qty.	p/n	Description
1	 0905-1175	Wash seal (PTFE)
1	 5062-2484	Gasket, seal wash (pack of 6)
1	 5063-6586	Sapphire piston, 100 μ L
1	 0905-1719	Metering Seal, 100 μ L

Preparations

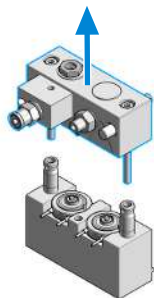
- Switch off the pump at the main power switch.
- Open the doors of the module.
- Use an optional solvent shutoff valve or lift up solvent filters for avoiding leakages.
- Remove the Pump Head Assembly.
- Remove the wash solvent tubings from the support ring inlet and outlet.

- 1 Place the pump head on a flat surface. Loosen the lock screw (two turns).

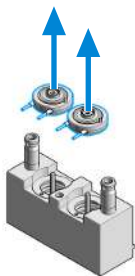


Maintenance**Maintenance of a Pump Head with Seal Wash**

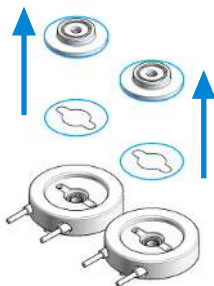
- 2 While holding the lower half of the assembly (piston housing), carefully pull the pump housing away from the piston housing.



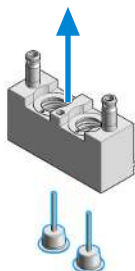
- 3 Remove the seal holder and the seal wash support rings from the piston housing.



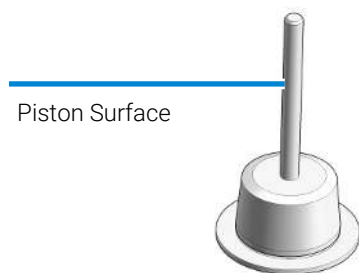
- 4 Remove the seal holder from the support ring assembly.



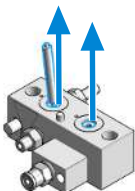
- 5 Lift the housing away from the pistons.



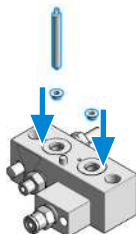
- 6 Check the piston surface and remove any deposits or layers: clean the piston surface with abrasive paper and rinse with 2-propanol. Replace piston if scratched.



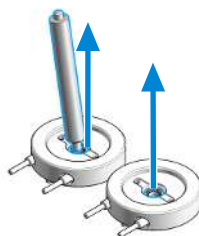
- 7 Using the steel side of the insert tool, carefully remove the seal from the pump housing.



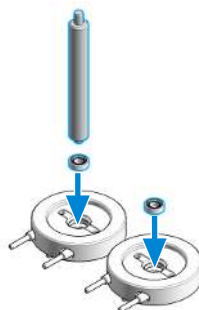
- 8 Using the plastic side of the insert tool, insert new seals into the pump head.



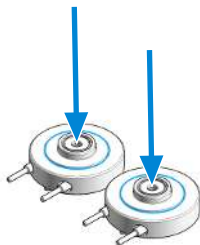
- 9 Using the steel side of the insert tool, remove the seal wash gasket and the wash seal from the support ring.



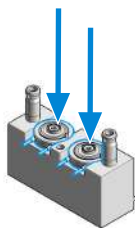
- 10 Using the plastic side of the insert tool, press the new wash seal (spring pointing upwards) into the recess of the support ring.



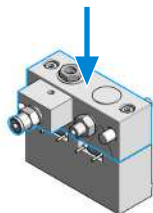
- 11 Place a seal wash gasket in the recess of the support ring. Use a matching orientation of gasket and support ring. Put the seal holder on top of the gasket.



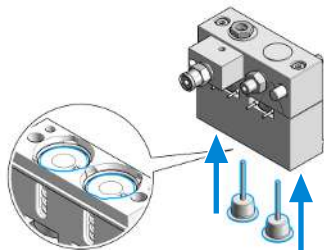
- 12 Place the support rings on the piston housing. Note the correct position of the pins on the support ring.



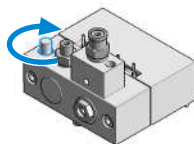
- 13 Place the pump housing onto the piston housing.



14 Insert the pistons and carefully press them into the seals.



15 Tighten the lock screw.



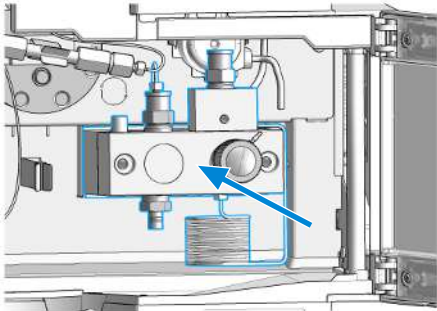
16 Install the pump head, see [Reinstall the Pump Head Assembly](#) on page 140.

Reinstall the Pump Head Assembly

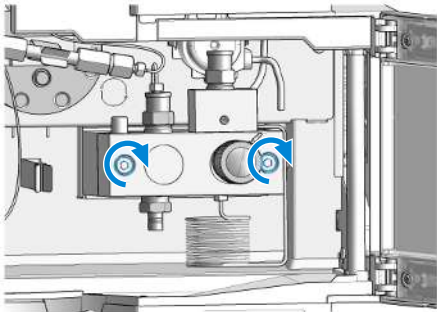
When • When reassembling the pump

Tools required	Qty.	p/n	Description
	1	 8710-0510	Open-end wrench 1/4-5/16 inch
	1		Hexagonal key, 4 mm
	1	 5023-0240	Hex driver, 1/4", slitted

- 1 Slide the pump head assembly onto the pump drive.

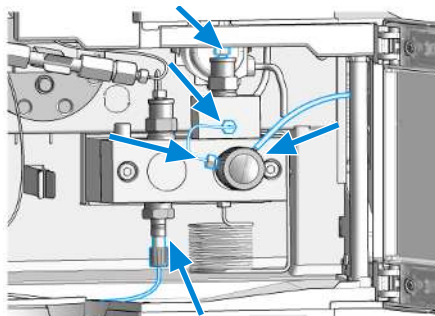


- 2 Using a 4 mm hexagonal key, tighten the pump head screws stepwise with increasing torque.

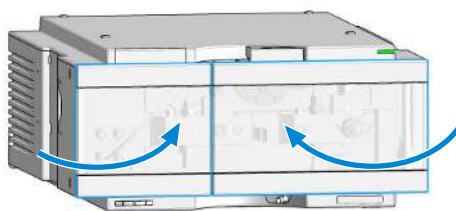


Maintenance**Reinstall the Pump Head Assembly**

- 3 Reconnect all tubings and capillaries.



- 4 Close the doors.



Seal Wear-in Procedure

NOTE

Procedure can be executed automatically in LabAdvisor

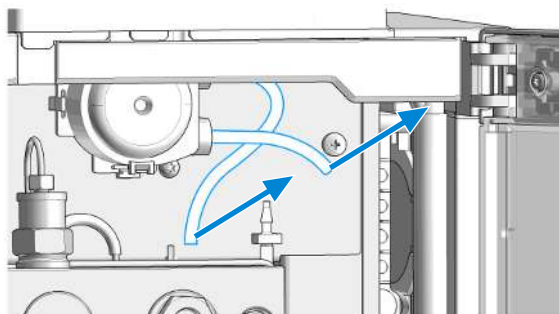
- 1 Put a bottle with 100 ml of isopropanol in the solvent cabinet and place the solvent intake filter of the pump head you want to wear in into this bottle.
- 2 Screw the 0100-1847 (PEEK adapter 1/4-28 to 10-32) onto the active inlet valve and connect the inlet tube from the bottle head directly to it.
- 3 Connect the 5022-2159 (Restriction capillary, SST 0.12 mm ID, 2 m long) to the purge valve. Connect its other end to a waste container.
- 4 Open the purge valve and purge the system for 5 min with isopropanol at a flow rate of 2 mL/min.
- 5 Close the purge valve, set the flow to a value that gives a pressure of 580 bar. Pump 15 min at this pressure to wear the seals in. The pressure can be monitored with the Instant Pilot, chromatographic data system or any other controlling device connected to your pump.
- 6 Turn OFF the pump, slowly open the purge valve to release the pressure from the system, disconnect the restriction capillary and reconnect the outlet capillary to the purge valve. Reconnect the intake tubing to the solvent selection valve and the connecting tube from the solvent selection valve (if installed) to the AIV.
- 7 Purge your system with the solvent used for your next application.

Exchange the Seal Wash Cartridge

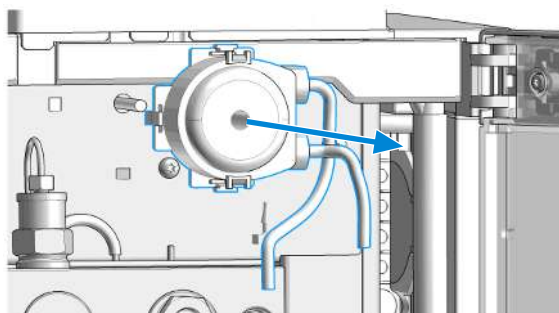
Parts required	Qty.	p/n	Description
	1	 5065-4445	Peristaltic pump cartridge

Preparations

- Switch off pump at the main power switch.
 - Open the doors.
- 1 Remove the wash solvent tubings from the support ring outlet and from the adapter leading to the waste bottle.

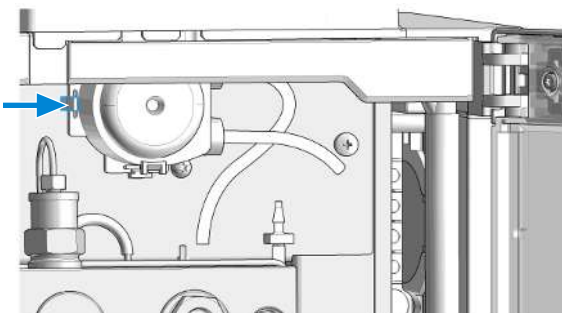
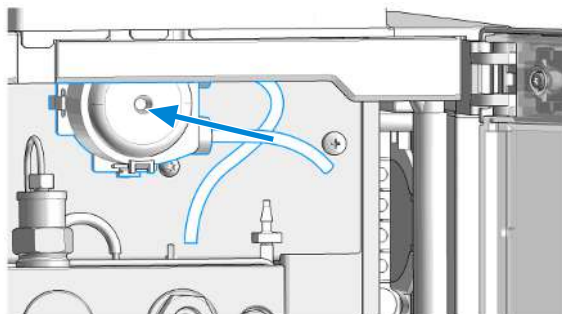


- 2 Unclip the peristaltic pump cartridge from the module housing and remove it.

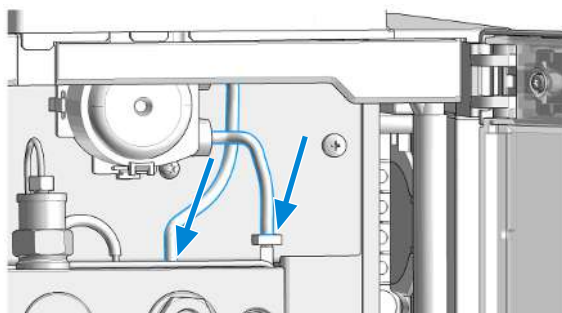


Exchange the Seal Wash Cartridge

- 3 Put the new peristaltic pump cartridge onto the rod of the pump motor and push the plastic clips into the module housing.



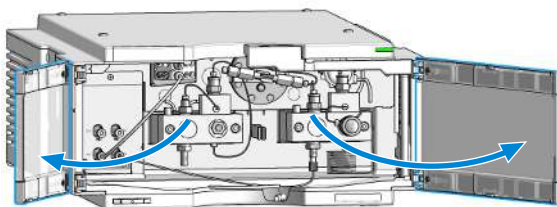
- 4 Connect the peristaltic pump tubes to the support rings outlet and to the adapter leading to the waste bottle.



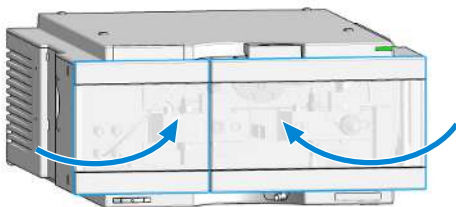
Replace Leak Handling System Parts

Parts required**Qty.****p/n****Description**1  5063-6527Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/
9 mm
approximately 85 mm required

- 1 Open the doors.





- 2 Press the Leak Adapter down and remove it together with the tubing.
- 3 Install the Leak Adapter by pressing it into the Main Cover.
- 4 Insert the Tubing (approximately 85 mm required for replacement) between Leak Adapter outlet and Leak Panel.
- 5 Close the doors.



Exchange the Inlet Valve

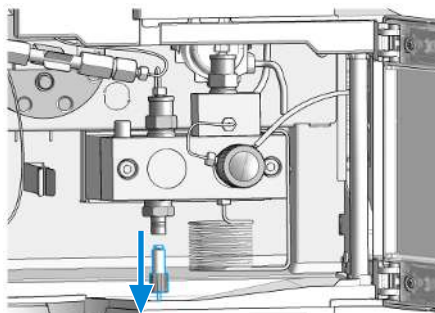
When • If internally leaking (backflow)

Tools required	Qty.	p/n	Description
	1	 8710-1924	Open-end wrench 14 mm

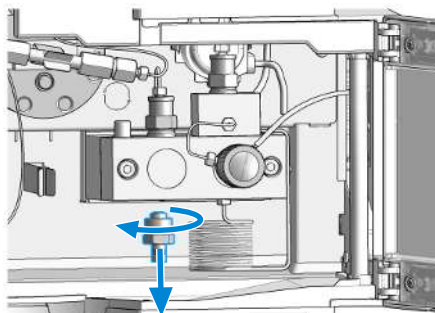
Parts required	Qty.	p/n	Description
	1	 G1312-60066	Passive inlet valve 1220/1260
	1	 G4302-60066	Inlet Valve SFC

Preparations • Switch off the pump at the main power switch

- 1 Disconnect the solvent inlet tube at the inlet valve (beware of leaking solvents).

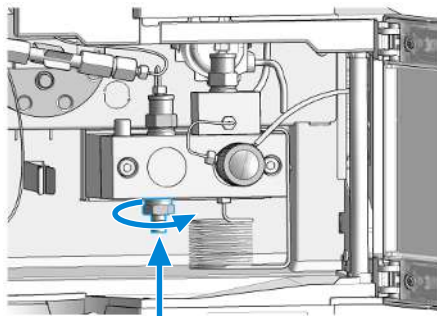


- 2 Using a 14 mm wrench, loosen the active inlet valve and remove the valve from the pump head.

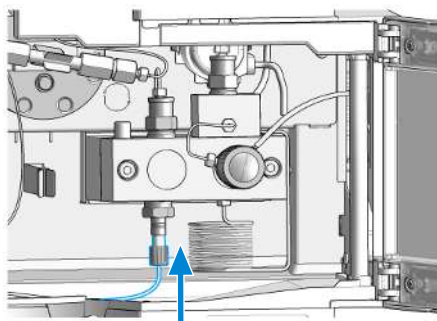


Exchange the Inlet Valve

- 3 Screw the new valve into the pump head and tighten using the 14 mm wrench. Do not overtighten.



- 4 Reconnect the inlet tube to the valve.






Exchange the Outlet Valve

When

- if leaking internally

Tools required

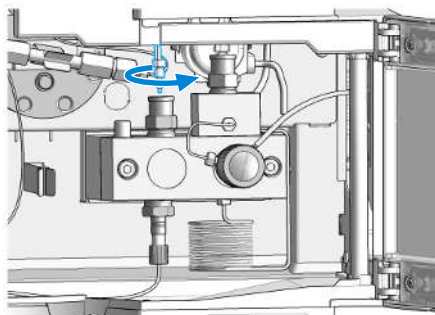
Qty.	p/n	Description
1	 8710-1924	Open-end wrench 14 mm
1	 8710-0510	Open-end wrench 1/4-5/16 inch
1	 5067-5688	Torque wrench 1 - 25 Nm with 14 mm wrench

Parts required

Qty.	p/n	Description
1	 G1312-60067	Outlet valve , OR
1	 G1312-60167	Outlet Valve Type N/SFC

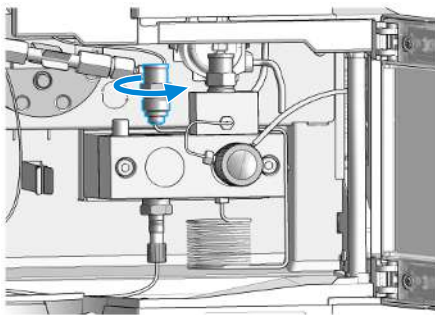
Preparations

- Switch off the pump at the main power switch
- 1 Using a ¼ inch wrench disconnect the absorber capillary from the outlet valve.



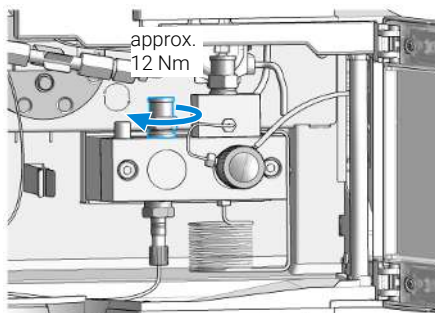
Exchange the Outlet Valve

- 2 Unscrew the valve with the 14 mm wrench and remove it from the pump body.

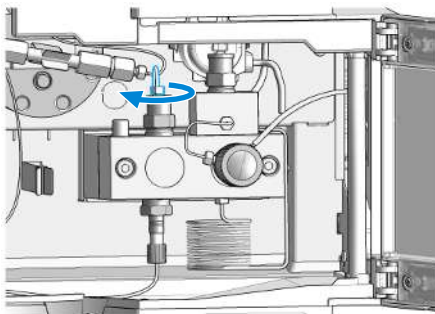
**NOTE**

Do not disassemble the outlet valve, as this can damage the valve.

- 3 Reinstall the outlet valve and tighten it using a torque wrench (approx. 12 Nm).



- 4 Reconnect the capillary.




Exchange the Solvent Selection Valve

When

- If leaking internally (crossflow between the ports), or if one of the channels is blocked

Tools required

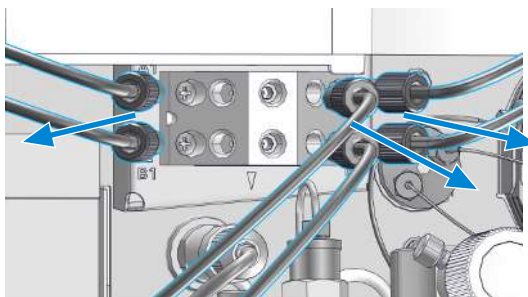
Qty.	p/n	Description
1	 8710-0899	Screwdriver Pozidrive Shaft

Parts required

Qty.	p/n	Description
1	 5067-5895	Solvent selection valve

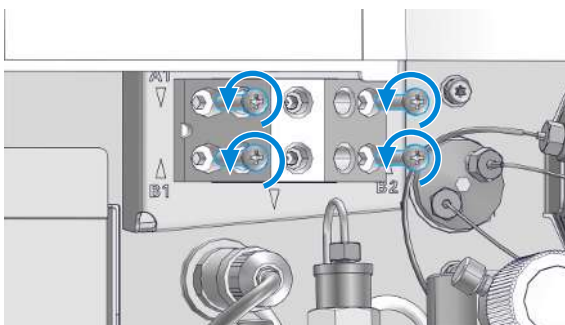
Preparations

- Switch off the pump at the main power switch
- 1 Lift solvent bottles out of the solvent cabinet and place them on the table. Disconnect the solvent tubes from the solvent selection valve and empty the tubes into the bottles. Place the bottles back into the solvent cabinet.
 - 2 Disconnect all tubings from the solvent selection valve.

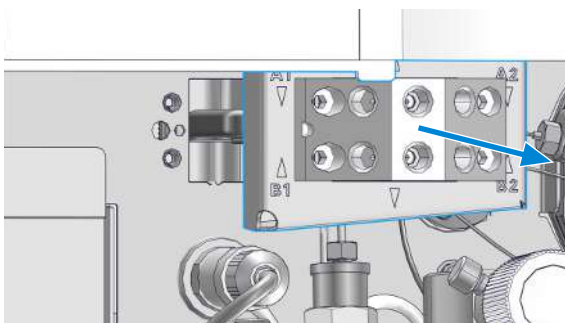


Exchange the Solvent Selection Valve

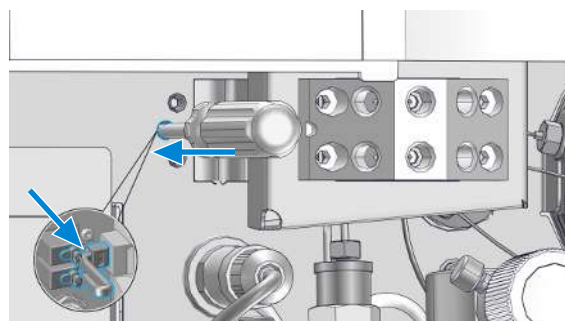
- 3 Using a screwdriver, loosen the holding screws of the valve holder.



- 4 Carefully pull the valve holder out.



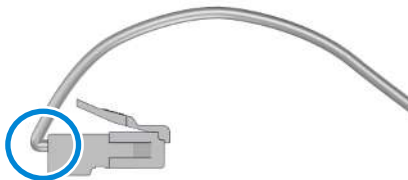
- 5 Disconnect the valve cable from the internal connection socket by pressing the fixing clip through the left side middle hole with the screwdriver.



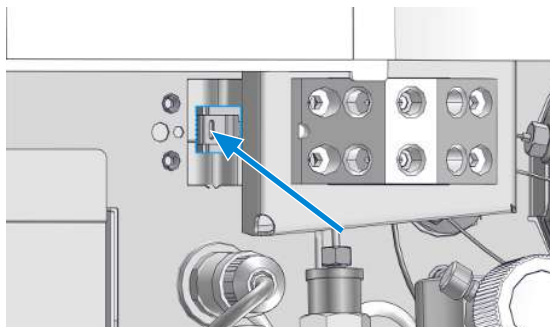
- 6 Completely remove the old valve.

Maintenance**Exchange the Solvent Selection Valve**

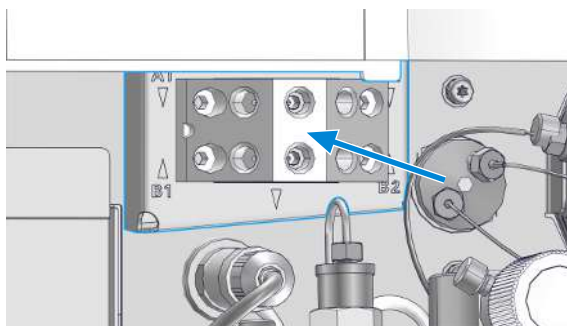
- 7 Bend the cable at the connector of the new valve.



- 8 Guide the cable and connector into the hole and push the connector into the socket.

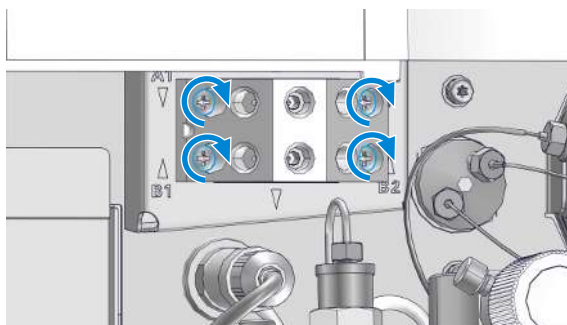


- 9 Exchange the defective solvent selection valve.

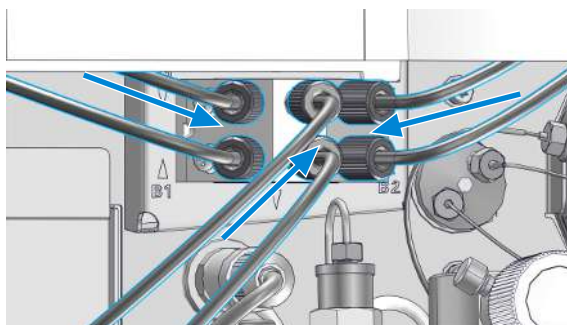


Exchange the Solvent Selection Valve

10 Tighten the screws of the valve holder.



11 Reconnect all tubings to the solvent selection valve.

**NOTE**

After an exchange of the valve it may be required to pump several mL of solvent before the flow stabilizes at a pressure ripple as low as it used to be when the system was still working properly.

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

There is no specific information for this module.

9

Parts and Materials for Maintenance

This chapter provides information on parts for maintenance.

Accessory Kit G7111-68755 156

Pump Head Assembly Channel B 157

Pump Head Assembly Channel A 159

Outlet Valve 161

Purge Valve Assembly 162

Active Seal Wash Option 163

HPLC System Tool Kit 164

Solvent Cabinet 165












Bottle Head Assembly 166

Hydraulic Path with Solvent Selection Valve 167

Cover Parts 169

Accessory Kit G7111-68755

The G7111-68755 (Accessory Kit) contains the following items:

Qty.	p/n	Description
2	 5043-1013	Tubing Clip
1	 5181-1519	CAN cable, Agilent module to module, 1 m
1	 5500-1246	Capillary ST 0.17 mm x 500 mm SI/SI
1	 5500-1217	Capillary, ST, 0.17 mm x 900 mm SI/SX
3	 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm
1	 G1311-90107	Algae note
3	 5500-1169	Tubing connector, Y-shaped, ID 6.4 mm
3	 5500-1155	Tube Connector, 90 degree, ID 6.4
1	 5043-1372	Tubing Connector Leak 3-1
2	 5043-1373	Tubing Connector Leak Cap
2	 0890-1195	Flexible sleeving 1.45 mm/2.5 mm, PTFE

Pump Head Assembly Channel B

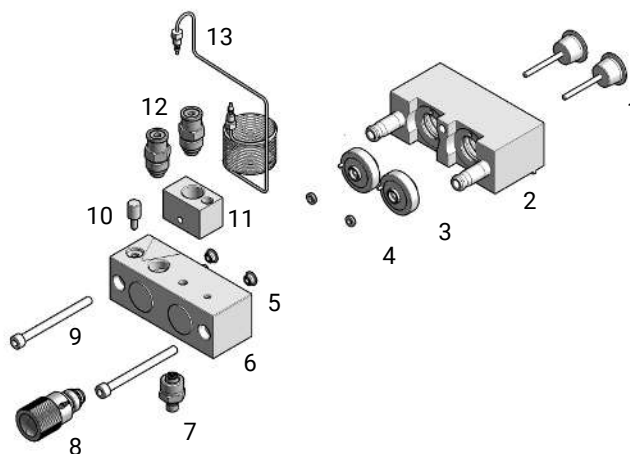





Figure 10: Pump Head Assembly Channel B

#	Qty.	p/n	Description
1	1	5067-4695	Sapphire piston (Bio-inert)
2	1	G1312-60062	Adapter, integrated, 1260
3	1	G4220-60015	Support ring including backup ring
4	1		Backup ring (part of G4220-60015)
5	1	0905-1503	Metering Seal, 100 µL
6	1	G1312-25260	Pump head body, 1260
7	1	G1312-60066	Passive inlet valve 1220/1260
8	1	G7111-60061	Purge valve
9	1	0515-2118	Screw, ST, M5 x 0.8 , 60 mm, Hex 4 mm
10	1	5042-1303	Lock screw
11	1	G4302-20000	Adapter OV SFC

Parts and Materials for Maintenance**Pump Head Assembly Channel B**

#	Qty.	p/n	Description
	1	 0515-0175	Mounting screw for manual purge valve holder, M4, 20 mm long
12	2	 G1312-60067	Outlet valve
13	1	 G1312-87300	Absorber capillary

Pump Head Assembly Channel A

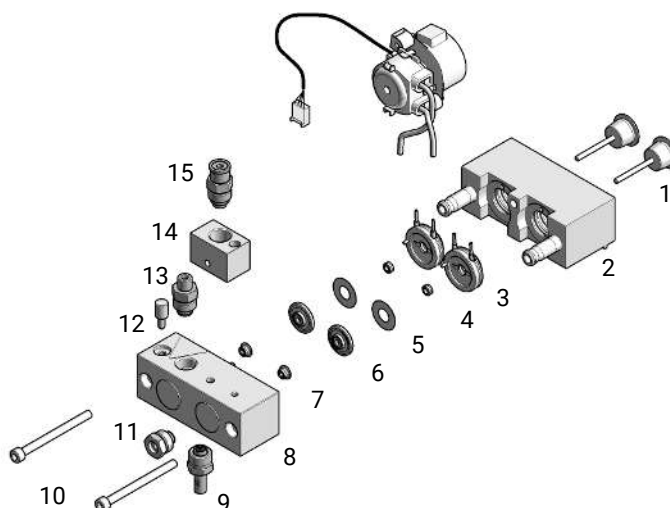










Figure 11: Pump Head Assembly Channel A

#	Qty.	p/n	Description
1	1	5067-4695	Sapphire piston (Bio-inert)
2	1	G1312-60062	Adapter, integrated, 1260
3	1	G4220-63010	Support Ring (Seal Wash)
4	1	0905-1718	Wash Seal PE
5	1	5062-2484	Gasket, seal wash (pack of 6)
6	1	G4220-60016	Seal holder including backup ring
7	1	0905-1719	Metering Seal, 100 µL
8	1	G1312-25260	Pump head body, 1260
9	1	G4302-60066	Inlet Valve SFC
10	1	0515-2118	Screw, ST, M5 x 0.8, 60 mm, Hex 4 mm
11	1	G1312-60001	Pump head adapter assembly, material: stainless steel

Parts and Materials for Maintenance

Pump Head Assembly Channel A

#	Qty.	p/n	Description
12	1	 5042-1303	Lock screw
13	1	 G4280-60026	High Pressure Filter Assembly (secondary pump head)
14	1	 G4302-20000	Adapter OV SFC
	1	 0515-0175	Mounting screw for manual purge valve holder, M4, 20 mm long
15	1	 G1312-60167	Outlet Valve Type N/SFC
	1	 5065-9978	Silicone tubing, 1 mm i.d., 3 mm o.d., 5 m, re-order number
	1	 G1316-87300	Capillary ST 0.17 mm x 90 mm S/S
	1	 G1313-87305	Restriction Capillary

Outlet Valve

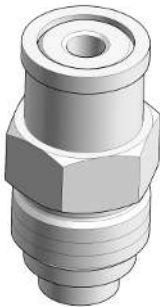




Figure 12: Outlet valve

p/n	Description
 G1312-60067	Outlet valve
 G1312-60167	Outlet Valve Type N/SFC

Purge Valve Assembly

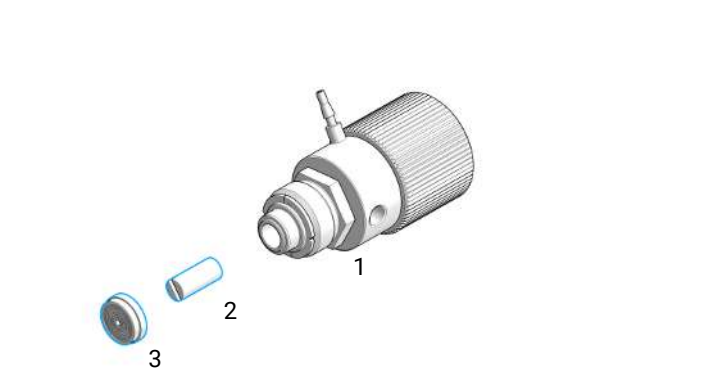


Figure 13: Purge Valve Assembly

#	Qty.	p/n	Description
1	1	G7111-60061	Purge valve
2	1	01018-22707	PTFE Frit (5/Pk)
3	1	5067-4728	Seal cap assembly
	1	5067-6595	1260 PV O-ring FKM 5/pack




















Active Seal Wash Option

The G1399B (Active Seal Wash Option kit) contains the following parts:

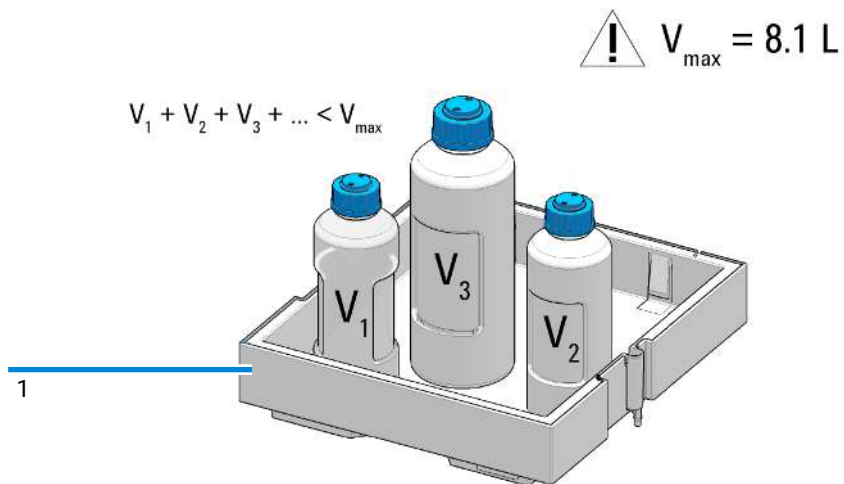
Qty.	p/n	Description
1	 5062-2484	Gasket, seal wash (pack of 6)
1	 01018-23702	Insert tool
4	 01018-60027	Support ring seal wash
2	 0515-1508	Screws for Seal Wash Pump Motor
1	 5065-9978	Silicone tubing, 1 mm i.d., 3 mm o.d., 5 m, re-order number
4	 0905-1175	Wash seal (PTFE)
1	 5063-6589	PTFE seal (pack of 2)
1	 1460-2763	Compression Spring ST
2	 1520-0260	Shock mount
1	 1540-0455	Edge protector
1	 5041-2120	Folding box
1	 5065-4445	Peristaltic pump cartridge
1	 5042-6422	Seal wash Pump Fittings x2
1	 5065-9943	Stepper Motor for the Peristaltic Pump
1	 G3010-01203	RFI Strip 10x30


HPLC System Tool Kit

G7120-68708 (InfinityLab System Tool kit) contains the following items:

p/n	Description
 9301-0411	Syringe, Plastic
 9301-1337	Syringe adapter
 0100-1710	Mounting Tool for Tubing Connections
 8710-0510	Open-end wrench 1/4-5/16 inch
 8710-1924	Open-end wrench 14 mm
 01018-23702	Insert tool
 0100-1681	Adapter syringe/seal wash tube
 8710-2394	Hex key 9/64 inch 15 cm long T-handle
 8710-1534	Wrench, open end, 4 mm
 8710-2409	Wrench open end, 5/16 – 3/8 inch
 8710-0899	Screwdriver Pozidrive Shaft
 5023-2500	Spanner double open ended SW-5
 5023-2504	Hex driver SW-4 slitted
 5810-0000	Hex driver SW-5 slitted
 5023-2502	Hex driver SW-6.35, slitted
 5023-2501	Screwdriver Torx-T10
 8720-0025	Wrench, 1/2 inch& 9/16 inch
p/n	Description
 5023-2524	Hex Key Set
p/n	Description
 01080-83202	Blank nut

Solvent Cabinet








#	p/n	Description
1	 5067-6871	Solvent Cabinet Kit

For details refer to: 01200-90150 (Usage Guideline for the Solvent Cabinet)

Bottle Head Assembly

The G7120-60007 (Bottle Head Assembly) contains:

p/n	Description
 5063-6598	Tefzel ferrules and SSL lock rings, 1/8 inch, 10/pck
 5063-6599	PPS nuts, 1/8 inch, 1/4-28 thread, 10/pck
	Wire marker
 5062-2483	Solvent tubing, 3.1 mm OD, 5 m
 5062-8517	Inlet filter adapter (4/Pk)
 5041-2168	Glass filter, solvent inlet, 20 µm

Hydraulic Path with Solvent Selection Valve

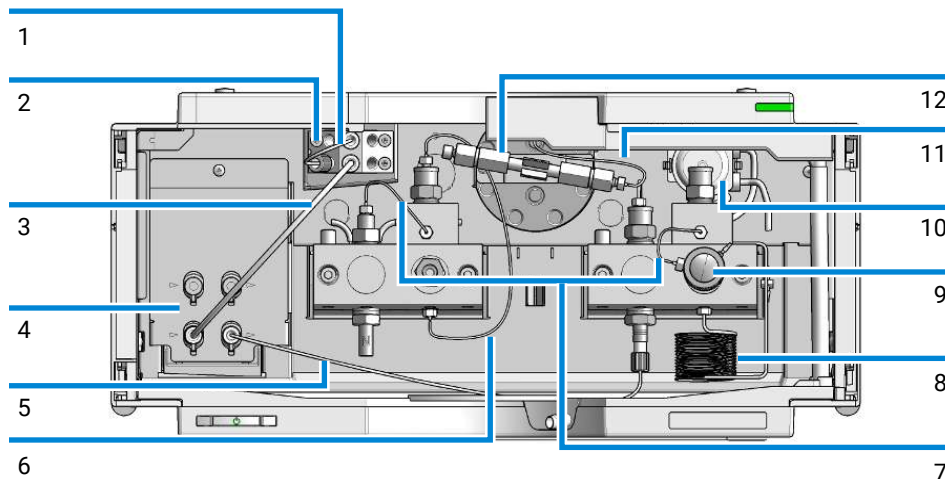











Figure 14: Hydraulic Path with Solvent Selection Valve

#	Qty.	p/n	Description
1	1	5067-4697	Solvent selection valve bridge tubing
2	1	5067-5895	Solvent selection valve
	1	5041-8365	Blank plug for unused SSV channels
3	2	G7111-60100	Solvent Tubes including labels Solvent selection valve to degasser
4	1	G7112-60070	Degasser Unit 2 Channels
5	1	G7112-67300	Connecting Tube Degasser to Channel B
6	1	G1312-87305	Capillary SSL, 0.17 x 150 mm (pressure sensor to damper)
7	2	G1316-87300	Capillary ST 0.17 mm x 90 mm S/S
8	1	G1312-87300	Absorber capillary
9	1	G7111-60061	Purge valve

Parts and Materials for Maintenance

Hydraulic Path with Solvent Selection Valve

#	Qty.	p/n	Description
10	1	 5064-5444	Peristaltic pump cartridge, silicone tubing
11	1	 G1312-87306	Capillary SSL, 0.17 x 105 mm (connections to solvent mixer)
	1	 G1312-04100	Bracket for solvent mixer
12	1	 G1312-87330	Mixer (capillary pump only)
	1	 5500-1246	Capillary ST 0.17 mm x 500 mm SI/SI
	1	 5500-1217	Capillary, ST, 0.17 mm x 900 mm SI/SX
	1	 5065-9978	Silicone tubing, 1 mm i.d., 3 mm o.d., 5 m, re-order number for seal wash option
	1	 5062-2461	Waste tube, 5 m (reorder pack)
	1	 G4301-60560	Cross and Cap. Kit for Aux Pres.Sensor

Cover Parts

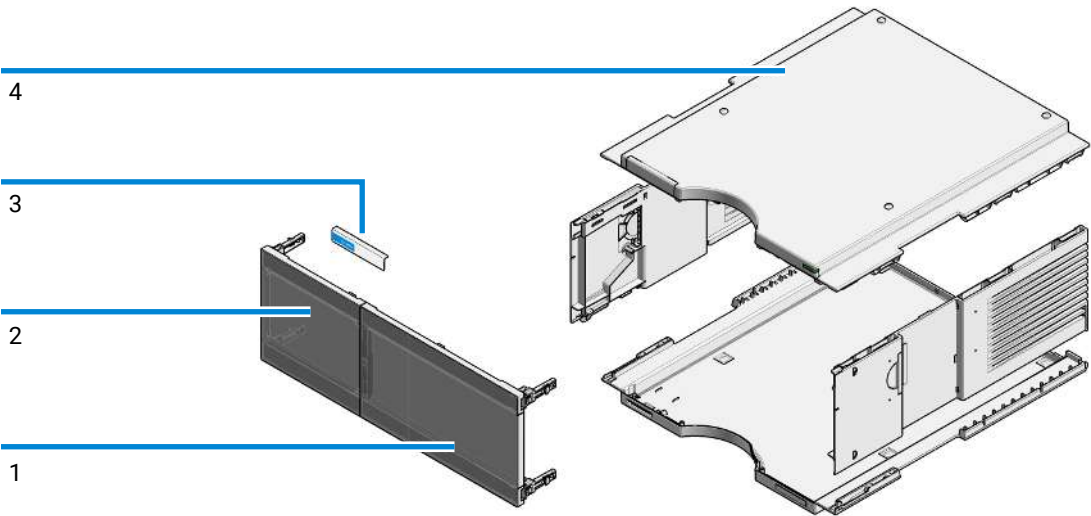






Figure 15: Cover Parts

#	p/n	Description
1	 5360-0018	Door 180mm right Infinity III
2	 5360-0017	Door 180mm left Infinity III
3	 5431-0117	Name Plate Infinity III 1260
4	 G7104-68713	Infinity II & III Cabinet Kit 180 (includes sides, bottom, top, leak adapter top and status indicator insert)

10

Identifying Cables

This chapter provides information on cables used with the modules.

Cable Overview 171

Analog Cables 173

Remote Cables 175

BCD Cables 179

CAN/LAN Cables 181

RS-232 Cables 182

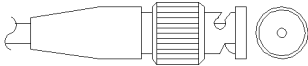
USB 183

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)

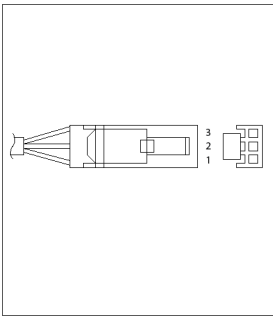
RS-232 cables	p/n	Description
	RS232-61601	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It is also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3, 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

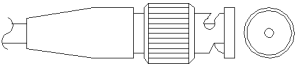


One end of these cables provides a BNC connector to be connected to Agilent modules. The other end depends on the instrument to which connection is being made.

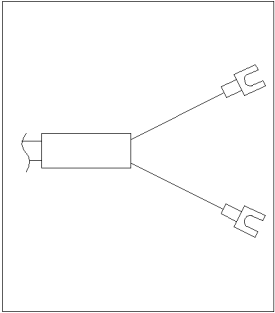
Agilent Module to 35900 A/D converters

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 +

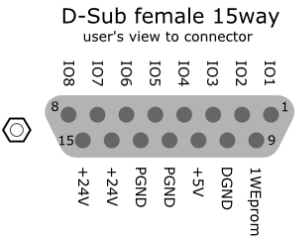
Agilent Module to General Purpose

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))

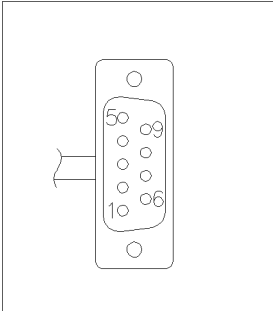
Table 12: 5188-8057 ERI to APG and RJ45

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		

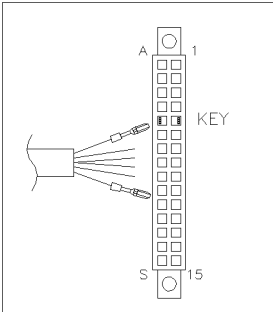


One end of these cables provides an Agilent Technologies APG (Analytical Products Group) remote connector to be connected to Agilent modules. The other end depends on the instrument to be connected to.

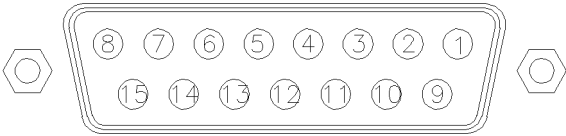
Agilent Module to Agilent 35900 A/D Converters

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

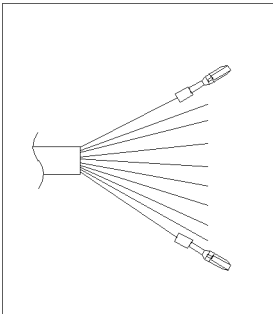
p/n 01046-60201	Wire Color	Pin Agilent module	Signal Name	Active (TTL)
	White	1	Digital ground	
	Brown	2	Prepare run	Low
	Gray	3	Start	Low
	Blue	4	Shut down	Low
	Pink	5	Not connected	
	Yellow	6	Power on	High
	Red	7	Ready	High
	Green	8	Stop	Low
	Black	9	Start request	Low

BCD Cables

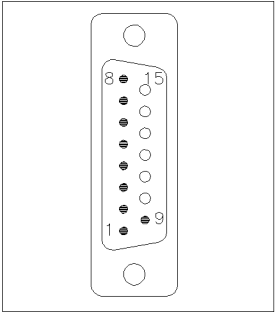


One end of these cables provides a 15-pin BCD connector to be connected to the Agilent modules. The other end depends on the instrument to be connected to

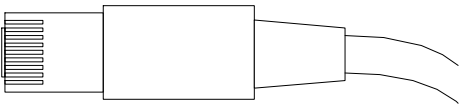
Agilent Module to General Purpose

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 185

Firmware Description 185

Electrical Connections 187

Interfaces 190

Instrument Layout 197

Early Maintenance Feedback (EMF) 197

Module-Specific Hardware Information 199

Setting the 6-bit Configuration Switch 199

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.

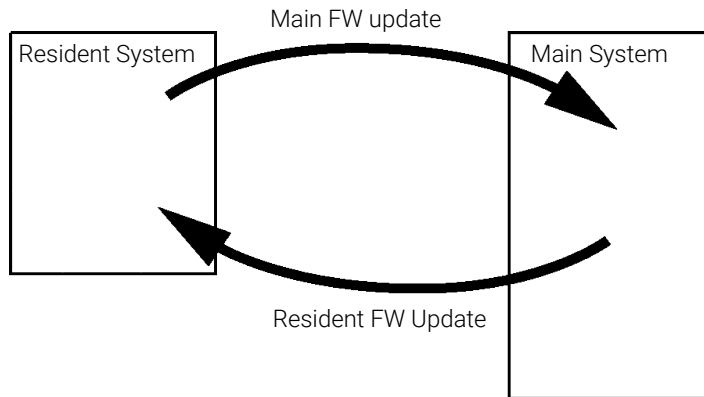


Figure 16: Firmware update mechanism

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.
- 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

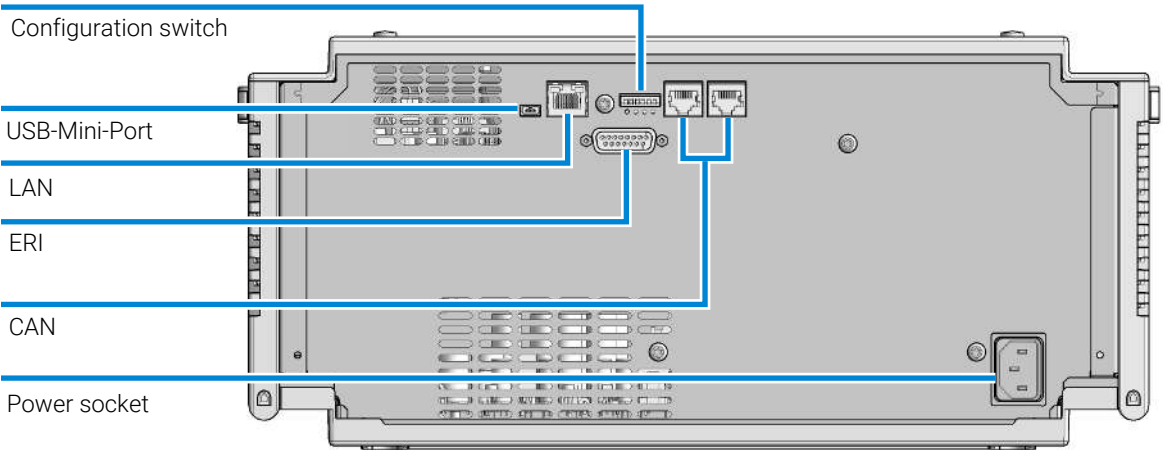


Figure 17: Rear view of the pump – electrical connections and label

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 13: 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	
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	

Module	CAN	USB	LAN (on-board)	RS-232	Analog	APG (A) / ERI (E)	Special
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 196.

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 14: 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.

Pin	Signal	Description
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.

ERI

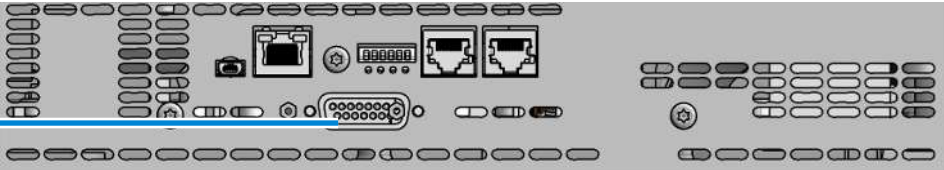


Figure 18: Location of the ERI interface

	Pin	Enhanced Remote
<div><p>D-Sub female 15way user's view to connector</p><p>IO1 IO2 IO3 IO4 IO5 IO6 IO7 IO8</p><p>1 8 15 9</p><p>1WEprom DGND +5V PGND PGND +24V +24V</p></div>	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
	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.

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.

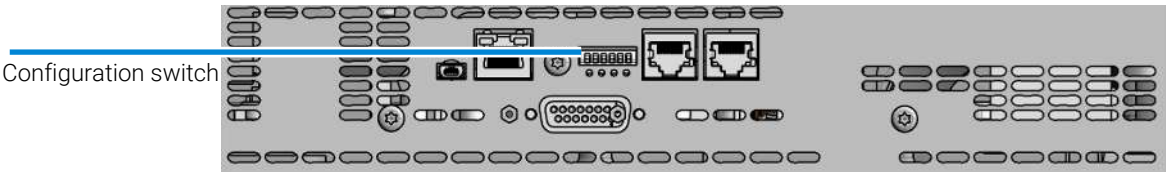


Figure 19: Location of configuration switch

Table 15: 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.

12

LAN Configuration

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

What You Have to Do First 203

TCP/IP Parameter Configuration 204

Configuration Switch 205

Initialization Mode Selection 206

Dynamic Host Configuration Protocol (DHCP) 208

General Information (DHCP) 208

Setup (DHCP) 208

Manual Configuration 211

With Telnet 211

PC and User Interface Software Setup 215

PC Setup for Local Configuration 215

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 21](#) on page 203).

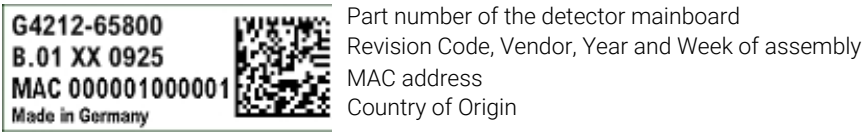


Figure 20: MAC label

- 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 21: 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 208
- 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 16](#) on page 206.

Configuration Switch

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

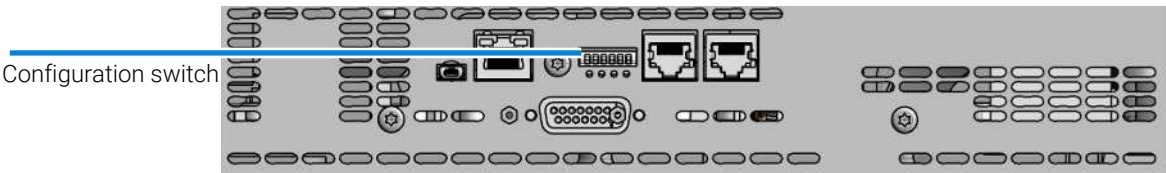


Figure 22: Location of configuration switch

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 16: 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.

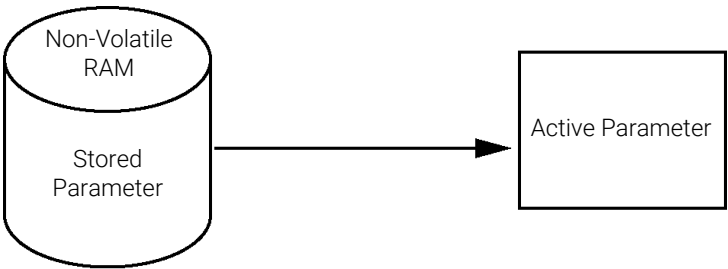


Figure 23: Using Stored (principle)

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 17](#) on page 207.

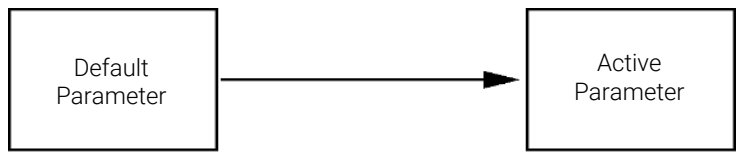


Figure 24: Using Default (principle)

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 17: 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).

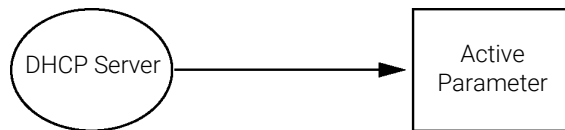


Figure 25: DHCP (principle)

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.

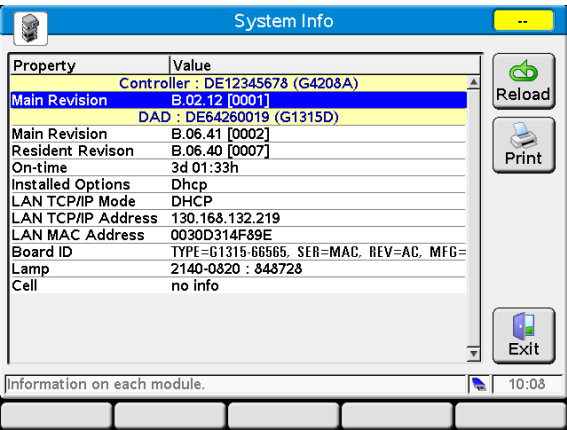


Figure 26: LAN setting on Instant Pilot

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

Table 18: 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 19: 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 208).

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>`

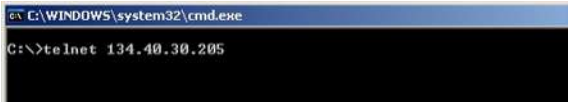


Figure 27: Telnet - Starting a session

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 205).

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

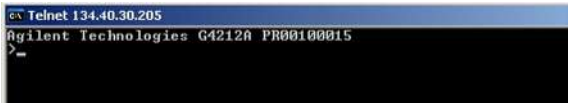


Figure 28: A connection to the module is made

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

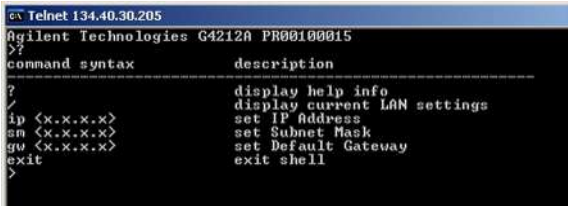
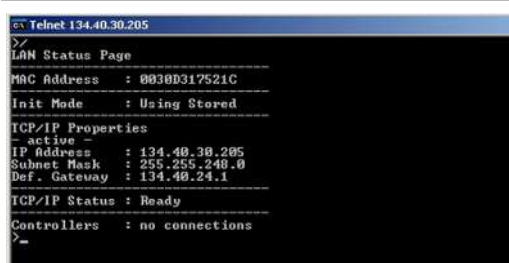


Figure 29: Telnet commands

Table 20: 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.

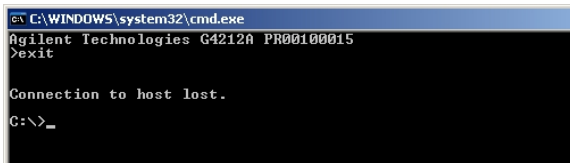


Figure 30: Closing the Telnet session

NOTE

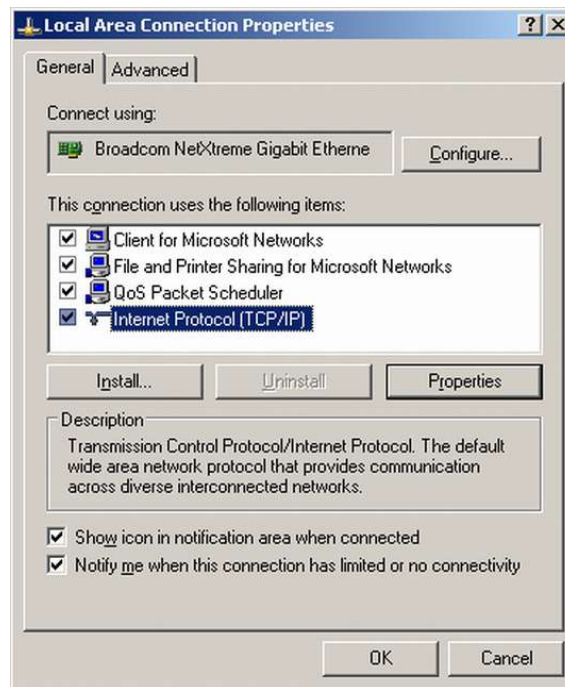
If the Initialization Mode Switch is changed now to “Using Stored” mode, the instrument will take the stored settings when the module is re-booted. In the example above it would be 192.168.254.12.

PC and User Interface Software Setup

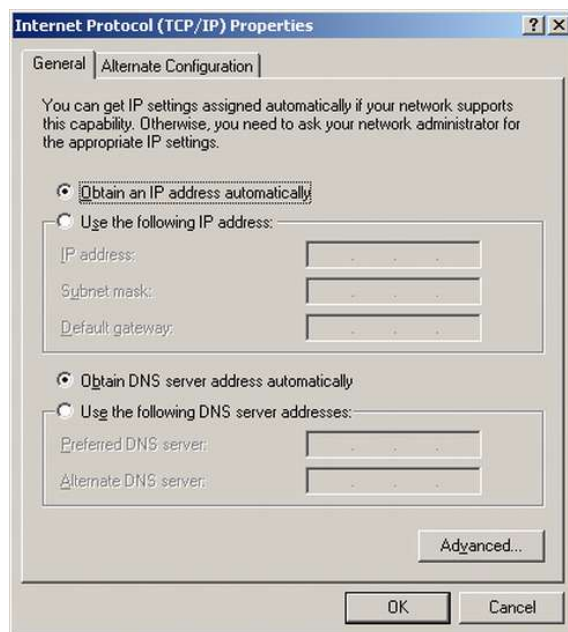
PC Setup for Local Configuration

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 17](#) on page 207).

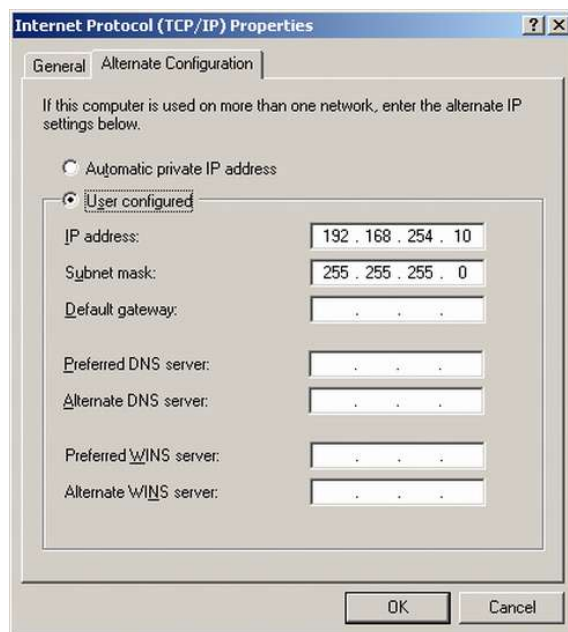
- 1 Open the Local Area Connection Properties and select **Internet Protocol (TCP/IP)**. Then click on **Properties**.



- 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 219

Safety Standards 219

General 219

Before Applying Power 220

Ground the Instrument 220

Do Not Operate in an Explosive Atmosphere 221

Do Not Remove the Instrument Cover 221

Do Not Modify the Instrument 221

In Case of Damage 221

Solvent Information 222

Algae Growth in HPLC Systems 224

Magnets 225

Safety Symbols 225

Material Information 227

Materials in Flow Path 227

General Information About Solvent/Material Compatibility 228

At-a-Glance Details About Agilent Capillaries 234**Waste Electrical and Electronic Equipment (WEEE) Directive 238****Radio Interference 239****Sound Emission 240****Agilent Technologies on Internet 241**

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 225.

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.

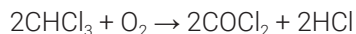
NOTE

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 [Algae Growth in HPLC Systems](#) on page 224
- 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.

Algae Growth in HPLC Systems

The presence of algae in HPLC systems can cause many problems that may be incorrectly diagnosed as instrument or application problems. Algae grow in aqueous media, preferably in a pH range from 4 to 8. Their growth is accelerated by buffers, for example phosphate or acetate. Since algae grow through photosynthesis, light will also stimulate their growth. Even in distilled water small-sized algae grow after some time.

Instrumental Problems Associated With Algae

Algae deposit and grow everywhere within the HPLC system, causing the following problems:

- Blocked solvent filters, or deposits on inlet or outlet valves, resulting in unstable flow, composition or gradient problems, or a complete failure of the pump.
- Plugging of small-pore, high-pressure solvent filters, usually placed before the injector, resulting in high system pressure.
- Blockage of PTFE frits, leading to increased system pressure.
- Plugging of column filters, giving high system pressure.
- Dirty flow cell windows of detectors, resulting in higher noise levels (since the detector is the last module in the flow path, this problem is less common).

How to Prevent and/or Reduce the Algae Problem

- Always use freshly prepared solvents, especially use demineralized water, which was filtered through 0.2 μm filters.
- Never leave mobile phase in the instrument for several days without flow.
- Always discard old mobile phase.
- Use the amber solvent bottle (9301-6526 (Solvent bottle, amber, 1000 mL)) supplied with the instrument for your aqueous mobile phase.
- If possible add a few mg/L sodium azide or a few percent organic solvent to the aqueous mobile phase.

Magnets

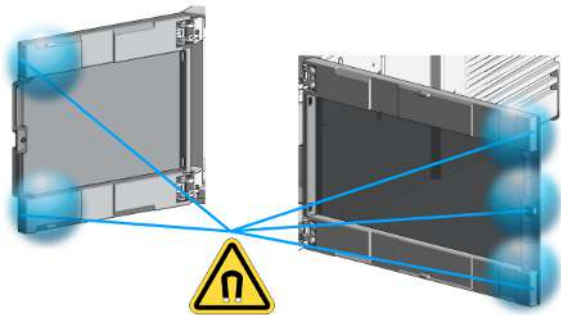








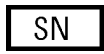







Figure 31: Magnets in doors of pumps, autosamplers, detectors, and fraction collectors

Safety Symbols

Table 21: 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.
	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.

Materials in Flow Path

Following materials are used in the flow path of this module:

Table 22: Materials in flow path

Part	Materials
Degasser chamber	TFE/PDD Copolymer, PFA (internal tubings), PEEK (inlets), FEP (tubings), ETFE (fittings)
SSV	PEEK, FFKM
Active inlet valve	SST, sapphire, ruby, ceramic, PTFE
Outlet valve	SST, gold, ruby, ZrO ₂ -based ceramic, tantalum
Adapter	SST, gold
Pump head (body)	SST
Pistons	Sapphire
Piston seals/wash seals	PTFE, SST (reversed phase) or UHMW-PE, SST (normal phase)
Pressure sensor	SST
Purge valve	SST, gold, PTFE, ceramic
Damping unit	SST, gold
Capillaries/fittings	SST
Tubings	PTFE

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)

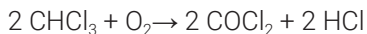
Tantalum is inert to most common HPLC solvents and almost all acids except fluoric acid and acids with free sulfur trioxide. It can be corroded by strong bases (e.g. hydroxide solutions > 10 %, diethylamine). It is not recommended for the use with fluoric acid and fluorides.

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 23: Example for a capillary description



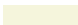










Code provided with the part	Meaing of the code
Color code: 	Material of the product is MP35N, the inner diameter is 0.20 or 0.25 mm
Capillary	The part is a connection capillary
MP35N	Material of the part is MP35N
0.25 x 80 mm	The part has an inner diameter of 0.25 mm and a length of 80 mm
SI/SI	Left 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 24](#) on page 235
- Type: [Table 25](#) on page 235
- Material: [Table 26](#) on page 236
- Dimension: [Table 27](#) on page 236
- Fittings: [Table 28](#) on page 237

Color Coding Guide

Table 24: 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 25: 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 26: 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 27: 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 28: 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.

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

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 1260 Infinity III SFC Binary Pump (G4782A).

The manual describes the following:

- Introduction
- Site Requirements and Specifications
- Using the Pump
- Optimizing Performance
- Troubleshooting and Diagnostics
- Error Information
- Test Functions and Calibration
- Maintenance
- Parts and Materials
- Cables
- Hardware Information
- LAN Configuration
- Appendix

www.agilent.com

© Agilent Technologies Inc. 2017 - 2024
Edition: 10/2024

Document No: SD-29000180 Rev. B

