

Agilent InfinityLab

Quick Change Valves G4234A/C

Instructions

This technical note describes the installation and application of the Agilent InfinityLab Quick Change 6 Column Selector Valve Heads G4234A and G4234C in a 1290 Infinity II Multicolumn Thermostat (MCT).

Contents

Typical Applications	2
Multicolumn Selection	2
Method Development	3
Delivery Checklist	4
Specifications	6
Install the Valve Heads	7
Remove the Transportation Lock and the Valve Dummy	7
Install the Valve Head and Connect Capillaries	9
Install the Capillaries	13
Valve Parts	14
Replacement Parts	14
Valve Head Parts	15



Typical Applications

Multicolumn Selection

Advantages:

- Increase productivity
- Higher instrument up-time

Quickly change between up to six different stationary phases for different applications, or use identical stationary phases in columns with different dimensions for either faster run-times (short columns) or higher resolution (long columns) or for loading studies with different internal diameters.

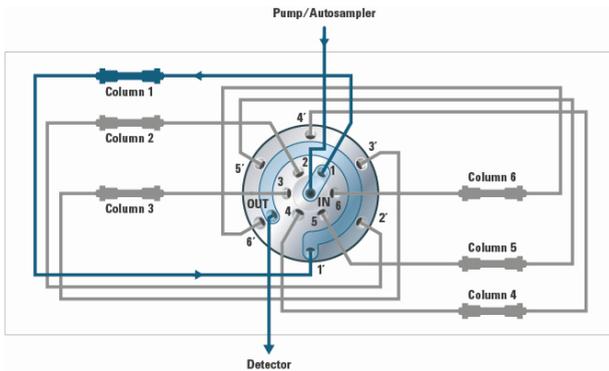


Figure 1 Multiple column selection, pos. 1 (example of schematic setup for 6-column selector)

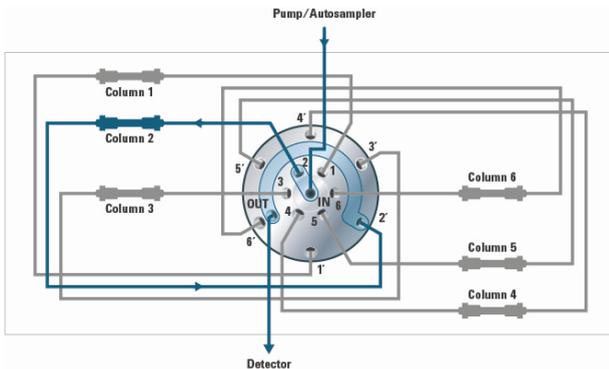


Figure 2 Multiple column selection, pos. 2 (example of schematic setup for 6-column selector)

Method Development

Advantages:

- Faster method development
- Automated method development possible

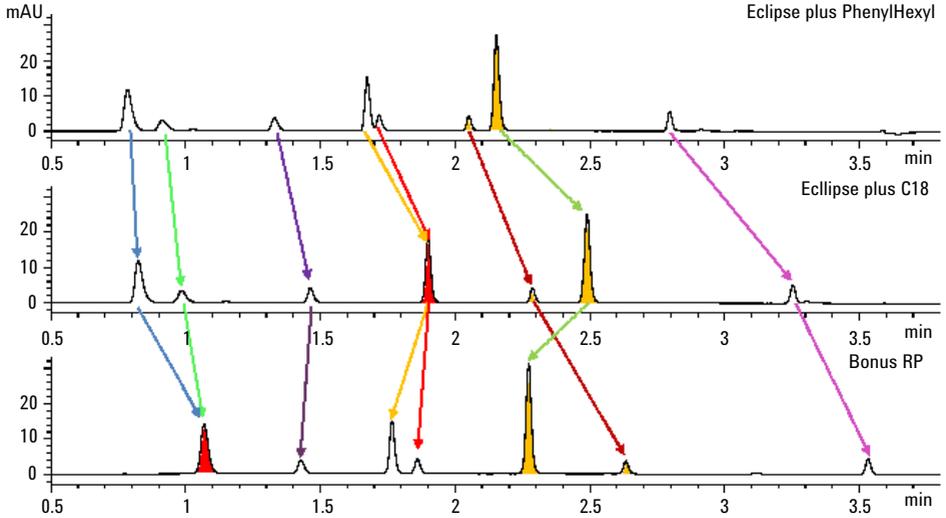


Figure 3 Totally different chromatographic results by using the same sample but three different stationary phases

Delivery Checklist

Check the content of the delivery. You should have received the following:

G4234A:

Item	p/n	Description
1	5067-4284	6-column selector valve head, 800 bar
2	5067-4270	6-Column Selector Capillary Kit 0.12 mm ID, MCT (OPTIONAL)

G4234C:

Item	p/n	Description
1	5067-4273	6-column selector valve head, 1300 bar
2	5067-4270	6-Column Selector Capillary Kit 0.12 mm ID, MCT (OPTIONAL)

6-Column Selector Capillary Kit 0.12 mm ID, MCT (5067-4270), contains:

#	p/n	Description
6	G7116-60015	Quick Connect Heat Exchanger Standard
6	5500-1201	Capillary ST 0.12 mm x 105 mm SL-- PS-LS Quick Connect Heat Exchanger to column
6	G7167-68703	Fitting Intermediate Kit fitting for column, capillary 5500-1201
6	5500-1199	Capillary ST 0.12 mm x 130 mm M4-SL PS-PS Valve to Quick Connect Heat Exchanger
6	5500-1200	Quick Turn Capillary ST 0.12 mm x 130 mm SL/M Column to valve
1	5063-6591	PEEK Fittings 10/PK Column outlet fitting for capillary 5500-1200
1	5500-1202	Capillary ST 0.12 mm x 500 mm M4-SL PS-PS Autosampler to valve
1	5500-1203	Capillary ST 0.12 mm x 280 mm M4-SL PS-PS Valve to detector
1	5500-1204	Capillary ST 0.12 mm x 150 mm M4-M4 PS-PS Valve to valve (bypass)
1	G1375-87326	Waste tube Waste tube incl. fitting
2	5067-6141	M4 Blank nut Block unused ports
1	5023-2504	Hex driver SW-4 slitted Tool for M4 fittings
6	G7116-68004	Column Holder Clamp (2/PK) for Infinity II
1	5043-0915	Fitting mounting tool for bio-inert capillaries
1	5067-6654	Number Kit 1-8 colored

Specifications

Table 1 G4234A (5067-4284), 6-column selector valve head, 800 bar

Type	Specification
Maximum pressure	800 bar
Typical application	6-column selection
Port size	Accepts M4 male threaded fittings
Liquid contacts	PEEK, Stainless Steel
pH range	0 – 14*

* incompatible with some mineral acids. For more information see Solvent Information.

Table 2 G4234C (5067-4273), 6-column selector valve head, 1300 bar

Type	Specification
Maximum pressure	1300 bar
Typical application	6-column selection
Port size	Accepts M4 male threaded fittings
Liquid contacts	PEEK, Stainless steel
pH range	0 – 14*

* incompatible with some mineral acids. For more information see Solvent Information.

Install the Valve Heads

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 7. 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.

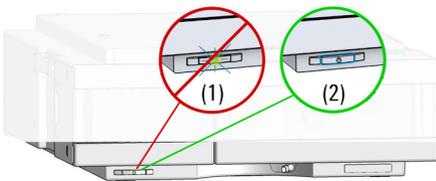
When all capillaries are installed, push the valve back into its housing, see [“Install the Valve Head and Connect Capillaries”](#) on page 9.

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).

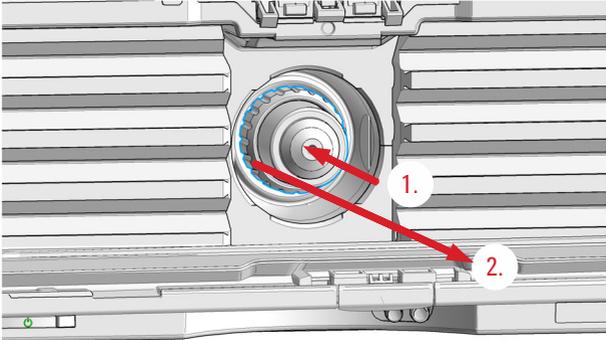
For the installation of a valve head to a G1170A Infinity Valve Drive you can ignore the steps that describe the MCT features of the transportation lock and spring loaded valve drive.

- 1 Switch off the module.

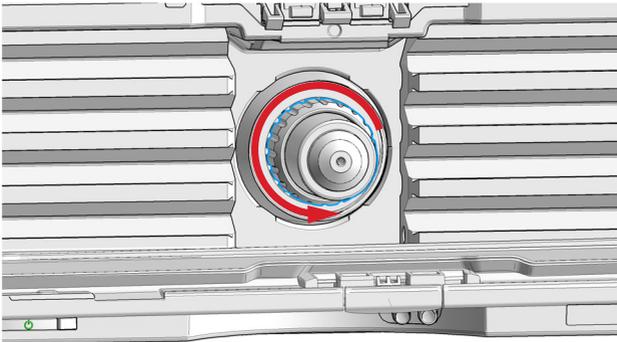


- 2 When unscrewing the transportation lock, push it back until the last screw is removed - the valve rail is spring-loaded.

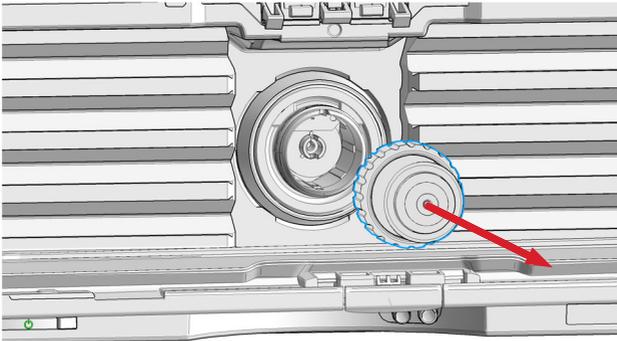
3 Press on the valve dummy (1.) to release it (2.) (spring-loaded valve rail).



4



5



Install the Valve Head and Connect Capillaries



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

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.
-

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

CAUTION

Wrong combination of fitting with valve

The InfinityLab Quick Turn Fitting (5067-5966) is not compatible with the G5639A Bio-inert 4-Column Selector Valve. Misuse can lead to extra dead volume and leaks.

- As fitting, use UHP fitting (5067-5403) instead.

NOTE

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

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

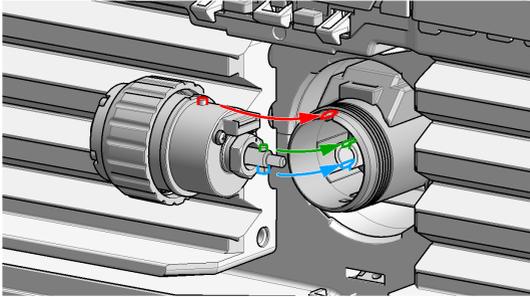
To allow correct valve identification, power off the valve drive for at least 10 s.

NOTE

For firmware requirements see Information on new RFID Tag Assembly Version Technical Note (01200-90133) which is included to each valve head.

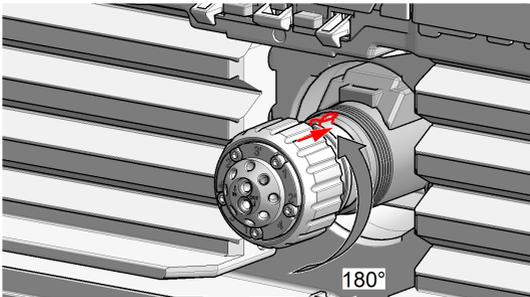
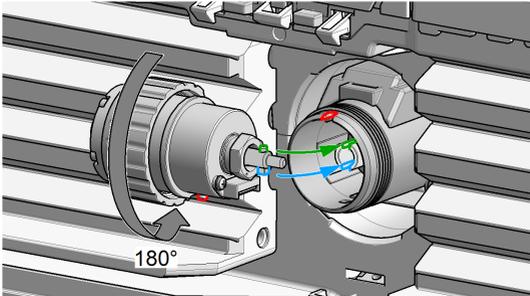
The following procedure shows the valve head installation with an G7116A/B (MCT) module as an example. For other modules it is similar.

- 1 Insert the valve head into the valve shaft.

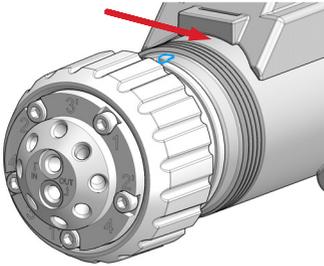


OR

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.



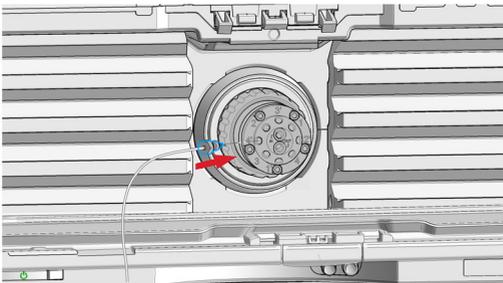
- 2** When the outer pin is locked into the groove, manually screw the nut onto the valve head.



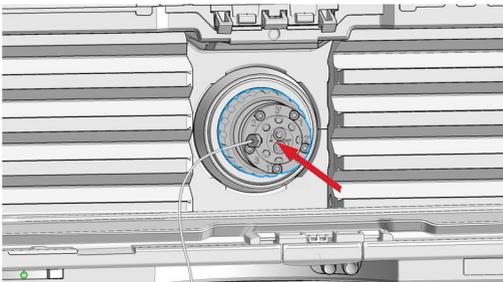
NOTE

Fasten the nut manually. Do not use any tools.

- 3** Install all required capillary connections to the valve.



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

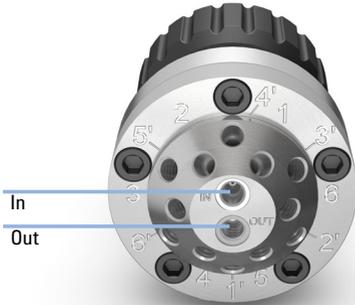


- 5** Power on or power-cycle your module, so the valve head gets recognized during module initialization.

Install the Capillaries

1 Install the in and out connectors.

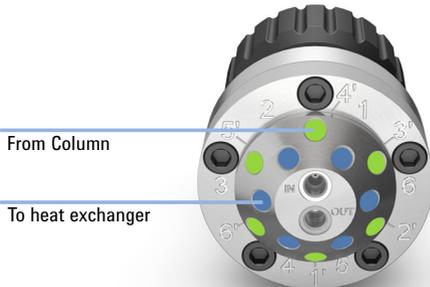
- from sampler to the valve (Capillary ST 0.12 mm x 500 mm M4-SL PS-PS (5500-1202))
- from valve to the detector (Capillary ST 0.12 mm x 280 mm M4-SL PS-PS (5500-1203))



The *In* port is hydraulically connected to the column inlet ports 1-6 on the inner ring while the *Out* port connects to the column outlet ports 1'-6' on the outer ring.

2 Install the column inlet and outlet connections.

- ports 1-6 for connections from valve to the heat exchanger (Capillary ST 0.12 mm x 130 mm M4-SL PS-PS (5500-1199)) or waste line (Waste tube (G1375-87326))
- ports 1'-6' for connections from column outlet to valve (Quick Turn Capillary ST 0.12 mm x 130 mm SL/M (5500-1200)), use fingertight PEEK fittings for connecting the column outlet



Valve Parts

Replacement Parts

Table 3 Replacement parts

Valve	Rotor Seal	Stator Head	Stator Screws (pack of 10)	Bearing Ring
5067-4284 6-column selector valve head, 800 bar	5068-0298 (PEEK)	5068-0241	5068-0089	1535-4045
5067-4273 6-column selector valve head, 1300 bar	5068-0242 (PEEK)	5068-0241	5068-0089	1535-4045

Valve Head Parts

NOTE

The figure below illustrates replacement parts for the valve heads, with the 12ps/13pt Selector valve as an example. The valves can vary in their appearance and do not necessarily include all of the illustrated parts. Neither, every spare part is available for each flavor of the valve.

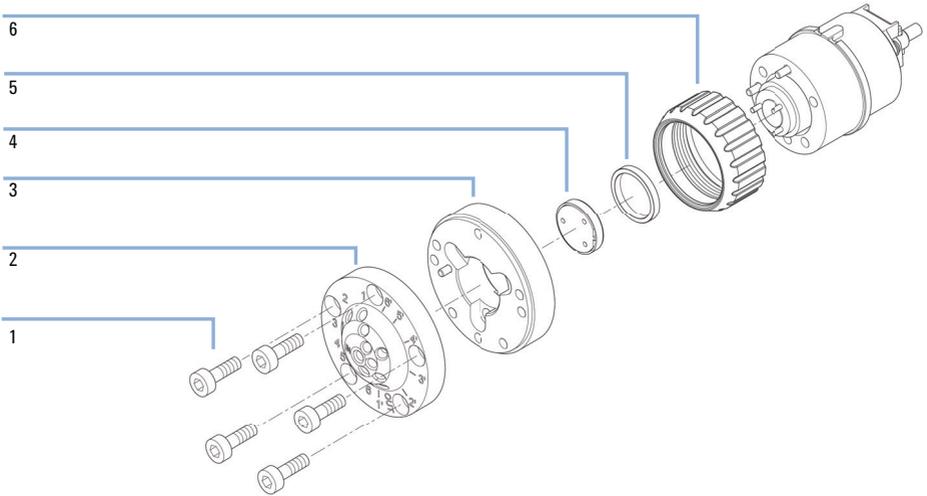


Figure 4 Valve Head Parts (example)

-
- | | |
|---|--|
| 1 | Stator screws |
| 2 | Stator head assembly |
| 3 | Stator ring (available for service only) |
| 4 | Rotor seal |
| 5 | Bearing ring |
| 6 | Spanner nut (available for service only) |
-



G4234-90010

Part Number:
G4234-90010 Rev. B

Edition: 01/2019
Printed in Germany

© Agilent Technologies, Inc
2017-2019

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