

Agilent InfinityLab LC Series

1290 Infinity III Valve Drive and Valve Heads

User Manual



Notices

Document Information

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

Document No: SD-29000412 Rev. D
Edition: 10/2024

Copyright

© Agilent Technologies, Inc.
2011-2024

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 5

1 Introduction 6

Overview of the Valve Drive 7

Overview of the Agilent InfinityLab Quick Change Valves 11

Typical Applications 13

2 Site Requirements and Specifications 19

Site Requirements 20

Specifications of the 1290 Infinity III Valve Drive and Valve Heads (G1170A) 23

3 Installation 25

Unpacking the Module 26

Installing the Valve Drive 28

Installing the Valve Heads 48

Handling Leak and Waste 55

4 Using the Module 65

General Information 66

Software Configuration 70

5 Diagnostics and Troubleshooting 74

Diagnostic Features 75

Overview of the Module's Indicators and Test Functions 76

6 Error Information 77

What Are Error Messages 78

General Error Messages 79

Module Specific Error Messages 88

7 Maintenance 93

Safety Information Related to Maintenance 94

Introduction to Maintenance 96

Overview of Maintenance	97
Cleaning the Module	98
Replacing Parts of the Valve Head	99
Replacing the Fuses of the Valve Drive	101
Replace Valve Heads of the Valve Drive	105
Replace the Module Firmware	110

8 Parts and Materials for Maintenance 111

Valve Drive Parts	112
Valve Options Overview	115
Accessory Kit	117

9 Identifying Cables 118

Cable Overview	119
Analog Cables	121
Remote Cables	123
BCD Cables	127
CAN/LAN Cables	129
RS-232 Cables	130
USB	131

10 Hardware Information 132

General Hardware Information	133
Module-Specific Hardware Information	146

11 Appendix 148

General Safety Information	149
Material Information	156
At-a-Glance Details About Agilent Capillaries	162
Waste Electrical and Electronic Equipment (WEEE) Directive	166
Radio Interference	167
RFID Statement	168
Sound Emission	170
Agilent Technologies on Internet	171



In This Book

This manual covers the following Agilent InfinityLab LC Series modules:

- Agilent 1290 Infinity III Valve Drive (G1170A)
- Agilent InfinityLab Quick Change Valves



1

Introduction

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

Overview of the Valve Drive 7

Overview of the Agilent InfinityLab Quick Change Valves 11

Typical Applications 13

Dual and Multiple Column Selection 13

Sample Enrichment and Sample Cleanup (2-Position/6-Port or 2-Position/10-Port Valves) 15

Alternating Column Regeneration (2-Position/10-Port Valves Only) 16

Solvent Selection 17

Overview of the Valve Drive

The Agilent 1290 Infinity III Valve Drive (G1170A) is a part of the InfinityLab LC Series. This is a valve drive delivered without a specific valve head. To be used in a HPLC system it is required to add a valve head to this device. It is compatible with all currently available Agilent InfinityLab Quick Change Valves (see [Table 1](#) on page 11).

It includes the following features:

- Built-in power supply
- Leak handling with a leak pane and a leak sensor underneath the valve head
- User exchangeable valve heads (Agilent InfinityLab Quick Change Valves)
- Valve type and pressure detected by RFID tag
- Flexible mounting bracket, for left- or right-side mounting on LC stacks or at the Agilent Column Organizer (G1383A).

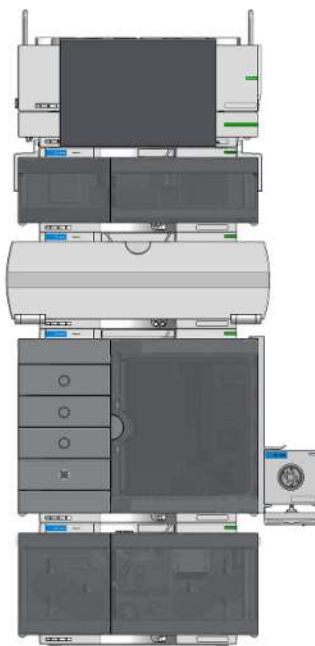


Figure 1: Mounting Examples for 1290 Infinity III Valve Drive

Introduction

Overview of the Valve Drive

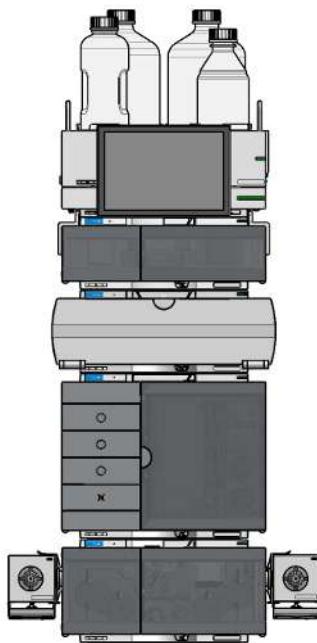


Figure 2: Mounting Examples for 1290 Infinity III Valve Drive

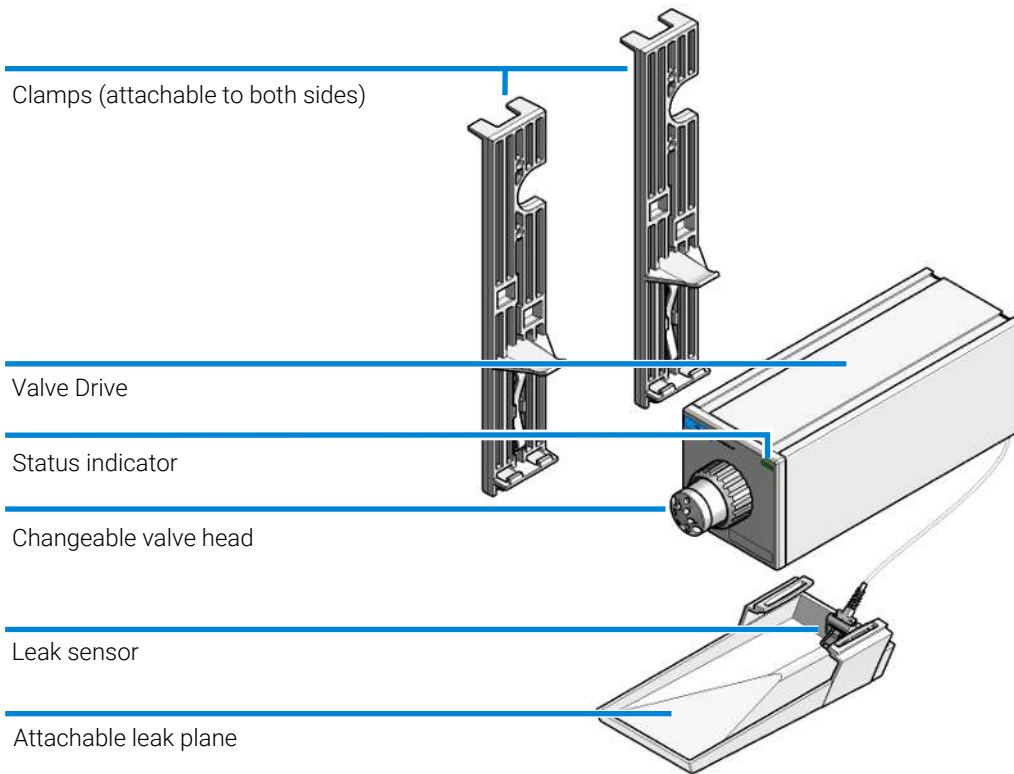


Figure 3: Overview of 1290 Infinity III Valve Drive

Introduction

Overview of the Valve Drive

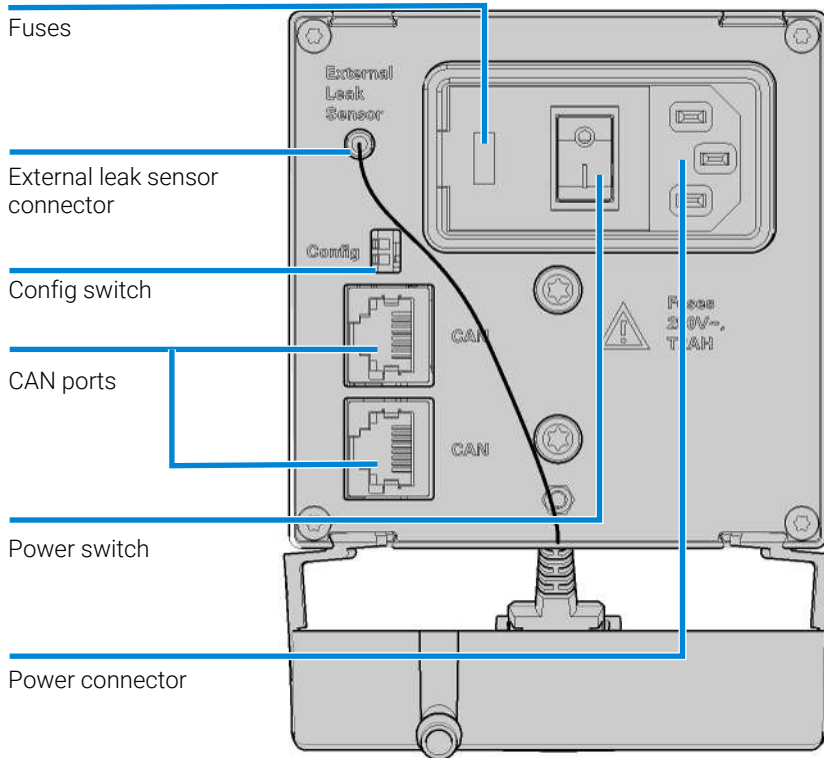


Figure 4: Rear View of 1290 Infinity III Valve Drive

Overview of the Agilent InfinityLab Quick Change Valves

The Agilent InfinityLab Quick Change Valves can be installed in the Multicolumn Thermostat (G7116A/B), the 1290 Infinity Flexible Cube (G4227A) and in the 1290 Infinity III Valve Drive (G1170A).

For bio-inert modules use bio-inert parts only!

Do not mix with bio / biocompatible parts.

For biocompatible modules use bio / biocompatible parts only!

Do not mix with bio-inert parts.



Table 1: Overview of the available Agilent Quick Change Valve Kits

Valve Kit PN	Pressure [bar]	Positions	Ports	Typical Application
G4231A	800	2	6	Any two-way switching, e.g. between two detectors, between waste and detector, between two columns
G4231C	1300	2	6	
G4232C	800	2	10	Anything a 2-position/6-port valve can do plus alternating column regeneration.
G4232D	1300	2	10	
G4234A	800	6	14	6 column selection
G4234C	1300	6	14	
G4235A	210	12	13	Solvent selection or fractionation
G4237A	800	4	10	4 column selection
G4239C	1300	8	18	8 column selection
G5631A	600	2	6	Bio-inert, same as G4231A
G5632A	600	2	10	Bio-inert, anything a 2-position/6-port valve can do plus alternation column regeneration

Introduction
Overview of the Agilent InfinityLab Quick Change Valves

Valve Kit PN	Pressure [bar]	Positions	Ports	Typical Application
G5639A	600	4	10	Bio-inert 4 column selection
G5641A	1300	2	10	Bio, anything a 2-position/6-port valve can do plus alternation column regeneration

Typical Applications

Agilent InfinityLab Quick Change Valves support a variety of valve applications. They can be mounted to an external 1290 Infinity III Valve Drive (G1170A) or a Multicolumn Thermostat (G7116A/B).

Examples of typical applications are:

- Dual and multiple column selection
- Sample enrichment and sample cleanup
- Alternating column regeneration
- Solvent selection

Dual and Multiple Column Selection

Dual Column Selection

Advantages:

- Increase productivity
- Higher instrument up-time
- Faster method scouting

Quickly change between two different stationary phases to check your separation selectivity, or use two identical stationary phases to have the second column immediately available after the first one loses efficiency, for example with complex matrices.

Multiple Column Selection

With the 6-column selection valve (G4234A/C) and the capillary kit for column selection you can set up your system for use with up to 6 columns as displayed in [Multiple Column Selection](#) on page 14. Or you can use the system with

5 columns and one flow path for flushing the system. This setup allows you to switch between columns for faster method development. The multicolumn setup might also be used, if several operators are sharing the system.

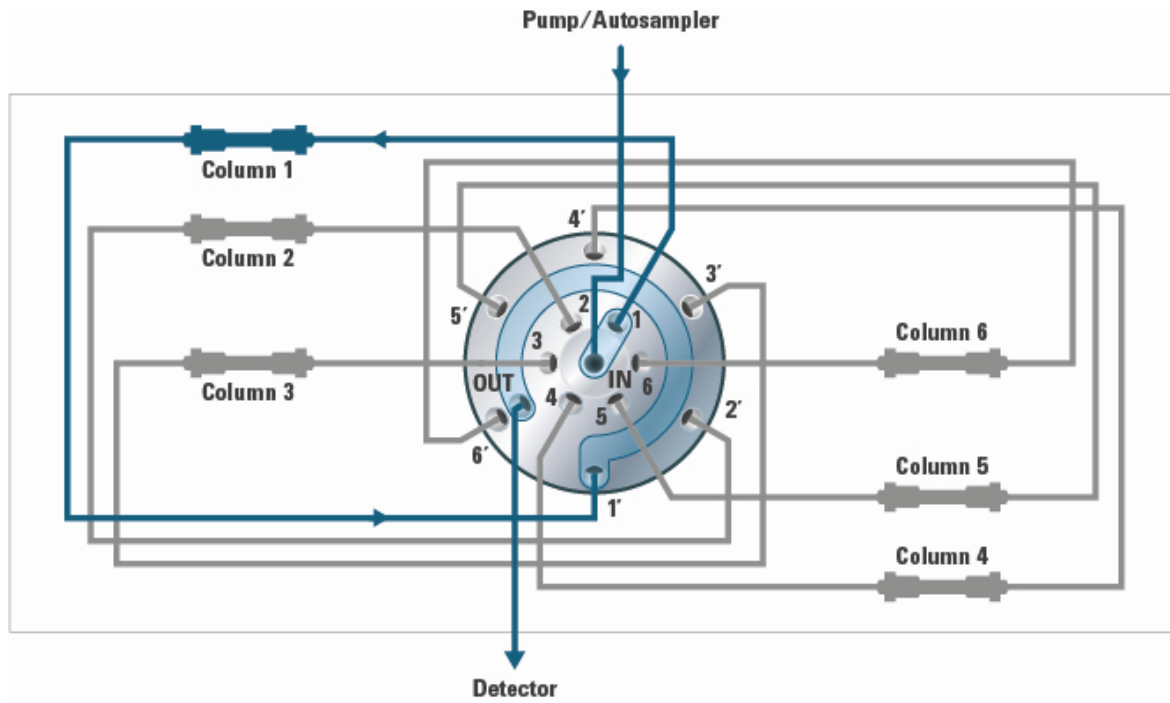
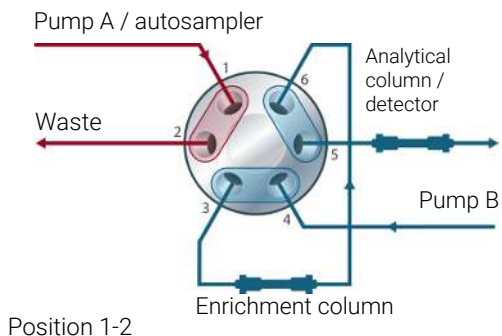
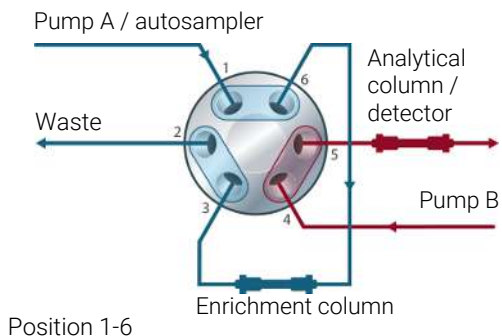


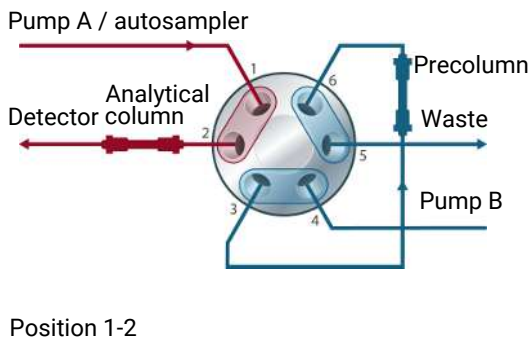
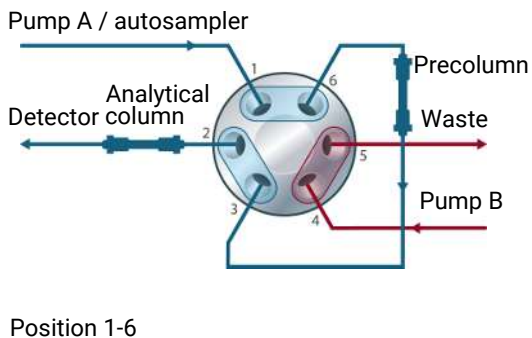
Figure 5: Multiple Column Selection (Example of Schematic Setup for 6-Column Selector)

Sample Enrichment and Sample Cleanup (2-Position/6-Port or 2-Position/10-Port Valves)

Sample Enrichment



Sample Cleanup



Advantages:

- Easy automation of sample preparation
- Higher reproducibility
- Increased productivity and sensitivity

Introduction

Typical Applications

Sample cleanup is essential for samples with complex matrices, such as food extracts and waste water. Before injection into a LC or LC/MS system, the sample matrix must be separated from the analytes of interest. Otherwise, contaminants can disrupt separation and detection or even damage the analytical column.

Enrichment Methods

Enrichment methods are the techniques of choice to obtain highest sensitivity and to remove the sample matrix. The analytes are retained and concentrated onto the precolumn, while the sample matrix is passed to waste. After the valve switch, a second pump backflushes the analytes out of the precolumn onto the separation column. This allows injection of large volumes onto the precolumn, significantly expanding sensitivity in the range of ten to several thousands.

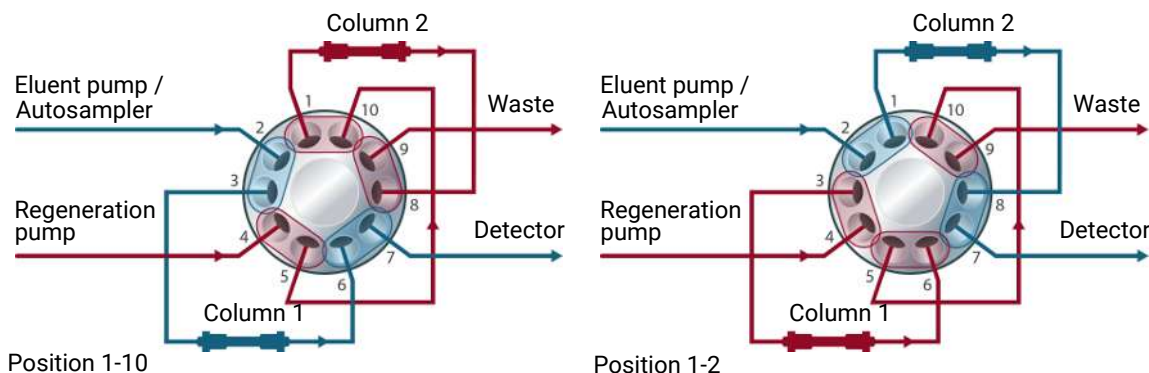
Sample Cleanup

Cleanup methods handle analytes and matrices in the opposite way to enrichment methods. Matrix components are retained on the precolumn while the analytes pass through to the separation column. After the valve switches, an extra pump backflushes the matrix components out of the precolumn to waste, while the analytes are separated on the main column. Backflushing prepares the precolumn for the next injection.

Alternating Column Regeneration (2-Position/10-Port Valves Only)

Advantages:

- High sample throughput
- Increased productivity
- High efficiency



Gradient elution is frequently used for fast separation of complex samples in LC. Since the gradient elution requires the column to regenerate before subsequent runs, an automated column regeneration system saves valuable analysis time. Agilent's InfinityLab Quick Change 2-position/10-port valve enables the simultaneous analysis of one sample on one LC column while an extra regeneration pump flushes and equilibrates a second, identical column. At the end of the run, the valve switches to the second position and the next sample is separated on the previously flushed and equilibrated column. Meanwhile, the regeneration pump flushes and equilibrates the first column. Up to 50 % of analysis time is often required to equilibrate columns. Using alternating column regeneration saves time and provides higher sample throughput.

Solvent Selection

The G4235A (12-position/13-port solvent selection valve) can be used for solvent selection (flow rate < 10 mL/min) as illustrated in [Solvent Selection](#) on page 18. It offers automated access to 12 different eluents.

Introduction

Typical Applications

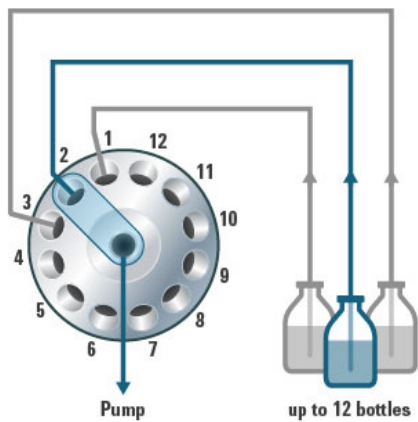


Figure 6: Solvent Selection (Schematic Setup)



2

Site Requirements and Specifications

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

Site Requirements 20

Power Considerations 20

Power Cords 21

Bench Space 22

Condensation 22

Specifications of the 1290 Infinity III Valve Drive and Valve Heads (G1170A) 23

Site Requirements

A suitable environment is important to ensure optimum 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 2](#) on page 23. 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

Inaccessible power plug.

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

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

WARNING

Hazard of electrical shock or damage of your instrumentation

can result, if the devices are connected to a line voltage higher than specified.

- Connect your instrument to the specified line voltage only.

WARNING

Electrical shock hazard

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

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

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

Power Cords

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

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

WARNING

Unintended use of power cords

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

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

WARNING

Absence of ground connection

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

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

WARNING

Electrical shock hazard

Solvents may damage electrical cables.

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

Bench Space

The module dimensions and weight (see [Table 2](#) on page 23) 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 1290 Infinity III Valve Drive and Valve Heads (G1170A)

Table 2: Physical Specifications of the 1290 Infinity III Valve Drive (G1170A)

Parameter Name	External Value	Comment
Weight	1.9 kg (4.3 lbs)	
Dimensions (height x width x depth)	90 x 90 x 300 mm (3.54 x 3.54 x 11.8 inches)	
Line voltage	100–240 V~, ±10%	Wide-ranging capability
Line frequency	50 or 60 Hz, ±5%	
Power consumption	20 VA, 4 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 (9843 ft)	
Safety standards: IEC, EN, CSA, UL	Overvoltage category II, Pollution degree 2	For indoor use only
ISM classification	ISM Group 1 Class B	According to CISPR 11

Table 3: Specifications of the Agilent InfinityLab Quick Change Valves

Kit PN	Valve PN	Description	max. Pressure [bar]	Fittings	Liquid Contact	pH Range
G5641A	5067-6682	2-position/10-port valve head, bio	1300	10-32	PEEK, MP35N	0-14 ¹
G4231A	5067-4282	2-position/6-port valve head	800	10-32	PEEK, ST	0-14 ¹
G4231C	5067-4241	2-position/6-port valve head	1300	10-32	PEEK, ST	0-14 ¹
G4232C	5067-4283	2-position/10-port valve head	800	10-32	PEEK, ST	0-14 ¹
G4232D	5067-4240	2-position/10-port valve head	1300	10-32	PEEK, ST	0-14 ¹
G4234A	5067-4284	6-position/14-port valve head	800	M4	PEEK, ST	0-14 ¹
G4234C	5067-4273	6-position/14-port valve head	1300	M4	PEEK, ST	0-14 ¹
G4235A	5067-4159	12-position/13-port valve head, bio-inert	210	10-32	PEEK, Ceramic	0-14 ¹
G4237A	5067-4279	4-position/10-port valve head	800	M4	PEEK, ST	0-14 ¹
G4239C	5067-4233	8-position/18-port valve head	1300	M4	PEEK, ST	0-14 ¹
G5631A	5067-4148	2-position/6-port valve head, bio-inert	600	10-32	PEEK, Ceramic	0-14 ¹
G5632A	5067-4132	2-position/10-port valve head, bio-inert	600	10-32	PEEK, Ceramic	0-14 ¹
G5639A	5067-4134	4-position/10-port valve head for columns selection, bio-inert	600	10-32	PEEK, Ceramic	0-14 ¹

Shipping and Storage Temperature: -40 – 80 °C

NOTE

Shipping, storing or operating this valve below 0 °C with water in the fluid passages may cause failure of the sealing surfaces.

NOTE

Operating valves above their maximum pressure limit will harm them. Please set a corresponding max. pressure limit in your Chromatographic Data System if a valve is positioned in the high pressure path.

¹ Incompatible with some mineral acids. For more information, see Solvent Information_all.

3 Installation

This chapter gives information about the installation of the valve drive and the valve heads.

Unpacking the Module 26

Damaged Packaging 26

Delivery Checklist 26

Accessory Kit 27

Installing the Valve Drive 28

Assembling and Installing the Leak Tray and Leak Sensor 28

Mounting the Valve Drive to an Instrument by Using a Valve Rail 32

Mounting the Valve Drive to a G1383A Column Organizer 36

Mounting the Valve Drive as a Part of Online Sample Manager 37

Mounting the Valve Drive to an Infinity III Module (except Multicolumn Thermostat) 39

Mounting the Valve Drive to a G7116A/B Multicolumn Thermostat 43

Installing the Valve Heads 48

Install the Valve Head and Connect Capillaries 48

Remove the Transportation Lock and the Valve Dummy 52

Handling Leak and Waste 55

Drain Connectors Installation 58

Waste Concept 63

Waste Guidance 63

Leak Sensor 64

Unpacking the Module

Damaged Packaging

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

CAUTION

"Defective on arrival" problems

If there are signs of damage, please do not attempt to install the module. Inspection by Agilent is required to evaluate if the instrument is in good condition or damaged.

- Notify your Agilent sales and service office about the damage.
- An Agilent service representative will inspect the instrument at your site and initiate appropriate actions.

Delivery Checklist

Ensure all parts and materials have been delivered with your module. The delivery checklist is shown below. For parts identification please check the illustrated parts breakdown in [Valve Drive Parts](#) on page 112. Please report any missing or damaged parts to your local Agilent Technologies sales and service office.

Table 4: Delivery Checklist

Description	Quantity
Universal Valve Drive	1
5067-4634 (Valve Rail Kit)	Optional
5067-1510 (Rail assy for column organizer)	Optional
G4800-64500 (Agilent LC Hardware Documentation Media)	1
Power cable	1

Installation
Unpacking the Module

Description	Quantity
G1369-60012 (LAN Interface Card)	Optional (required if no proper hosting module is available, see Interfaces on page 139)
G1170-68705 (Accessory Kit)	1

Accessory Kit

The 1290 Infinity III Valve Drive is shipped with G1170-68705 (Accessory Kit) (see [Accessory Kit](#) on page 117).

Installing the Valve Drive

The Valve Drive can be installed in different ways.

It can be attached to either side of your instrument with the use of the Valve Rail (InfinityLab LC Series, 1260 and 1290 Series Pumps and Detectors, for older modules order new cover kits), or it can be mounted to a G1383A column organizer by using the optional column stand mount.

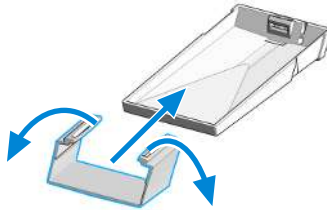
The Universal Valve Drive can also be installed on all Infinity II/III modules on either sides (5067-5685 Clamp Guide Kit). Installation of the Valve Drive on the G7116A/B requires either 5067-6138 Valve Holder Kit Right or 5067-6139 Valve Holder Kit Left.

Assembling and Installing the Leak Tray and Leak Sensor

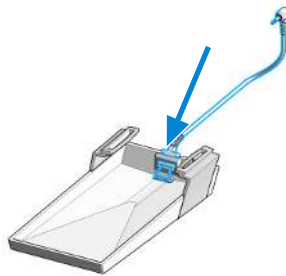
NOTE

Only connect or disconnect the leak sensor while the valve drive is powered off. The valve drive must be 'powered off' for at least 10 s to recognize any hardware changes correctly.

- 1 Attach the leak tray holder to the leak tray.



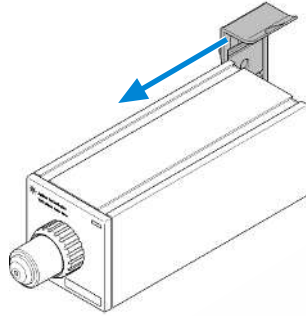
- 2 Push the leak sensor into its holder.



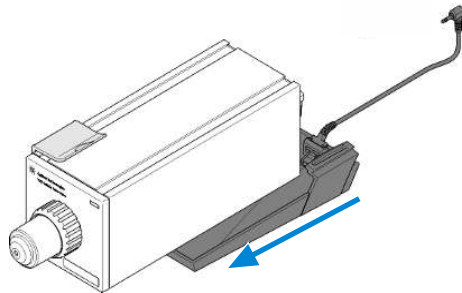
Installation

Installing the Valve Drive

- 3 Slide the first clamp on to one side of the valve drive housing.



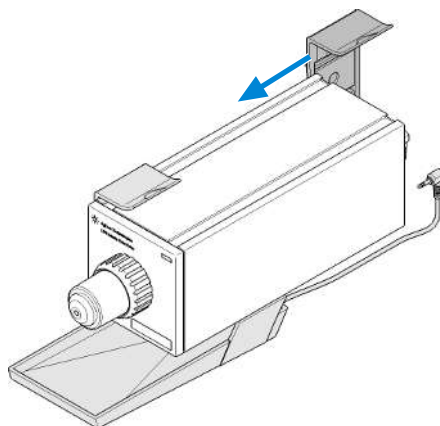
- 4 Slide the leak tray onto the bottom of the valve drive.



Installation

Installing the Valve Drive

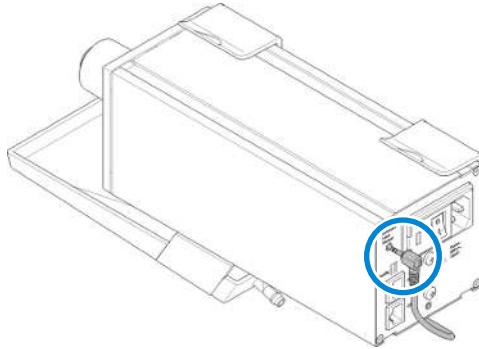
- 5 Slide the second clamp on to the same side as the first one.



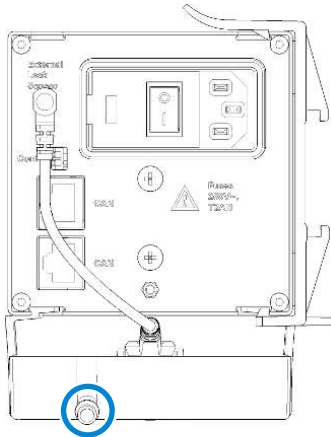
Installation

Installing the Valve Drive

- 6 Connect the leak sensor cable to the valve drive.



- 7 Attach the waste tubing to the leak plane and guide it to a proper waste container.



Mounting the Valve Drive to an Instrument by Using a Valve Rail

CAUTION

"Defective on arrival" problems

If there are signs of damage, please do not attempt to install the module. Inspection by Agilent is required to evaluate if the instrument is in good condition or damaged.

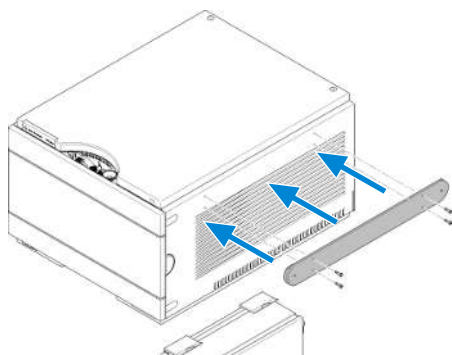
- Notify your Agilent sales and service office about the damage.
- An Agilent service representative will inspect the instrument at your site and initiate appropriate actions.

NOTE

The Installation of the leak tray and the external leak sensor is optional. If you want to attach the leak tray to the valve drive you need to slide the leak tray on to the valve drive before having both clamps attached to the valve drive.

Mounting the valve to an instrument by using a valve rail

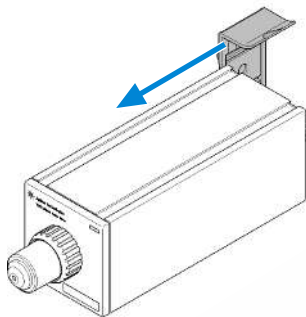
- 1 Ensure the power switch at the rear of the module is OFF and the power connector is unplugged.
- 2 Identify the marks for the screws on the side panel of the cabinet assy and screw the valve rail to the side panel of the cabinet.



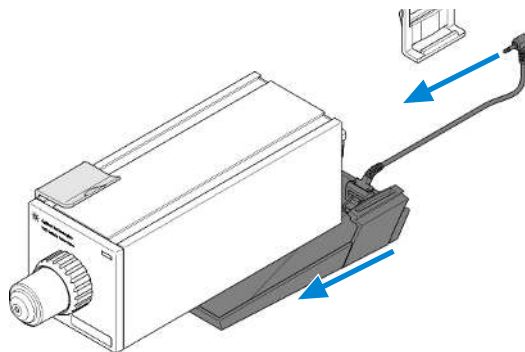
Installation

Installing the Valve Drive

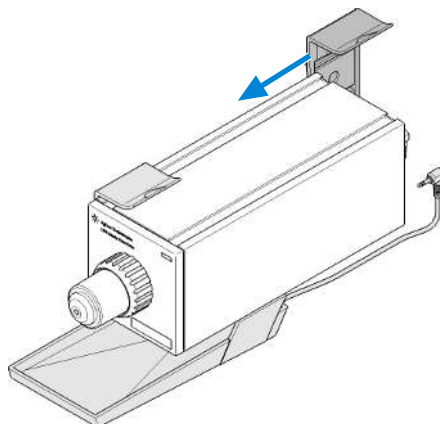
- 3 Slide the first clamp on to one side of the valve drive housing.



- 4 Install the leak tray and leak sensor to the valve.



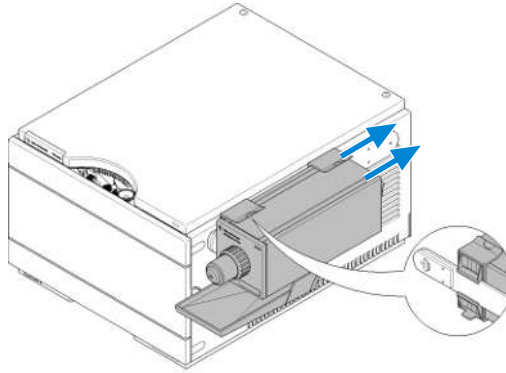
- 5 Slide the second clamp on to the same side as the first one.



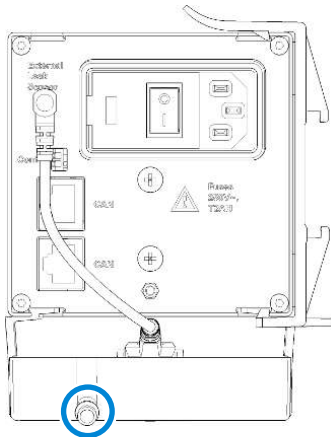
Installation

Installing the Valve Drive

- 6 Slide the valve on to the valve rail.



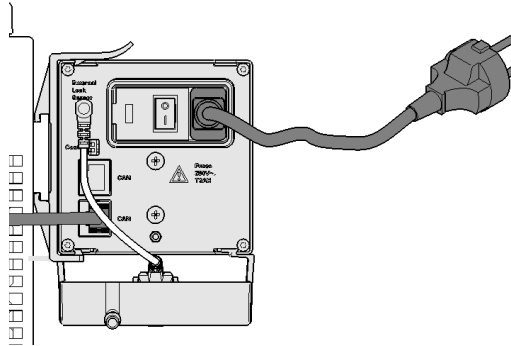
- 7 Attach the waste tubing to the leak plane and guide it to a proper waste container.



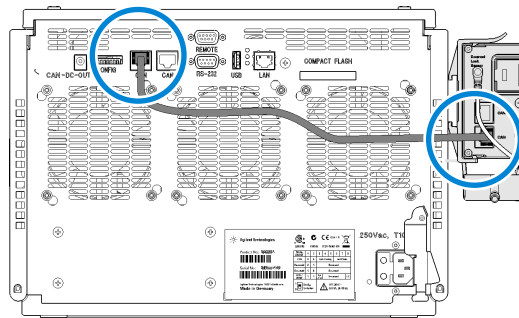
Installation

Installing the Valve Drive

- 8 Connect the power cable to the power connector at the rear of the module.



- 9 Connect the CAN interface connection.



- 10 Power on the module by switching the Power button at the rear of the module if a valve head has been already installed. Otherwise continue with [G7116B_Installation of the Valve Heads](#) on page 48.




Mounting the Valve Drive to a G1383A Column Organizer

NOTE

For more information about the column organizer read the G1383-90011 Column Organizer Quick Reference Guide

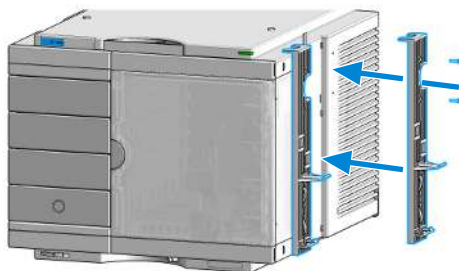
- 1 Attach the column stand mount to the column organizer.
- 2 Attach both clamps to one side of the valve drive.
- 3 Install the leak tray and leak sensor to the valve.
- 4 Slide the valve onto the column stand mount.
- 5 Attach the waste tubing to the leak plane and guide it to a proper waste container.
- 6 Connect the power cable to the power connector at the rear of the module.
- 7 Connect the CAN interface connection to one of the CAN ports at the rear of the module.
- 8 Power on the module by switching the Power button at the rear of the module if a valve head has been already installed. Otherwise continue with [G7116B_Installation of the Valve Heads](#) on page 48.

Mounting the Valve Drive as a Part of Online Sample Manager

Parts required	Qty.	p/n	Description
	2	 G3167-42000	Single Holder UVD Multi Function
	4	 0515-5869	Screw-Tapping Pan-HD Hexalobular-Recess
	1	 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm

Preparations

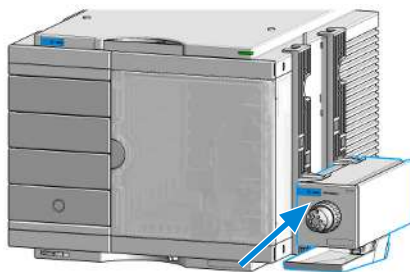
- Ensure the power switch at the rear of the module is off and the power connector is unplugged.
- 1 Identify the marks for the screws on the side panel of the Sampler and screw the clamps to the side panel of the module.



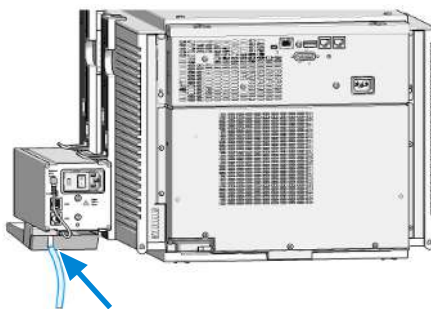
Installation

Installing the Valve Drive

- 2 Slide the valve drive into the clamp guides from front to back.



- 3 Attach the waste tubing to the leak plane and guide it to a proper waste container.



Mounting the Valve Drive to an Infinity III Module (except Multicolumn Thermostat)

Parts required**Qty.****p/n****Description**

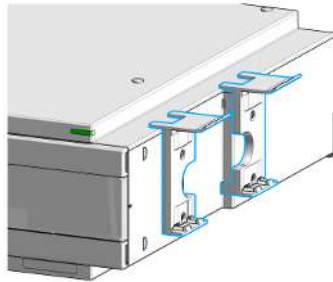
1



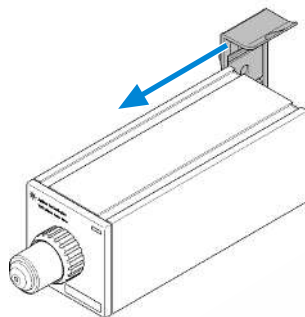
5067-5685

Infinity II & III Clamp Guide Kit

- 1 Ensure the power switch at the rear of the module is OFF and the power connector is unplugged.
- 2 Identify the marks for the screws on the side of the Infinity III module cabinet and screw the valve clamps to the side panel of the cabinet using the screws from the kit.



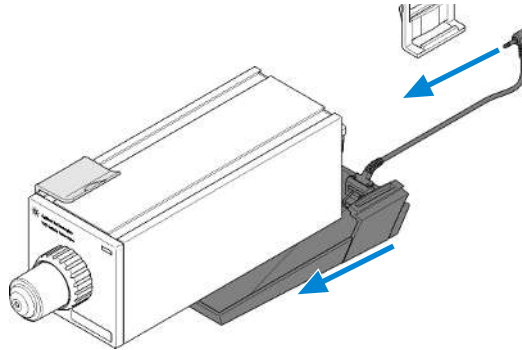
- 3 Slide the first clamp on to one side of the valve drive housing.



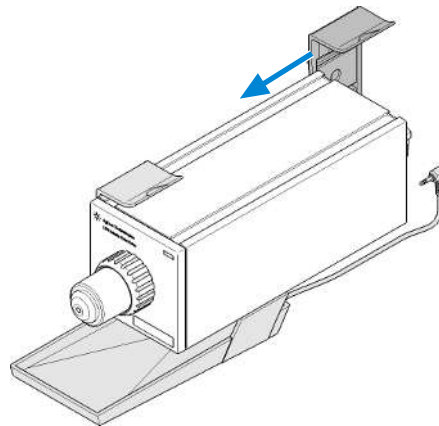
Installation

Installing the Valve Drive

- 4 Install the leak tray and leak sensor to the valve.



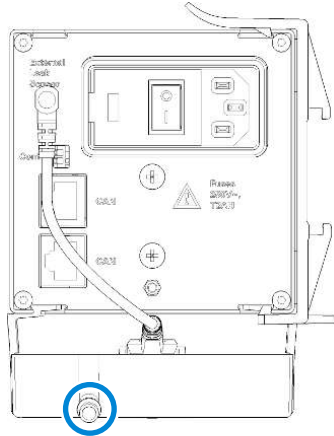
- 5 Slide the second clamp on to the same side as the first one.



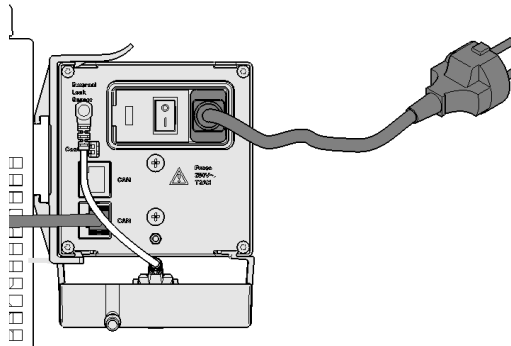
Installation

Installing the Valve Drive

- 6 Attach the waste tubing to the leak plane and guide it to a proper waste container.



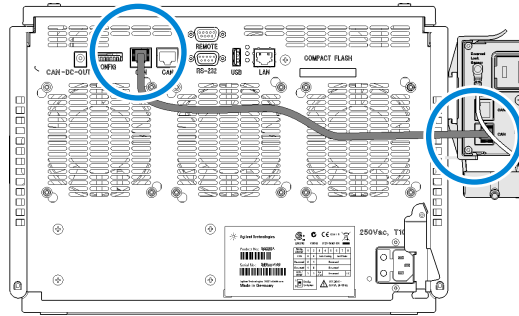
- 7 Connect the power cable to the power connector at the rear of the module.



Installation



Installing the Valve Drive

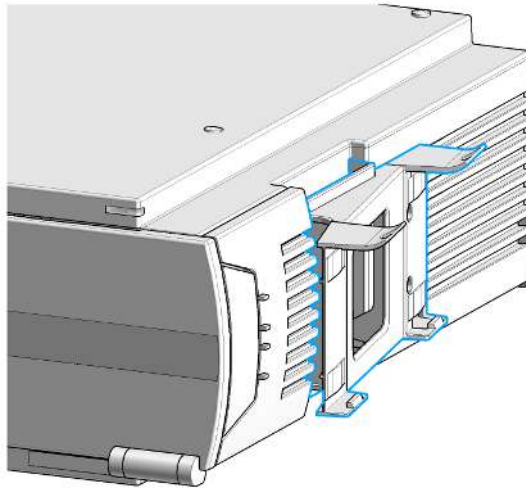
- 8 Connect the CAN interface connection.



- 9 Power on the module by switching the Power button at the rear of the module if a valve head has been already installed. Otherwise continue with [G7116B_Installation of the Valve Heads](#) on page 48.

Mounting the Valve Drive to a G7116A/B Multicolumn Thermostat

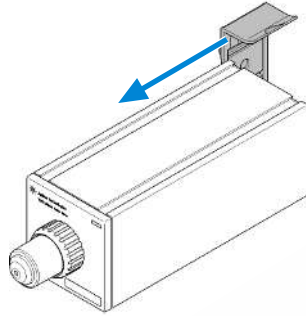
Parts required	Qty.	p/n	Description
	1	 5067-6138	Infinity II & III Valve Holder Kit Right , or
	1	 5067-6139	Infinity II & III Valve Holder Kit Left
1 Ensure the power switch at the rear of the module is OFF and the power connector is unplugged.			
2 Identify the marks for the screws on the side of the Infinity II/III MCT cabinet and screw the valve clamp to the side panel of the cabinet using the screws from the kit.			



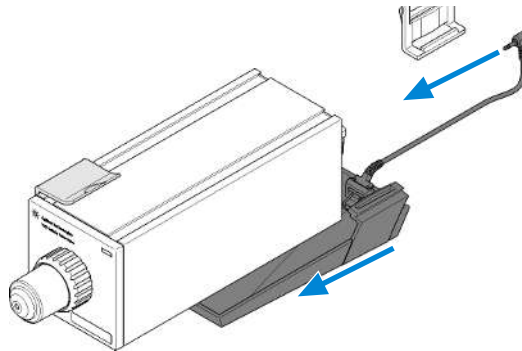
Installation

Installing the Valve Drive

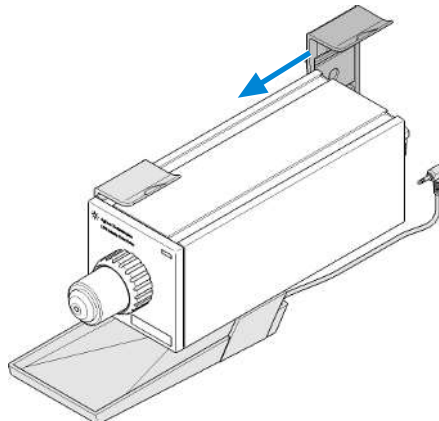
- 3 Slide the first clamp on to one side of the valve drive housing.



- 4 Install the leak tray and leak sensor to the valve.



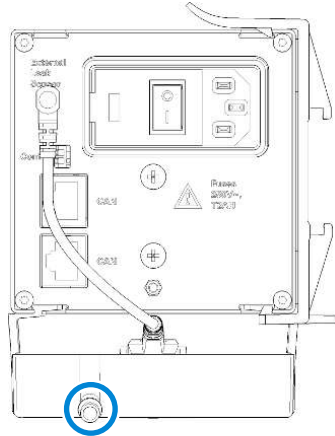
- 5 Slide the second clamp on to the same side as the first one.



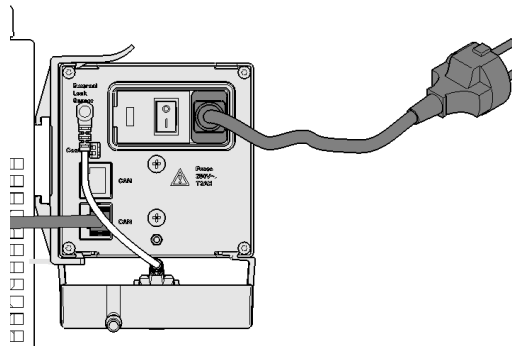
Installation

Installing the Valve Drive

- 6 Attach the waste tubing to the leak plane and guide it to a proper waste container.



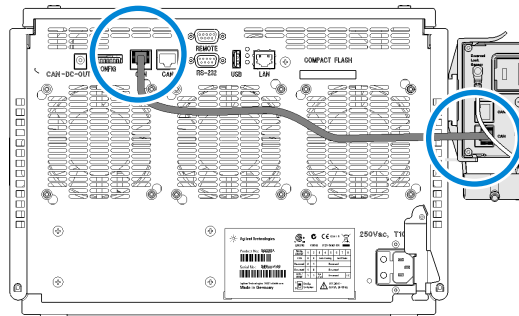
- 7 Connect the power cable to the power connector at the rear of the module.



Installation

Installing the Valve Drive

- 8 Connect the CAN interface connection.



- 9 Power on the module by switching the Power button at the rear of the module if a valve head has been already installed. Otherwise continue with [G7116B_Installation of the Valve Heads](#) on page 48.

Installing the Valve Heads

If ordered, the valve drives are factory-installed in the Multicolumn Thermostat. The valve heads are interchangeable and can be easily mounted.

At the first installation, the transportation lock and the dummy valve have to be removed, see [Remove the Transportation Lock and the Valve Dummy](#) on page 52. The valve heads can be installed by mounting the valve heads onto the valve drives and fastening the nut manually (do not use any tools).

Be sure that the guide pin snaps into the groove of the valve drive thread.

NOTE

The valves are mounted on pull-out rails to allow easy installation of capillaries. Push the valve gently into its housing until it snaps into the inner position, push it again and it slides out.

When all capillaries are installed, push the valve back into its housing, see [Install the Valve Head and Connect Capillaries](#) on page 48.

Install the Valve Head and Connect Capillaries



For bio-inert modules use bio-inert parts only!
Do not mix with bio / biocompatible parts.



For biocompatible modules use bio / biocompatible parts only!
Do not mix with bio-inert parts.

CAUTION

The valve actuator contains sensitive optical parts, which need to be protected from dust and other pollution. Pollution of these parts can impair the accurate selection of valve ports and therefore bias measurement results.

- Always install a valve head for operation and storage. For protecting the actuator, a dummy valve head can be used instead of a functional valve. Do not touch parts inside the actuator.

CAUTION

Column Damage or Bias Measurement Results

Switching the valve to a wrong position can damage the column or bias measurement results.

- Fit the lobe to the groove to make sure the valve is switched to the correct position.

CAUTION

Valve Damage

Using a low pressure valve on the high pressure side can damage the valve.

- When using multiple column compartments as part of a method development solution, make sure that the high pressure valve head is connected to the autosampler and the low pressure valve head is connected to the detector.

NOTE

For information about the compatibility mode of 800 bar valve heads see 01200-90134 (Information on RFID Tag Technical Note) .

NOTE

For a correct installation of the valve head, the outside pin (red) must completely fit into the outside groove on the valve drive's shaft (red). A correct installation is only possible if the two pins (green and blue) on the valve head fit into their corresponding grooves on the valve drive's actuator axis. Their match depends on the diameter of the pin and groove.

NOTE

The tag reader reads the valve head properties from the valve head RFID tag during initialization of the module. Valve properties will not be updated, if the valve head is replaced while the module is on. Selection of valve port positions can fail, if the instrument does not know the properties of the installed valve.

NOTE

To allow correct valve identification, power off the valve drive for at least 10 s.

NOTE

For firmware requirements see 01200-90134 (Information on RFID Tag Technical Note) which is included to each valve head.

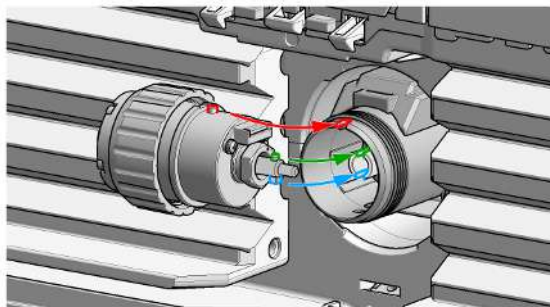
CAUTION

Sample degradation and contamination of the instrument

Metal parts in the flow path can interact with the bio-molecules in the sample leading to sample degradation and contamination.

- For bio applications, always use dedicated bio parts, which can be identified by the bio-inert symbol or other markers described in this manual.
- Do not mix bio, and non-bio modules or parts in a bio system.

- 1 Insert the valve head into the valve shaft.

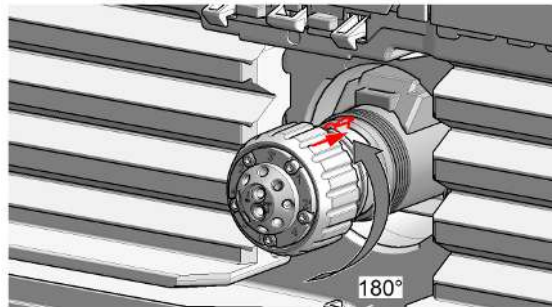
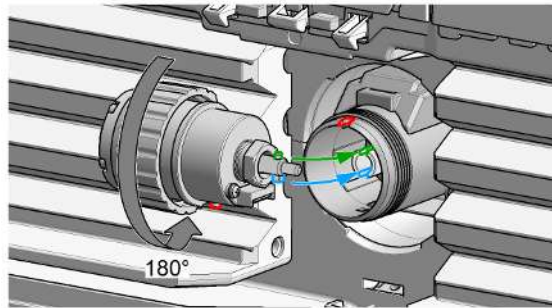


Installation

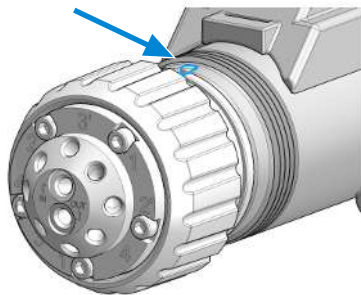
Installing the Valve Heads

OR

If the outside pin does not fit into the outside groove, you have to turn the valve head until you feel that the two pins snap into the grooves. Now you should feel additional resistance from the valve drive while continuously turning the valve head until the pin fits into the groove.



- 2 When the outer pin is locked into the groove, manually screw the nut onto the valve head.



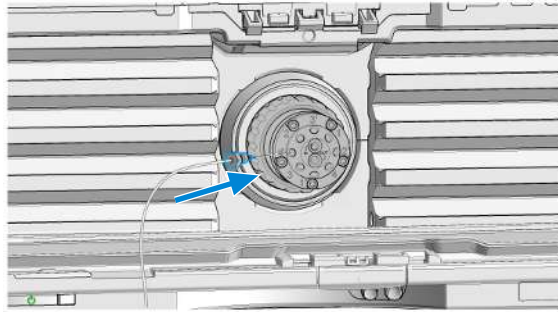
NOTE

Fasten the nut manually. Do not use any tools.

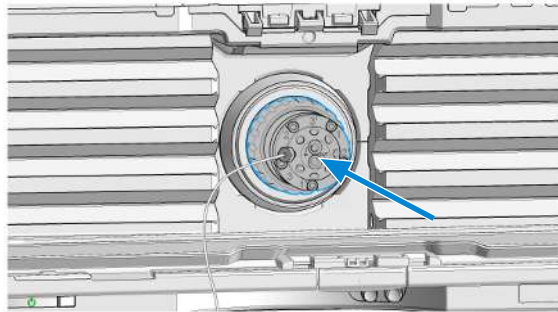
Installation

Installing the Valve Heads

- 3 Install all required capillary connections to the valve.



- 4 Push the valve head until it snaps in and stays in the rear position.



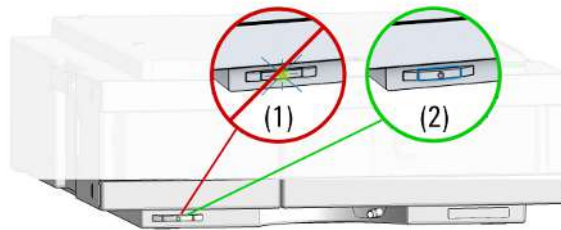
- 5 Power on or power-cycle your module, so the valve head gets recognized during module initialization.

Remove the Transportation Lock and the Valve Dummy

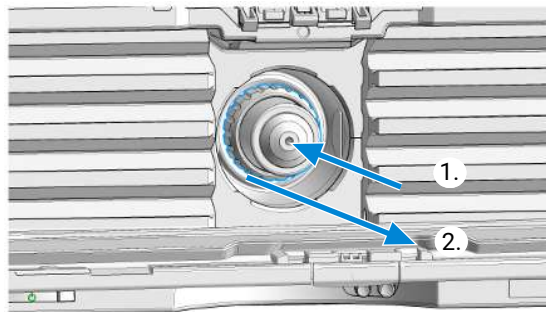
The following procedure demonstrates the necessary steps for installing the valve head to the valve drive of a Multicolumn Thermostat (MCT).

For the installation of a valve head to a G1170A Valve Drive you can ignore the steps that describe the MCT features of the transportation lock and spring loaded valve drive.

- 1 Switch off the module.

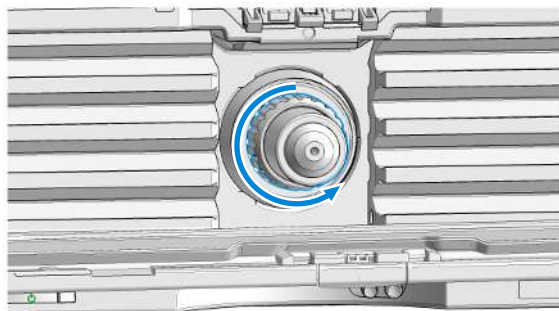


- 2 When unscrewing the transportation lock, push it back until the last screw is removed - the valve rail is spring-loaded.
- 3 Press on the valve dummy (1.) to release it (2.) (spring-loaded valve rail).

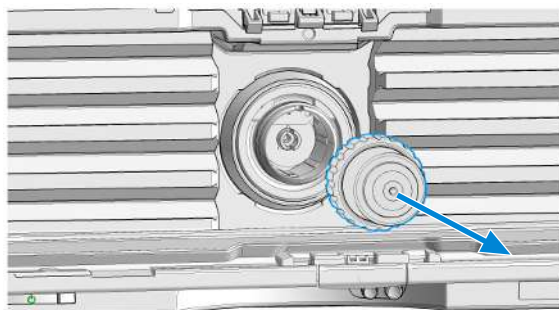


Installation**Installing the Valve Heads**

- 4 Unscrew the valve dummy.



- 5 Remove the valve dummy from the valve drive.



Handling Leak and Waste

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

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

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

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

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

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

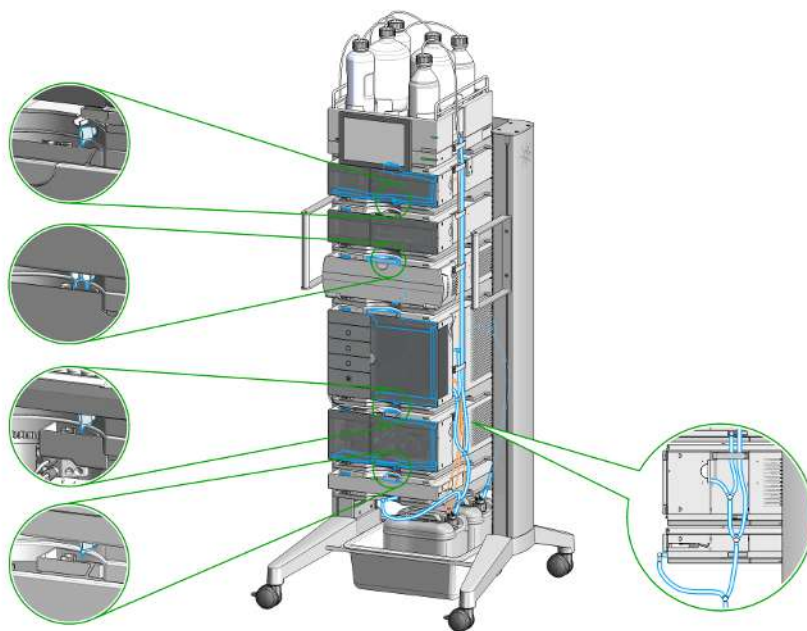


Figure 7: Infinity III Leak Waste Concept (Flex Bench installation)

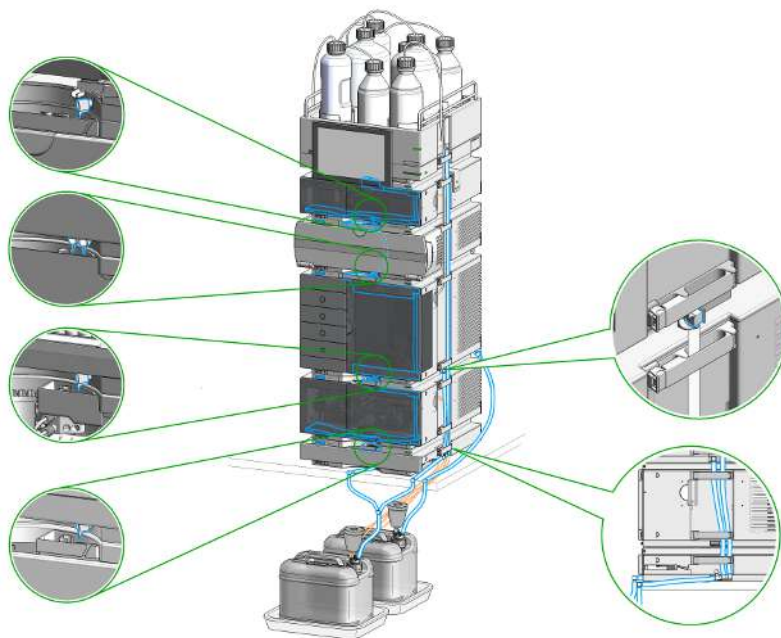


Figure 8: Infinity III Single Stack Leak Waste Concept (bench installation)

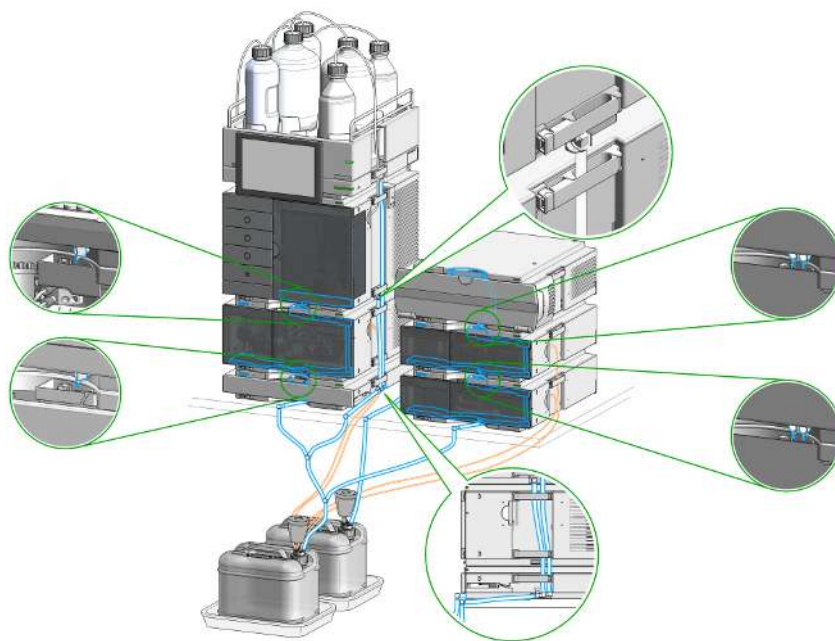


Figure 9: Infinity III Two Stack Leak Waste Concept (bench installation)

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

Drain Connectors Installation

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

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

For illustration, see [Handling Leak and Waste](#) on page 55.

Parts required

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

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

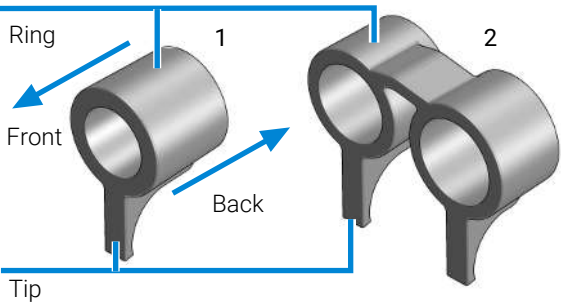


Figure 10: Overview of Drain Connectors: Single (left) and Double (right)



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

Table 5: Compatibility of drain connectors and modules

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

Preparations

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

NOTE

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

NOTE

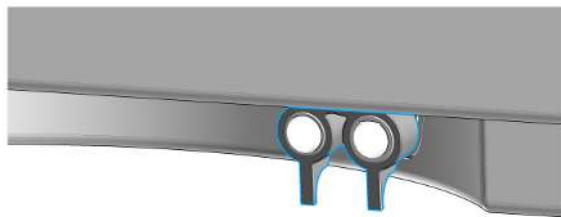
In case of incorrect installation, drain connectors cannot fully perform the intended function.

NOTE

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

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

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

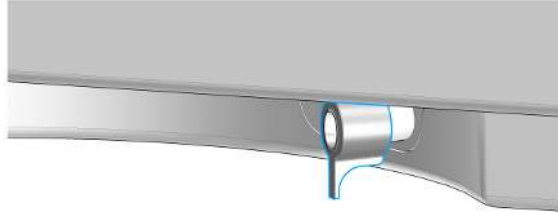


Install Single Drain Connectors on other modules in the LC stack

Installation

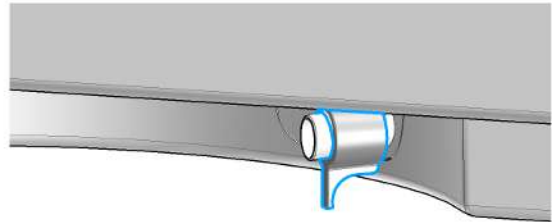
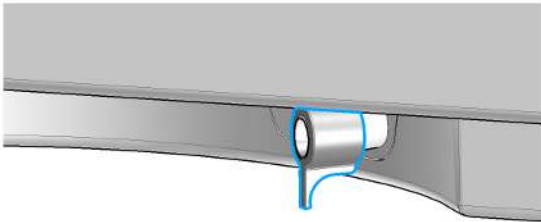
Handling Leak and Waste

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



Make sure that the following requirements are covered:

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



Waste Concept

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



Waste Guidance



NOTE

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

Leak Sensor

CAUTION

Solvent incompatibility

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

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



4 Using the Module

This chapter provides information on how to use the module.

General Information 66

Turn On/Off 66

Status Indicators 68

Software Configuration 70

Configuring the Valves in the Software 70

Assign a Pressure Limit Cluster to the Valve 70

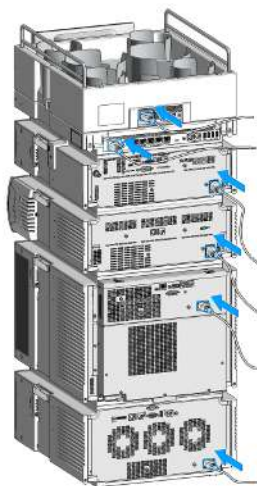
Possible Method Settings for the Valve 72

General Information

Turn On/Off

This procedure exemplarily shows an arbitrary LC stack configuration.

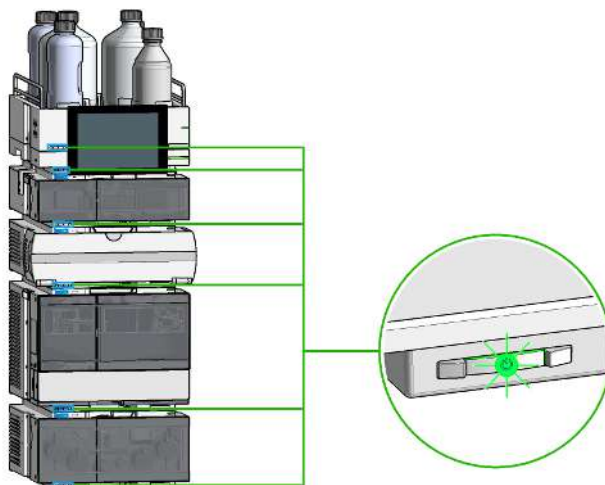
1



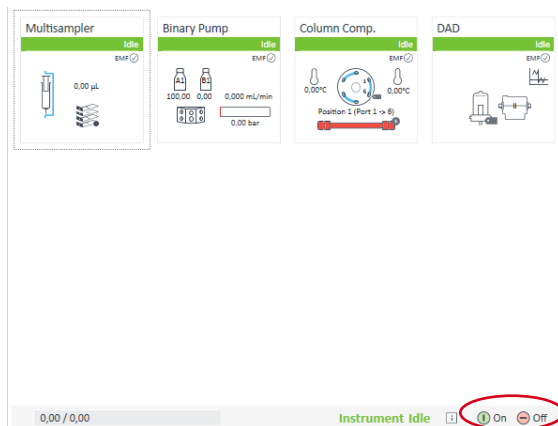
Using the Module

General Information

2 On/Off switch: On



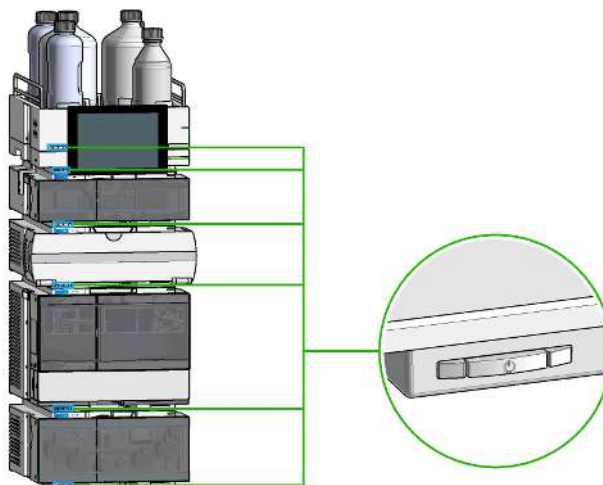
3 Turn instrument On/Off with the control software.



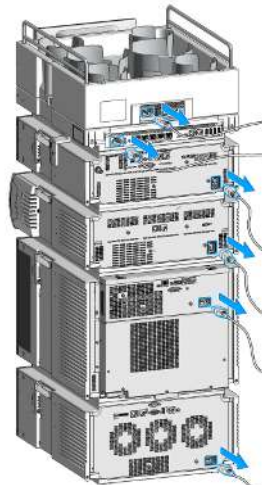
Using the Module

General Information

4 On/Off switch: Off



5



Status Indicators

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

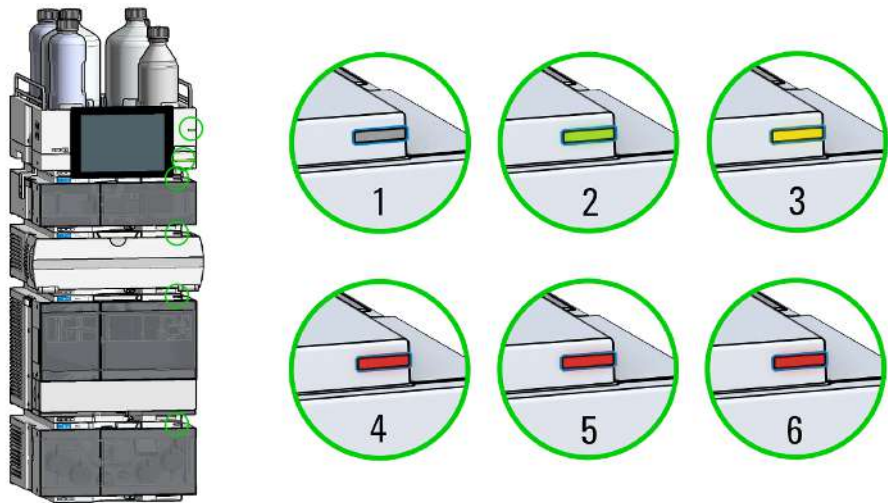


Figure 11: Arbitrary LC stack configuration (example)

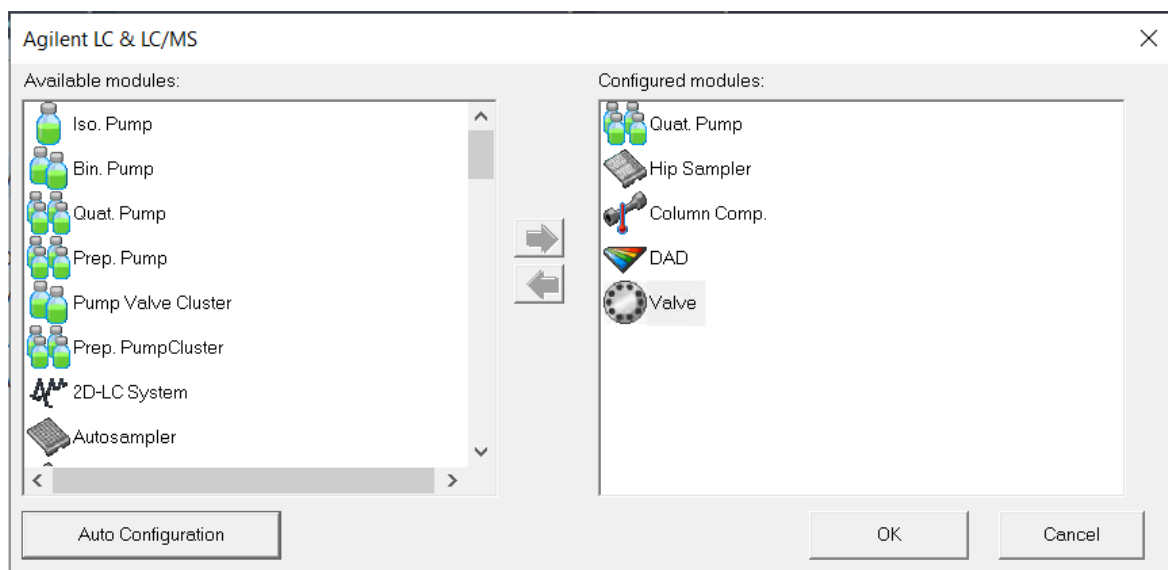
1	Idle
2	Run mode
3	Not-ready. Waiting for a specific pre-run condition to be reached or completed.
4	Error mode - interrupts the analysis and requires attention (for example, a leak or defective internal components).
5	Resident mode (blinking) - for example, during update of main firmware.
6	Bootloader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

InfinityLab Assist Hub Status Indicator

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

Software Configuration

Configuring the Valves in the Software

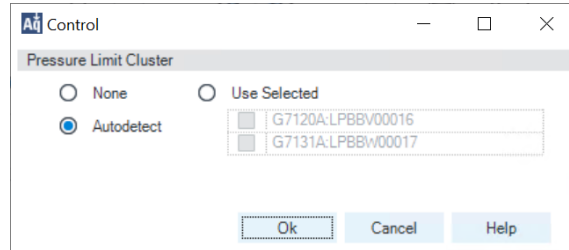


- 1 Configure your instrument by choosing between available modules.
- 2 Choose your valve drive(s) from the list of 'Configurable modules'(auto or manual configuration).

NOTE

A valve head needs to be detected from the module before you can configure it in the software. Please refer to [G7116B_Installation of the Valve Heads](#) on page 48 for details.

Assign a Pressure Limit Cluster to the Valve



The valve head's RFID Tag holds the information about its maximum pressure. An assigned pressure cluster partner assures that this limit is not exceeded. The maximum pressure information of the valve head will become the new maximum pressure setting of the assigned pressure cluster partner. This helps to avoid damage to the valve head if the maximum pressure of the pump is higher than the valve head's maximum pressure.

- 1 Open the valve's control menu (right click on the valve's GUI).
- 2 Assign the valve's pressure partner ('None' or several partners possible).

NOTE

Assigning a Pressure Limit Cluster only works in combination with a 1290 Infinity and 1290 Infinity II/III pump.

Possible Method Settings for the Valve

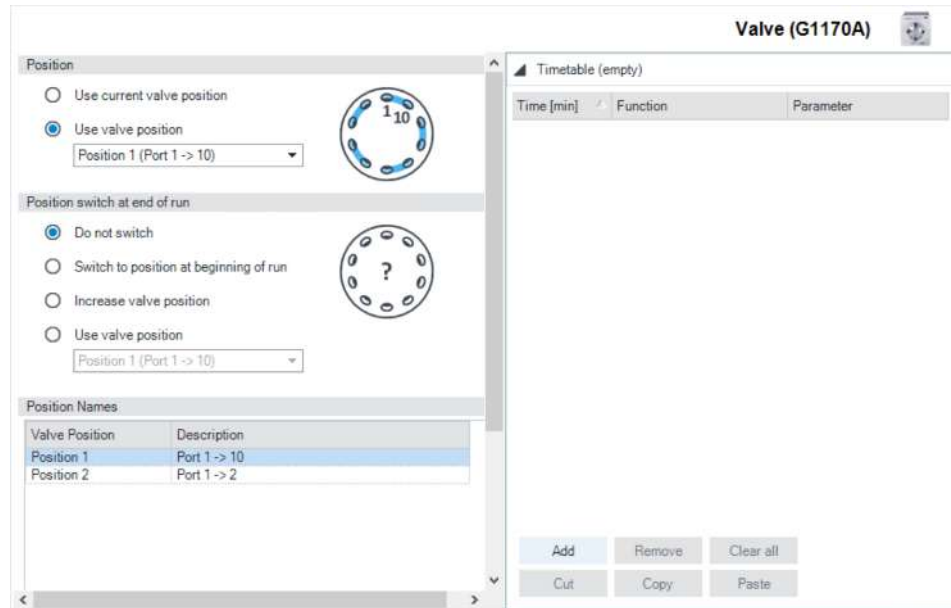


Figure 12: Method settings of a valve installed in the External Valve Drive

After the Valve Drives and Quick Change Valve Heads have been installed and configured, the valve parameters can be edited in the software. Choose a method by right clicking on the valve's GUI icon to open the valve method dialog box as displayed [Figure 12](#) on page 72.

Position

When loading a method the valve is switched to the position that is displayed in the Position dropdown box. If you select Use Current, the valve remains in the current position, when a new method is loaded.

If you select a new position in the dropdown box and click OK, the Setup Valve dialog box will be closed and the valve switches to the new position.

Position switch at end of run

Gives you the possibility to define a position the valve is switched to at the end of the run.

Position Names

Define the Position Names that is used for the method report and the instrument actuals. The Position Description is limited to 19 characters.

Time Table

The Time Table can be used to edit and run a valve program during a sequence of sample runs. The Time Table contains 3 columns (Line, Time and Position). The Time Table is limited to 20 lines. The number of selection in the Position column depends on the valve that has been configured.

Position X	Switches the valve to the selected position.
Next Position	Switches to the next available position. If the valve is on the highest position it will switch to position 1.
Insert	Inserts a line in the Time Table above the selected position.
Append	Appends a line at the end of the Time Table.
Cut	Cuts the selected line(s) out of the Time Table and saves it to the clipboard.
Copy	Copies the selected line(s) from the table to the clipboard.
Paste	Pastes line(s) from the Clipboard to the Time Table.



5

Diagnostics and Troubleshooting

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

Diagnostic Features 75

User Interfaces 75

Troubleshooting With HPLC Advisor 75

Overview of the Module's Indicators and Test Functions 76

Status Indicators 76

Error Messages 76

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.
- Screenshots used within these procedures are based on the Agilent Lab Advisor Software.

Troubleshooting With HPLC Advisor

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

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

Overview of 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).



6 Error Information

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 78

General Error Messages 79

- Timeout 79
- Shutdown 79
- Remote Timeout 80
- Lost CAN Partner 81
- Leak 82
- Leak Sensor Open 83
- Leak Sensor Short 84
- Compensation Sensor Open 85
- Compensation Sensor Short 86

Module Specific Error Messages 88

- Initialization of Valve Failed 88
- Valve Switching Failed 88
- Valve Tag Violation 89
- Pressure Cluster Partner Missing 90
- Position Cluster Partner Missing 91

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 Service Manual for the degasser or the pump that has the degasser built-in.

Remote Timeout

Error ID: 70

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

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

Lost CAN Partner

Error ID: 71

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

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

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

Leak

Error ID: 64

A leak was detected in the module.

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

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

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 mainboard.	• 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.

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.

Compensation Sensor Open

Error ID: 81

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

The resistance across the temperature compensation sensor (NTC) on the mainboard 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 mainboard.	<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 mainboard in the module has failed (open circuit).

The resistance across the temperature compensation sensor (NTC) on the mainboard 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 mainboard.	<ul style="list-style-type: none">• Please contact your Agilent service representative.

Module Specific Error Messages

Initialization of Valve Failed

Error ID: 24000

During the initialization process the motor of the valve drive moves to some special positions depending on the installed valve head. A failure in this process means either that the movement couldn't be performed properly or it was not noticed correctly by the sensor.

Probable cause		Suggested actions
1	Mechanical problems. Friction too high or blockages on the valve drive's motor or on the valve head.	<ul style="list-style-type: none">• Check valve head for correct installation.• Try to identify the source of trouble by installing a different valve head if possible.• Please contact your Agilent service representative.
2	Defect Sensor on the Valve Drive Motor.	<ul style="list-style-type: none">• Check valve head for correct installation.• Try to identify the source of trouble by installing a different valve head if possible.• Please contact your Agilent service representative.

Valve Switching Failed

Error ID: 24001

The valve drive was not able to operate the valve head correctly. Either due to mechanical reasons or the movement couldn't be detected correctly.

Probable cause		Suggested actions
1	Mechanical problems. Friction too high or blockages on the valve drive's motor or on the valve head.	<ul style="list-style-type: none">• Check valve head for correct installation.• Try to identify the source of trouble by installing a different valve head if possible.• Please contact your Agilent service representative.
2	Defect Sensor on the Valve Drive Motor.	<ul style="list-style-type: none">• Check valve head for correct installation.• Try to identify the source of trouble by installing a different valve head if possible.• Please contact your Agilent service representative.

Valve Tag Violation

Error ID: 24006

The valve drive identified a different valve head than it had identified during the last initialization.

NOTE

Soft power-down power supply of the valve drive.
Whenever you want to power cycle the valve drive for a re-boot, it needs to be powered off for at least 10 seconds.

Probable cause		Suggested actions
1	A valve head has been exchanged (hot-plugged) while the valve drive was still powered on.	<ul style="list-style-type: none">Change the valve head. It is important to have the valve switched off for at least 10 s after or before a new valve head has been installed.

Pressure Cluster Partner Missing

Error ID: 2523

The connection from the valve drive to a defined pressure cluster partner is lost.

Probable cause		Suggested actions
1	Communication issues.	<ul style="list-style-type: none">• Check the CAN cable connections of the modules.
2	Configuration mismatch.	<ul style="list-style-type: none">• Check and correct if necessary the valve configuration and presence of defined pressure cluster partner.

Position Cluster Partner Missing

Error ID: 4526

Probable cause		Suggested actions
1	Communication issues.	<ul style="list-style-type: none">• Check the CAN cable connections of the modules.
2	Configuration mismatch.	<ul style="list-style-type: none">• Check and correct if necessary the valve configuration and presence of defined position cluster partner.• If the module was moved to another LC stack, perform Firmware Declustering in Service & Diagnostic section of Lab Advisor.

This chapter gives you an overview and instructions about the possible maintenance and repair procedures that can be performed by the user.

Safety Information Related to Maintenance 94

Introduction to Maintenance 96

Overview of Maintenance 97

Cleaning the Module 98

Replacing Parts of the Valve Head 99

Replacing the Fuses of the Valve Drive 101

Replace Valve Heads of the Valve Drive 105

InfinityLab LC Method Development Solutions User Guide 000

Replace the Module Firmware 110

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.

CAUTION

Sample degradation and contamination of the instrument

Metal parts in the flow path can interact with the bio-molecules in the sample leading to sample degradation and contamination.

- For bio applications, always use dedicated bio parts, which can be identified by the bio-inert symbol or other markers described in this manual.
- Do not mix bio, and non-bio modules or parts in a bio system.

Introduction to Maintenance

The module is designed for easy maintenance. The most frequent maintenance such as maintaining valve heads can be done from the front with the modules in place in the system stack.

NOTE

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

Overview of Maintenance

The following pages describe maintenance procedures (simple repairs) that can be done without opening the main cover.

Table 6: Maintenance Procedures

Procedure	Typical Frequency	Notes
Cleaning the Module on page 98	If required	
Replacing Parts of the Valve Head on page 99	If the valve performance shows indication of leakage or wear	
Replacing the Fuses of the Valve Drive on page 101	When a fuse is defect	
Replace Valve Heads of the Valve Drive on page 105	If the valve performance shows indication of leakage or wear	
Replace the Module Firmware on page 110	If required	

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.

Replacing Parts of the Valve Head

For details about the needed parts and orientation please refer to [Valve Drive Parts](#) on page 112.

Disassembling and reassembling the valve head

For bio-inert modules use bio-inert parts only!
Do not mix with bio / biocompatible parts.



For biocompatible modules use bio / biocompatible parts only!
Do not mix with bio-inert parts.



When

- Stator head: Scratches and damage on the inner surface, blockages
- Stator face assy: When visibly scratched, or when the valve performance shows indication of leakage or wear
- Rotor seal assy: When the valve performance shows indication of leakage or wear

Tools required

Qty.	p/n	Description
1		Hex key

Parts required

Qty.	p/n	Description
1		Quick Change Valve Head

WARNING

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

- Be sure that no solvent can drop out of the solvent connections when removing them from your valve head.
- 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.

- 1 Use the Hex Key to open and remove the Stator Screws from the Stator Head.

- 2 Carefully disassemble the necessary Valve Head parts to gain access to the one you want to replace. While doing so please observe the orientation of the parts.
- 3 Independent of the part you want to replace always inspect all parts for signs of damage.
- 4 Replace the proposed part.

NOTE

Always mind the correct orientation of the parts and avoid to touch their surfaces.


- 5 Turn each of the screws an equal amount until they are finger-tight, then tighten them for another half turn.

Replacing the Fuses of the Valve Drive

When

- If the valve drive shows no reaction.

Tools required

Qty.	p/n	Description
	 5023-3138	Reversible Screwdriver + Blade 1,0 x 5,5

Parts required

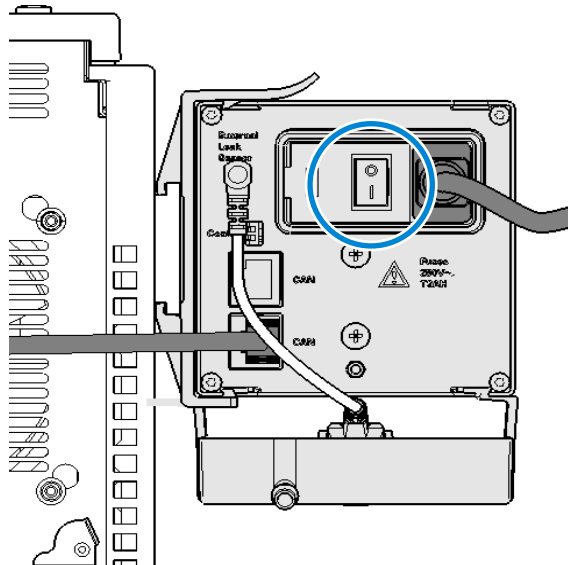
Qty.	p/n	Description
2	 2110-1486	Fuse 2 AT250 V

WARNING

Electrical shock

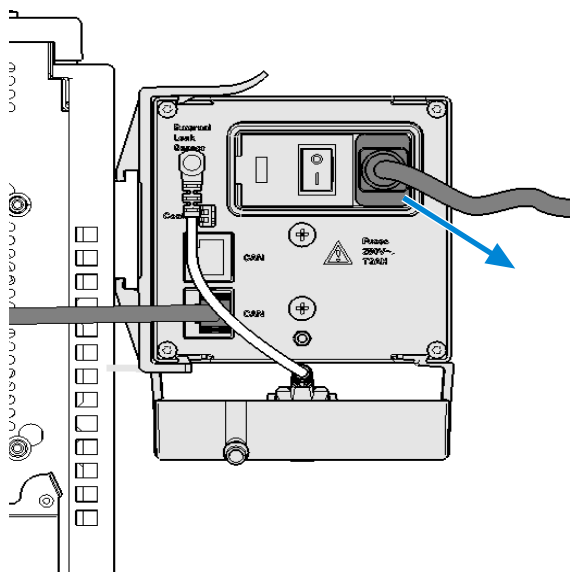
- Disconnect the valve drive from line power before changing a fuse or trying to open the hatch of the power input socket.
- Never re-connect the line power before having the power input socket closed.

- 1 Switch off the instrument. The power switch is located at the rear of the valve drive.

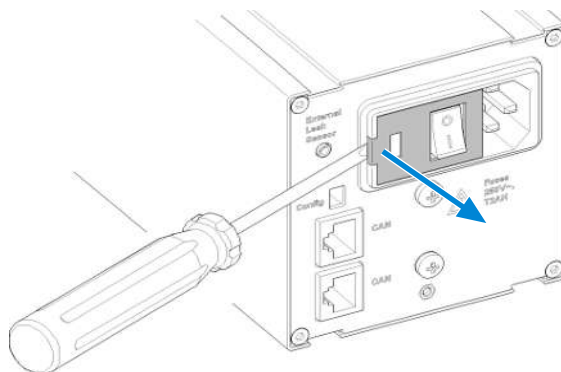


Replacing the Fuses of the Valve Drive

- 2 Disconnect the power cable from the power input socket.

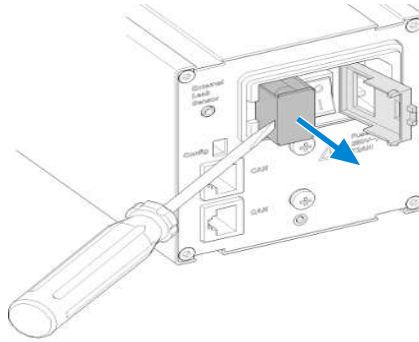


- 3 To access the fuse drawer, gently lift the outer plastic housing of the power inlet socket using a flat screwdriver.



Maintenance**Replacing the Fuses of the Valve Drive**

- 4 Pull out the fuse drawer as shown.

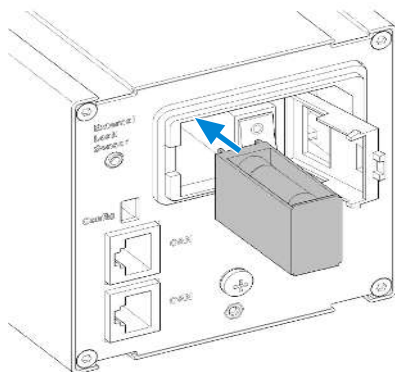


- 5 Replace the defect fuse(s).

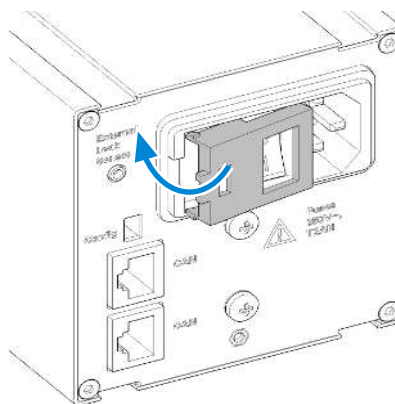
Maintenance

Replacing the Fuses of the Valve Drive

- 6 Slide in the fuse drawer and push till it fits tightly.



- 7 Finally, close the fuse drawer housing, reconnect the instrument to the power line and switch it on.



Replace Valve Heads of the Valve Drive

Several optional valve heads are available, which can be installed and exchanged easily.



For bio-inert modules use bio-inert parts only!

Do not mix with bio / biocompatible parts.



For biocompatible modules use bio / biocompatible parts only!

Do not mix with bio-inert parts.

Parts required	Qty.	p/n	Description
	1		Agilent Quick Change Valve Head For details, see Valve Options Overview on page 115

WARNING

Toxic, flammable and hazardous solvents, samples and reagents

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

- Be sure that no solvent can drop out of the solvent connections when removing them from your valve head.
- 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.

CAUTION

Valve Damage

Using a low pressure valve on the high pressure side can damage the valve.

- When using multiple column compartments as part of a method development solution, make sure that the high pressure valve head is connected to the autosampler and the low pressure valve head is connected to the detector.

CAUTION**Column Damage or Bias Measurement Results**

Switching the valve to a wrong position can damage the column or bias measurement results.

- Fit the lobe to the groove to make sure the valve is switched to the correct position.

CAUTION

The valve actuator contains sensitive optical parts, which need to be protected from dust and other pollution. Pollution of these parts can impair the accurate selection of valve ports and therefore bias measurement results.

- Always install a valve head for operation and storage. For protecting the actuator, a dummy valve head (part of G1316-67001 (Transportation Lock Kit)) can be used instead of a functional valve. Do not touch parts inside the actuator.

NOTE

For details, please refer to the *InfinityLab LC Method Development Solutions User Guide (InfinityLab-Method-Development-Solution-UseMa-en-SD-29000211.pdf, SD-29000211)*.

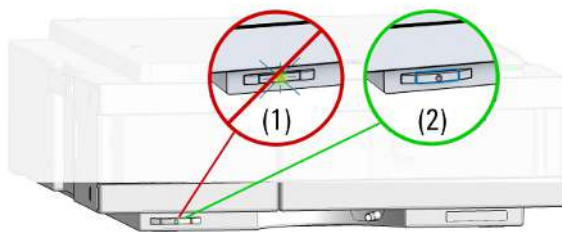
NOTE

The tag reader reads the valve head properties from the valve head RFID tag during initialization of the module. Valve properties will not be updated, if the valve head is replaced while the module is on. Selection of valve port positions can fail, if the instrument does not know the properties of the installed valve.

NOTE

To have the valve correctly recognized by the valve drive you must have the valve drive powered off for at least 10 seconds.

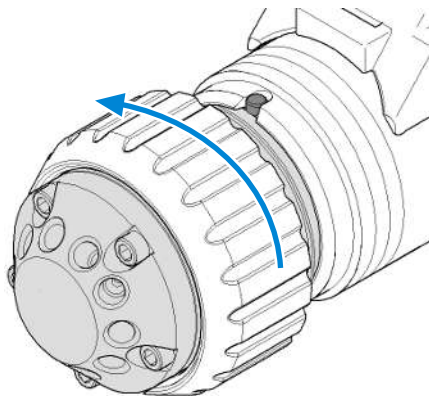
- 1 Switch off the module.



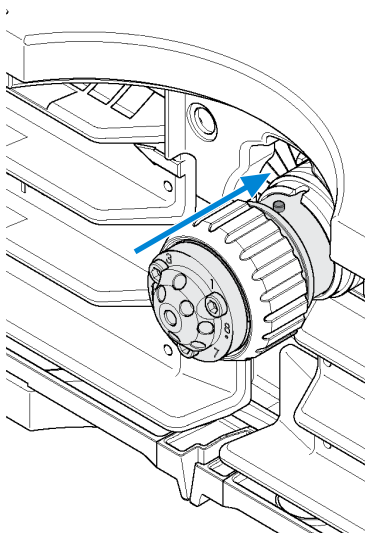
Maintenance

Replace Valve Heads of the Valve Drive

- 2 Push the valve head for bringing it to its outer position (column compartment only).
- 3 Remove all capillary connections from the valve head.
- 4 Unscrew the nut and remove the valve head.

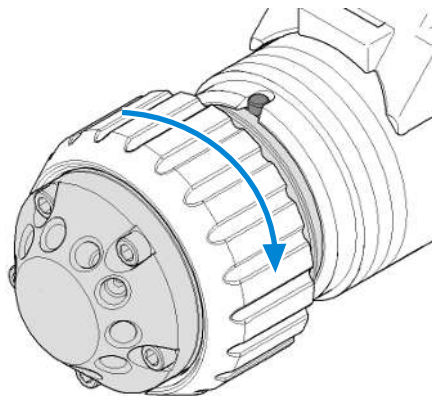


- 5 Put the new valve head onto the valve drive such that the lobe fits to the groove (see also [G7116B_Installation of the Valve Heads](#) on page 48).

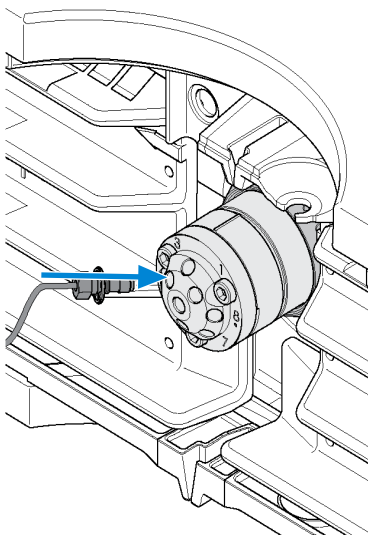


Maintenance**Replace Valve Heads of the Valve Drive**

- 6 Fasten the valve head onto the valve drive using the union nut (see also [G7116B_Installation of the Valve Heads](#) on page 48).

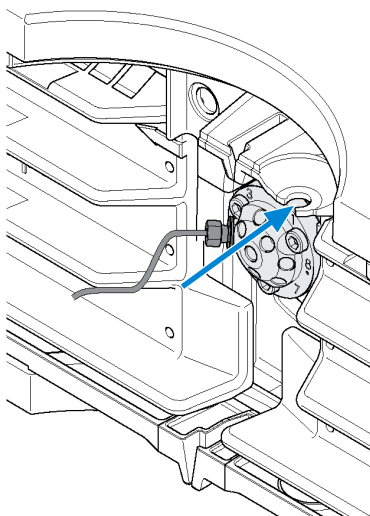


- 7 Install all required capillary connections to the valve head.

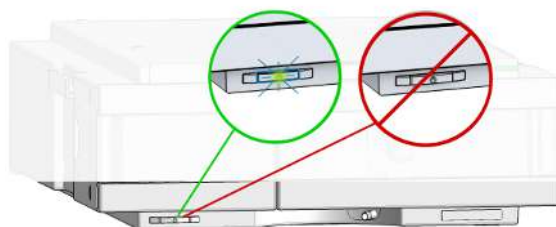


Maintenance**Replace Valve Heads of the Valve Drive**

- 8 Push the valve head until it snaps in and stays in the rear position (column compartment only).



- 9 Switch on the module.



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.



8 Parts and Materials for Maintenance

This chapter provides information on parts for maintenance.

Valve Drive Parts 112

1290 Infinity Valve Drive Parts 112

1290 Infinity II/III Valve Drive Parts 113

Valve Options Overview 115

Accessory Kit 117

Valve Drive Parts

1290 Infinity Valve Drive Parts

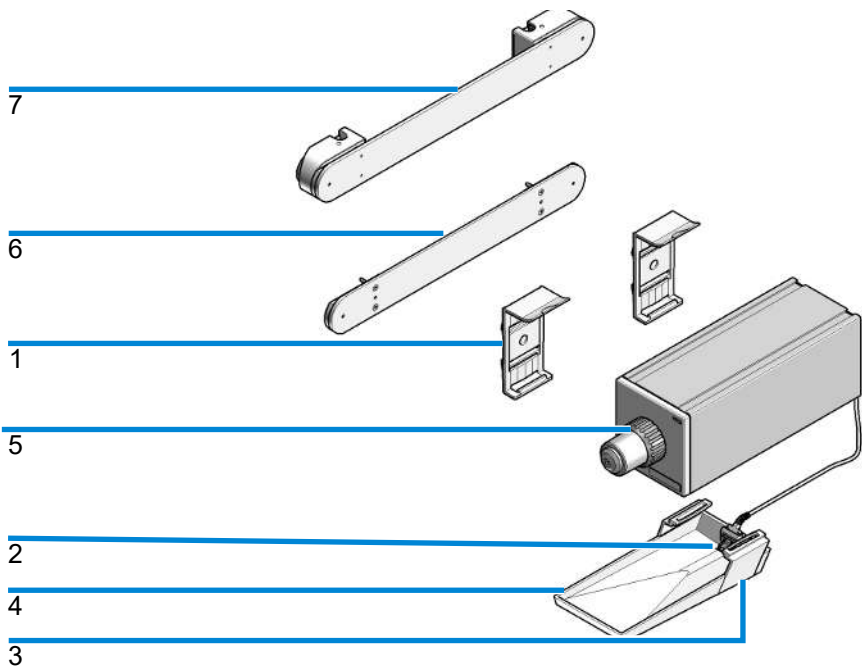










Figure 13: Parts for 1290 Infinity Valve Drive

#	p/n	Description
1	 5043-0275	Clamp guide For attaching the valve to a rail assembly
2	 5067-4792	Leak sensor assembly External leak sensor
3	 5043-0271	Holder leak plane
4	 5043-0270	Leak plane
5	 5068-0106	Spanner nut
	 2110-1486	Fuse 2 AT250 V

#	p/n	Description
6	 5067-4634	Valve Rail Kit
7	 5067-1510	Rail assy for column organizer

1290 Infinity II/III Valve Drive Parts

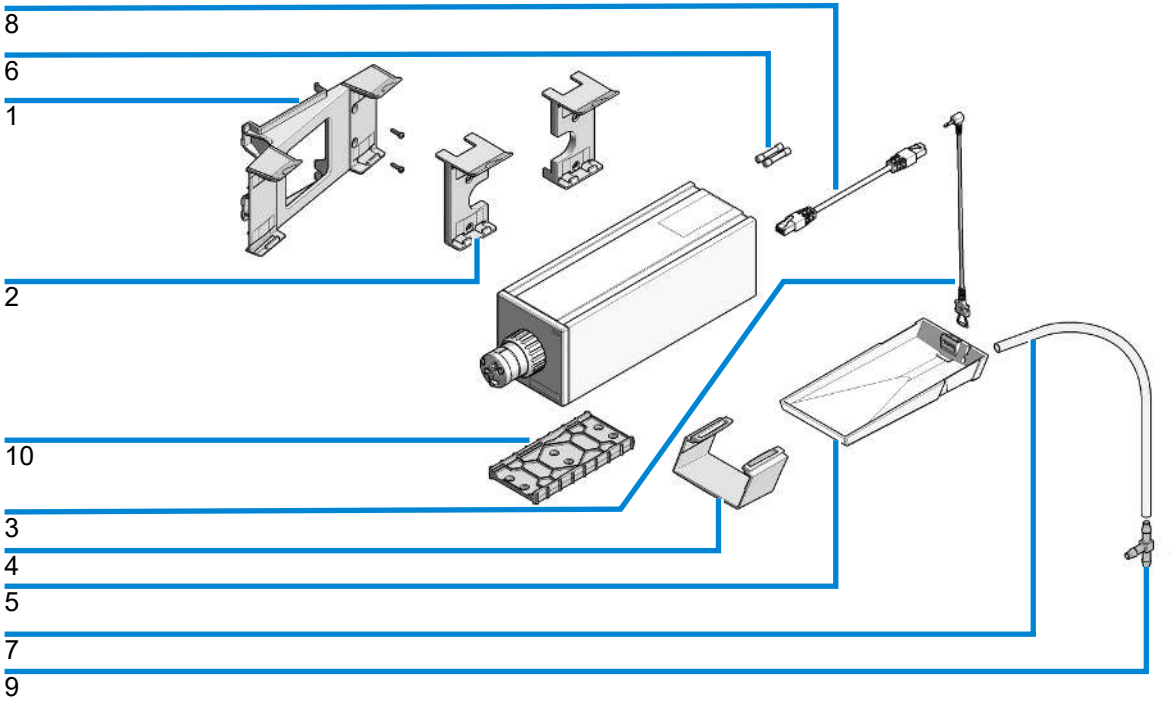













Figure 14: Parts for 1290 Infinity II/III Valve Drive

#	p/n	Description
1	 5067-6138	Infinity II & III Valve Holder Kit Right For G7116A/B
	 5067-6139	Infinity II & III Valve Holder Kit Left For G7116A/B (Not shown)
2	 5067-5685	Infinity II & III Clamp Guide Kit

#	p/n	Description
3	 5067-4792	Leak sensor assembly External leak plane
4	 5043-0271	Holder leak plane
5	 5043-0270	Leak plane
6	 2110-1486	Fuse 2 AT250 V
7	 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm
8	 5181-1519	CAN cable, Agilent module to module, 1 m
9	 5500-1156	T-Tube Connector ID6.4
10	 5043-0269	Adapter-profile For G1170A (Multiple valve drives can be connected with adapter profiles)

Valve Options Overview

This overview gives a summary of the main parts and assemblies. More details are available with each valve option in this chapter.










Table 7: Replacement parts standard valve heads

Valve Head	Rotor Seal	Stator Head	Stator Screws	Stator Ring	Stator Face	Other
5067-4233 8-Position/18-Port Valve 1300 bar	5068-0200 (P EEK)	5068-0199	5068-0089	n.a.		
5067-4241 2-Position/6-Port Valve 1300 bar	5068-0207 (P EEK)	5068-0006	1535-4857	5068-0120		
5067-4240 2-Position/10-Port Valve 1300 bar	5068-0205 (P EEK)	5068-0011	5068-0019	n.a.		
5067-4273 6-Position/14-Port Valve 1300 bar	5068-0242 (P EEK)	5068-0241	5068-0089	n.a.		
5067-4284 6-Position/14-Port Valve 800 bar	5068-0298 (P EEK)	5068-0241	5068-0089	n.a.		
5067-6682 2-Position/10-Port Valve Bio 1300 bar	5068-0205 (P EEK)	5068-0286	5068-0019	n.a.		Bearing ring: 1535-4045
5067-4279 4-Position/10-Port Valve 800 bar	5068-0264 (PEEK)	5068-0263	5068-0019	n.a.	n.a.	Bearing ring: 1535-4045
5067-4282 2-Position/6-Port Valve 800 bar	0101-1409 (PEEK)	0101-1417	1535-4857	5068-0120	n.a.	Bearing ring: 1535-4045

Valve Head	Rotor Seal	Stator Head	Stator Screws	Stator Ring	Stator Face	Other
5067-4148 2-Position/6-Port Bio-Inert Valve 600 bar	0101-1409 (PEEK)	5068-0060	5068-0020 (10/pack)	n.a.	0100-1851 (ceramic)	Bearing ring: 1535-4045
5067-4132 2-Position/10-Port Bio-Inert Valve 600 bar	5068-0041 (PEEK)	5068-0040	5068-0059	n.a.	5068-0095	Bearing ring: 1535-4045
5067-4283 2-Position/10-Port Valve 800 bar	0101-1415 (PEEK)	5068-0165	5068-0019	n.a.	n.a.	Bearing ring: 1535-4045
5067-4134 4-Position/10-Port Bio-Inert Valve 600 bar	5068-0045 (PEEK)	5068-0044	5068-0059	n.a.	5068-0093	Bearing ring: 1535-4045

Accessory Kit

G1170-68705 (Accessory Kit)

Qty.	p/n	Description
1	 5043-0270	Leak plane
1	 5043-0271	Holder leak plane
1	 5067-4792	Leak sensor assembly
1	 5181-1519	CAN cable, Agilent module to module, 1 m
2	 5043-0275	Clamp guide (for Infinity modules)
1	 5067-5685	Clamp Guide Kit-IF-II (for Infinity II modules except G7116A/B MCT)
2	 2110-1486	Fuse 2 AT250 V
1	 5063-6527	Tubing, Silicon Rubber, 1.2 m, ID/OD 6 mm/9 mm
1	 5500-1156	T-Tube Connector ID6.4 (not orderable separately)

9

Identifying Cables

This chapter provides information on cables used with the modules.

Cable Overview 119

Analog Cables 121

Remote Cables 123

BCD Cables 127

CAN/LAN Cables 129

RS-232 Cables 130

USB 131

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

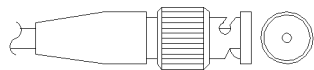
p/n	Description
5181-1516	CAN cable, Agilent module to module, 0.5 m
5181-1519	CAN cable, Agilent module to module, 1 m

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)

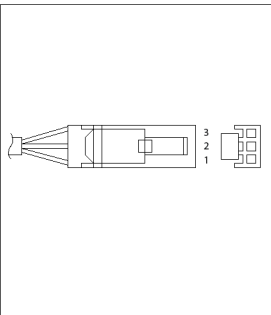
RS-232 cables	p/n	Description
	RS232-61601	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It is also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3,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

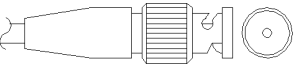


One end of these cables provides a BNC connector to be connected to Agilent modules. The other end depends on the instrument to which connection is being made.

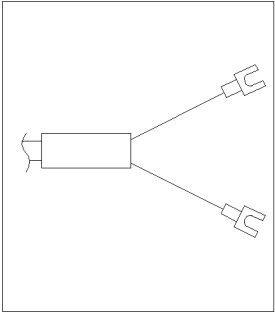
Agilent Module to 35900 A/D converters

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 +

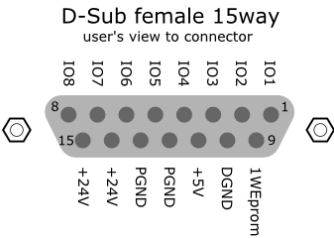
Agilent Module to General Purpose

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))

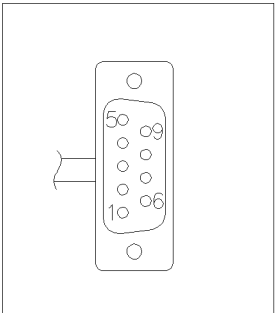
Table 8: 5188-8057 ERI to APG and RJ45

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		

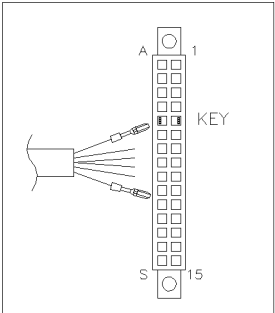


One end of these cables provides an Agilent Technologies APG (Analytical Products Group) remote connector to be connected to Agilent modules. The other end depends on the instrument to be connected to.

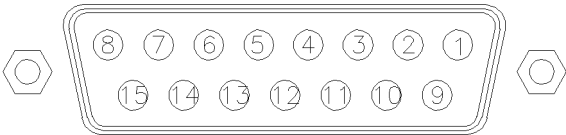
Agilent Module to Agilent 35900 A/D Converters

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

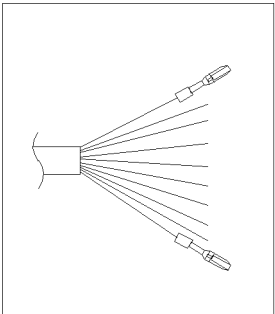
p/n 01046-60201	Wire Color	Pin Agilent module	Signal Name	Active (TTL)
	White	1	Digital ground	
	Brown	2	Prepare run	Low
	Gray	3	Start	Low
	Blue	4	Shut down	Low
	Pink	5	Not connected	
	Yellow	6	Power on	High
	Red	7	Ready	High
	Green	8	Stop	Low
	Black	9	Start request	Low

BCD Cables

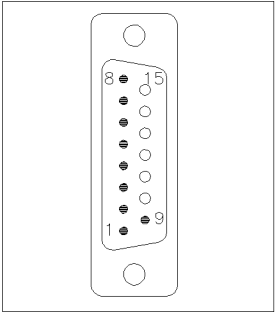


One end of these cables provides a 15-pin BCD connector to be connected to the Agilent modules. The other end depends on the instrument to be connected to

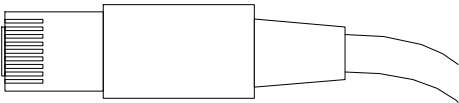
Agilent Module to General Purpose

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)

10

Hardware Information

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

General Hardware Information 133

Firmware Description 133

Electrical Connections 135

Interfaces 139

Instrument Layout 144

Early Maintenance Feedback (EMF) 144

Module-Specific Hardware Information 146

Config Switch Settings of the G1170A Valve Drive 146

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.

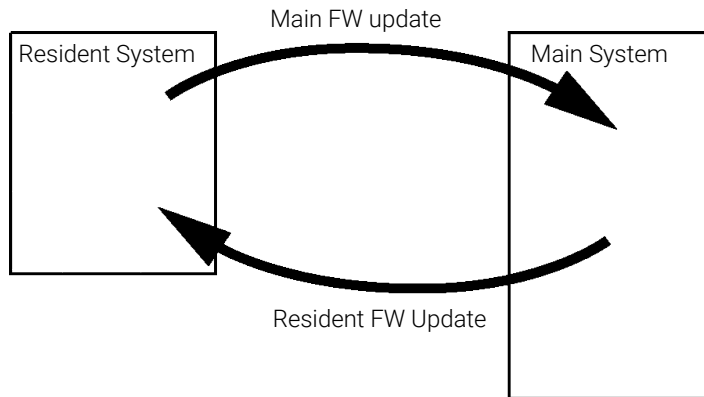


Figure 15: Firmware update mechanism

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.
- 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.

- 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.

Front and Rear View of the Module

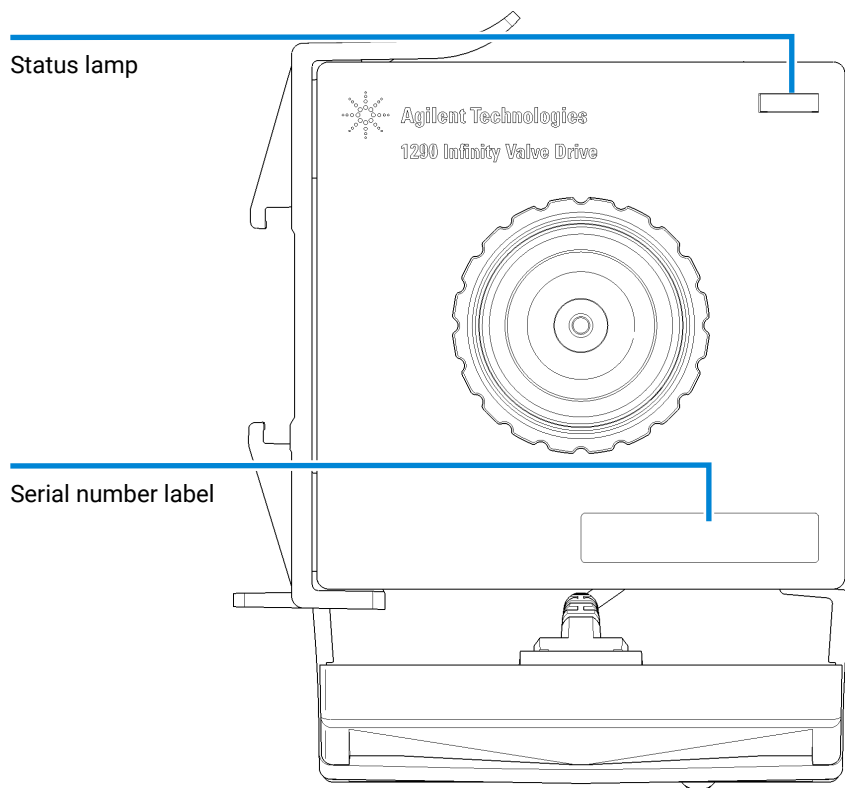


Figure 16: Front view of 1290 Infinity III Valve Drive

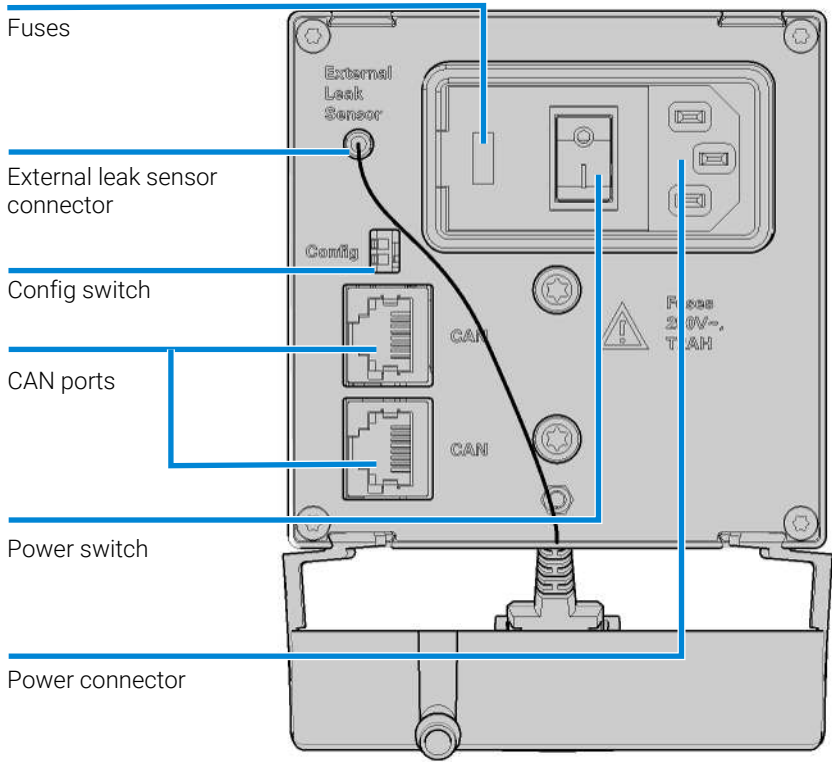


Figure 17: Rear view of 1290 Infinity III Valve Drive

Serial Number Information

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

CCYWWSSSSS	Format
CC	country of manufacturing <ul style="list-style-type: none">• DE = Germany• JP = Japan• CN = China
YWW	year and week of last major manufacturing change, e.g. 820 could be week 20 of 1998 or 2008
SSSSS	real serial number

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 9: Agilent InfinityLab LC Series interfaces

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

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

NOTE

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

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

Overview Interfaces

CAN

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

LAN

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

NOTE

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

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

USB

The USB interface replaces the RS-232 Serial interface in new generation modules. For details on USB refer to [USB \(Universal Serial Bus\)](#) on page 144.

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 10: ERI signal distribution

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

Special Interfaces

There is no special interface for this module.

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

Config Switch Settings of the G1170A Valve Drive

Configuration switch settings

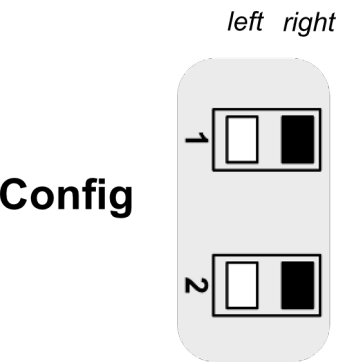


Figure 18: Config Switches

Table 11: Dip switches settings for G1170A

Mode select	1	2
Default	right	right
Coldstart	right	left
Boot resident	left	right
Not supported	left	left

Special Settings

Boot-Resident

Firmware update procedures may require this mode in case of firmware loading errors (main firmware part). If you use the following switch settings and power the instrument up again, the instrument firmware stays in the 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.

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.

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

General Safety Information 149

Safety Standards 149

General 149

Before Applying Power 150

Ground the Instrument 150

Do Not Operate in an Explosive Atmosphere 151

Do Not Remove the Instrument Cover 151

Do Not Modify the Instrument 151

In Case of Damage 151

Solvent Information 152

Safety Symbols 154

Material Information 156

General Information About Solvent/Material Compatibility 156

At-a-Glance Details About Agilent Capillaries 162

Waste Electrical and Electronic Equipment (WEEE) Directive 166

Radio Interference 167

RFID Statement 168

Sound Emission 170

Agilent Technologies on Internet 171

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 154.

Ground the Instrument

WARNING

Missing electrical ground

Electrical shock

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

Do Not Operate in an Explosive Atmosphere

WARNING

Presence of flammable gases or fumes

Explosion hazard

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

Do Not Remove the Instrument Cover

WARNING

Instrument covers removed

Electrical shock

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

Do Not Modify the Instrument

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

In Case of Damage

WARNING

Damage to the module

Personal injury (for example electrical shock, intoxication)

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

Solvent Information

WARNING

Toxic, flammable and hazardous solvents, samples and reagents

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

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

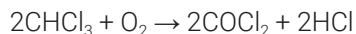
NOTE

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

Recommendations on the Use of Solvents

Observe the following recommendations on the use of solvents.

- Brown glass ware can avoid growth of algae.
- Follow the recommendations for avoiding the growth of algae, see the pump manuals.
- Small particles can permanently block capillaries and valves. Therefore, always filter solvents through 0.22 µm filters.
- Avoid or minimize the use of solvents that may corrode parts in the flow path. Consider specifications for the pH range given for different materials such as flow cells, valve materials etc. and recommendations in subsequent sections.
- Avoid the use of the following steel-corrosive solvents:
 - solutions of alkali halides and their respective acids (for example, lithium iodide, potassium chloride, and so on),
 - high concentrations of inorganic acids like sulfuric acid and nitric acid, especially at higher temperatures (if your chromatography method allows, replace by phosphoric acid or phosphate buffer which are less corrosive against stainless steel),
 - halogenated solvents or mixtures which form radicals and/or acids, for example:



This reaction, in which stainless steel probably acts as a catalyst, occurs quickly with dried chloroform if the drying process removes the stabilizing alcohol,

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

Recommended Wash Solvents

- water
- ethanol









- methanol
- water/acid (especially for basic compounds)
- water/base (especially for acidic compounds)
- water/acetonitrile

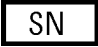





NOTE

For different wash solvents as mentioned above, verify that the wash solvent is suitable for the silicone wash tubing.

Safety Symbols

Table 12: 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. <ul style="list-style-type: none">– 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. <ul style="list-style-type: none">– Do not proceed beyond a caution until you have fully understood and met the indicated conditions.

Material Information

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

General Information About Solvent/Material Compatibility

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

Disclaimer

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

MP35N

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

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

Polyphenylene Sulfide (PPS)

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

PEEK

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

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

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

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

Polyimide

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

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

Polyethylene (PE)

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

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

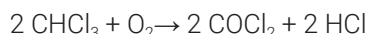
Tantalum (Ta)

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

Stainless Steel (SST)

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

- Solutions of alkali halides, their respective acids (for example, lithium iodide, potassium chloride) and aqueous solutions of halogens.
- High concentrations of inorganic acids like nitric acid, sulfuric acid, and organic solvents especially at higher temperatures (replace, if your chromatography method allows, by phosphoric acid or phosphate buffer, which are less corrosive against stainless steel).
- Halogenated solvents or mixtures, which form radicals and/or acids, for example:



This reaction, in which stainless steel probably acts as a catalyst, occurs quickly with dried chloroform if the drying process removes the stabilizing alcohol.

- Chromatographic grade ethers, which can contain peroxides (for example, THF, dioxane, diisopropyl ether). Such ethers should be filtered through dry aluminum oxide, which adsorbs the peroxides.
- Solutions of organic acids (acetic acid, formic acid, and so on) in organic solvents. For example, a 1 % solution of acetic acid in methanol will attack steel.
- Solutions containing strong complexing agents (for example, EDTA, ethylenediaminetetraacetic acid).
- Mixtures of carbon tetrachloride with isopropanol or THF.

Titanium (Ti)

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

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

Diamond-Like Carbon (DLC)

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

Fused Silica and Quartz (SiO₂)

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

Gold

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

Zirconium Oxide (ZrO₂)

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

Platinum/Iridium

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

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

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

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

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

Sapphire, Ruby, and Al₂O₃-Based Ceramics

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


At-a-Glance Details About Agilent Capillaries

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

Syntax for capillary description

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

Table 13: Example for a capillary description



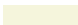










Code provided with the part	Meaing of the code
Color code: 	Material of the product is MP35N, the inner diameter is 0.20 or 0.25 mm
Capillary	The part is a connection capillary
MP35N	Material of the part is MP35N
0.25 x 80 mm	The part has an inner diameter of 0.25 mm and a length of 80 mm
SI/SI	Left 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 14](#) on page 163
- Type: [Table 15](#) on page 163
- Material: [Table 16](#) on page 164
- Dimension: [Table 17](#) on page 164
- Fittings: [Table 18](#) on page 165

Color Coding Guide

Table 14: 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 15: 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 16: 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 17: 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 18: Fitting left/fitting right (indicates which fitting is used on both ends of the capillary)

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

Waste Electrical and Electronic Equipment (WEEE) Directive

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

**NOTE**

Do not dispose of in domestic household waste
To return unwanted products, contact your local Agilent office, or see <https://www.agilent.com> for more information.

Radio Interference

Cables supplied by Agilent Technologies are screened to provide optimized protection against radio interference. All cables are in compliance with safety or EMC regulations.

Test and Measurement

If test and measurement equipment is operated with unscreened cables, or used for measurements on open set-ups, the user has to assure that under operating conditions the radio interference limits are still met within the premises.

RFID Statement

Brasil

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

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

Canada

Statement according to RSS GEN Issue 5:

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

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

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

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

Mexico

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

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

Thailand

เครื่องโทรคมนาคมและอุปกรณ์นี้มีความสอดคล้องตามมาตรฐานหรือข้อกำหนดทางเทคนิคของ กสทช.

This telecommunication equipment conforms to NTC/NBTC technical requirement.

USA

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

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

CAUTION

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

NOTE

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

Table 19: Operating frequencies and maximum power levels

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

Sound Emission

Sound Pressure

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

Schalldruckpegel

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

Agilent Technologies on Internet

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

<https://www.agilent.com>

In This Book

This manual contains technical reference information about the Agilent 1290 Infinity III Valve Drive (G1170A) and the Agilent InfinityLab Quick Change Valves.

The manual describes the following:

- introduction,
- specifications,
- configuration,
- optimizing,
- troubleshooting and diagnostics,
- maintenance,
- parts identification,
- hardware information,
- safety and related information.

www.agilent.com

© Agilent Technologies Inc. 2011-2024
Edition: 10/2024

Document No: SD-29000412 Rev. D

