

# 1220 Infinity II

## Active Inlet Valve Upgrade Kit

### Technical Note

This note describes the procedures to install an AIV upgrade kit to 1220 Infinity II LC. This upgrade kit only applies to 1220 Infinity II LC modules with minimum Firmware revision B.07.10. With the AIV upgrade, some applications (e.g. involving buffers) may work even better. This upgrade kit also gives customers the backward compatibility, if the analysis methods are transferred from other AIV systems.

## Contents

[Delivery Checklist](#) 2

[Install the Active Inlet Valve](#) 2



**Agilent Technologies**

# Delivery Checklist

p/n	Description
0100-2298	Adapter, PEEK int. 1/4-28 to ext. 10-32 (Used for G4286B Isocratic pump)
G1311-67304	Connecting tube
G1312-60020	Cartridge for active inlet valve 600 bar
G1312-60025	Active inlet valve without cartridge

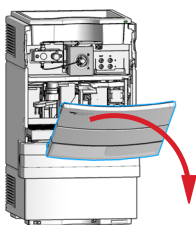
## Install the Active Inlet Valve

Tools required	Description
	Wrench, 14 mm

Parts required	p/n	Description
	G5699B	1220 Infinity II AIV Upgrade Kit

- Preparations**
- Switch off the instrument at the main power switch.
  - Use an optional solvent shutoff valve or lift up solvent filters in solvent reservoirs to avoid leakages.

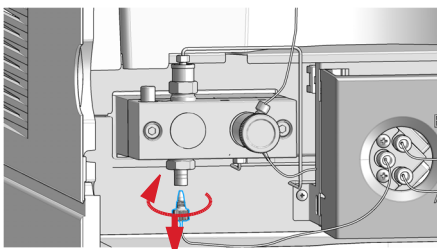
**1** Remove the upper front cover.



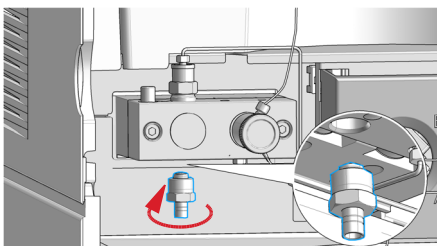
- 2 Disconnect the solvent inlet tube from the inlet valve.

## NOTE

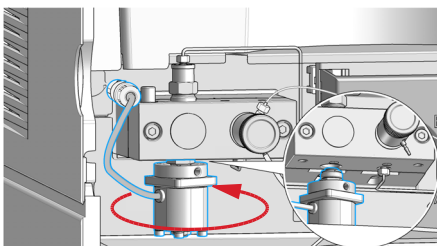
Beware of leaking solvents due to hydrostatic pressure.



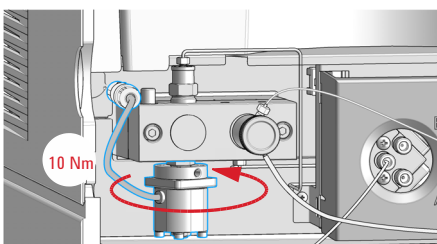
- 3 Using a 14 mm wrench, loosen the passive inlet valve and remove the valve from the pump head.



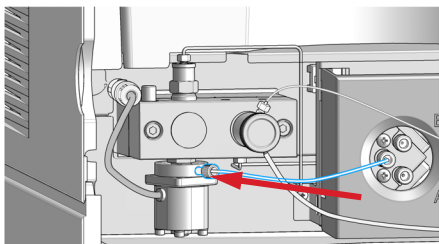
- 4 Insert the active inlet valve into the pump head. Using the 14 mm wrench, turn the nut until it is hand tight. Position the valve such that the solvent inlet tube connection points towards the front.



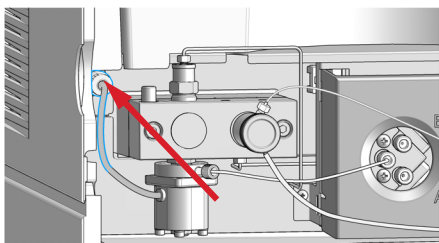
- 5 Tighten the active inlet valve using a 14 mm wrench (approx. 10 Nm).



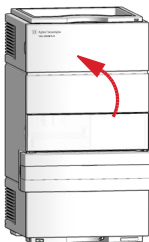
- 6** Use the new solvent tubing provided in update kit to connect the flow path between AIV and DCGV.



- 7** Connect the active inlet valve cable to the connector in the Z-panel.



- 8** Close the upper front cover.



Part Number:  
G5699-90000

Edition: 10/2016  
Printed in Germany

© Agilent Technologies, Inc 2016

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