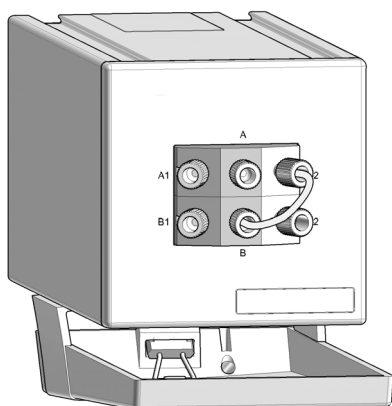




# Agilent 1260 Infinity II Fraction Collector Cluster Valve

## Technical Note

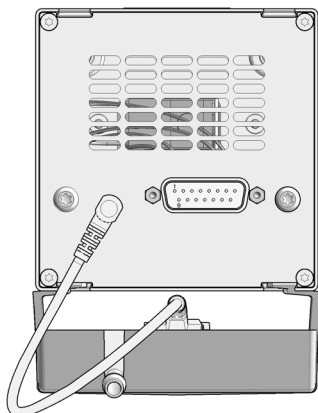
### Overview



**Figure 1** 1260 Infinity II Fraction Collector Cluster Valve



**Agilent Technologies**



**Figure 2** Rear view

For installation of the 1260 Infinity II Fraction Collector Cluster Valve, please refer to the Agilent 1260 Infinity II Fraction Collector Cluster Valve Installation Note (G9322-90120).

p/n	Description
G9322-64000	Fraction Collector Cluster Valve
G9322-68002	Bridging Capillary Set
5188-8059	ERI-Extension-Cable 1.2 m
5043-0271	Holder leak plane
5043-0270	Leak plane
5067-4792	Leak sensor assembly

# Specifications

## Physical Specifications

**Table 1**    Physical Specifications 1260 Infinity II Fraction Collector Clustering Valve (G9322A)

Type	Specification	Comment
Weight	0.6 kg (1.4 lbs)	
Dimensions (height x width x depth)	90 x 90 x 205 mm (3.5 x 3.5 x 8.1 inches)	
Line Voltage	24 V DC / 5 V DC	
Power consumption	3 W	
Ambient operating temperature	4 – 40 °C (39 – 104 °F)	
Ambient non-operating temperature	-40 – 70 °C (-40 – 158 °F)	
Humidity	<95 % r.h. at 40 °C (104 °F)	Non-condensing
Operating altitude	Up to 3000 m (9842 ft)	
Non-operating altitude	Up to 4600 m (15092 ft)	For storing the module
Safety standards: IEC, EN, CSA, UL	Pollution degree 2	For indoor use only
ISM Classification	ISM Group 1 Class B	According to CISPR 11

# Performance Specifications

**Table 2** Performance Specifications 1260 Infinity II Fraction Collector Clustering Valve (G9322A)

Type	Specification	Comment
External Leak Sensor	Yes	Via a 2.5 mm jack plug
Flow rate	0 – 200 mL/min	
Maximum Pressure	8 bar (116 psi)	
Valve Identification	Via EPROM	
Instrument Control	LC & CE Drivers A.02.19 or above Instrument Control Framework (ICF) A.02.04 or above Instant Pilot (G4208A) with firmware B.02.22 or above InfinityLab LC Companion (G7108A) Lab Advisor B.02.11 or above	For details about supported software versions refer to the compatibility matrix of your version of the LC & CE Drivers
Communications	ERI (1x)	G9322A is a hosted ERI module. The LC stack needs to contain an ERI connector.
Maintenance and safety-related features	Leak handling with external leak sensor and leak pane. Leak detection, safe leak handling, leak output signal for shutdown of pumping system, and low voltages in major maintenance areas Extensive diagnostics, error detection and display with Agilent Lab Advisor software	

**Table 2**    Performance Specifications 1260 Infinity II Fraction Collector Clustering Valve (G9322A)

Type	Specification	Comment
GLP features	Early maintenance feedback (EMF) for continuous tracking of instrument usage with user-settable limits and feedback messages. Electronic records of maintenance and errors	
Housing	All materials recyclable	

# Waste Electrical and Electronic Equipment Directive

## Abstract

The Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC), adopted by EU Commission on 13 February 2003, is introducing producer responsibility on all electric and electronic appliances starting with 13 August 2005.

### NOTE

This product complies with the WEEE Directive (2002/96/EC) marking requirements. The affixed label indicates that you must not discard this electrical/electronic product in domestic household waste.

Product Category:

With reference to the equipment types in the WEEE Directive Annex I, this product is classed as a Monitoring and Control Instrumentation product.

---



### NOTE

Do not dispose of in domestic household waste

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

---





G9322-90130

Part Number:  
G9322-90130 Rev. B

Edition: 12/2017  
Printed in Germany

© Agilent Technologies, Inc 2017

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