



# Agilent 1220 Infinity II LC

## Column Compartment Upgrade Kit

### Installation Note

This note describes the procedures to install an Column Compartment upgrade kit to an existing Agilent 1220 Infinity II LC without Column Compartment.

## Introduction

With the Column Compartment Upgrade Kit, the following instruments can be upgraded:

- G4286B
- G4288B
- G4288C

After installation, the Column Compartment can be heated from 5 °C above ambient temperature to 80 °C (see [Table 1](#) on page 1 for details).

**Table 1** Specifications

Type	Specification
Temperature range	<ul style="list-style-type: none"><li>• 5 °C above ambient to 80 °C</li><li>• 10 °C above ambient to 80 °C (G4294B with DAD)</li></ul>
Temperature stability	±0.15 °C, constant composition and flow rate
Temperature accuracy	±0.8 °C
Column capacity	one 25 cm column
Internal volume	6 µL

## Delivery Checklist

p/n	Description
G4280-60017	Heater door assembly
G4280-60040	Complete column compartment assembly
G1316-87300	Capillary ST 0.17 mm x 90 mm S/S
01090-87304	Capillary ST 0.17 mm x 280 mm S/S

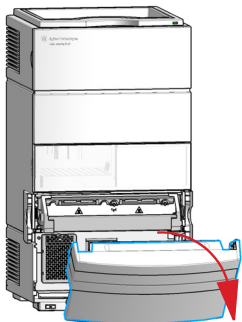
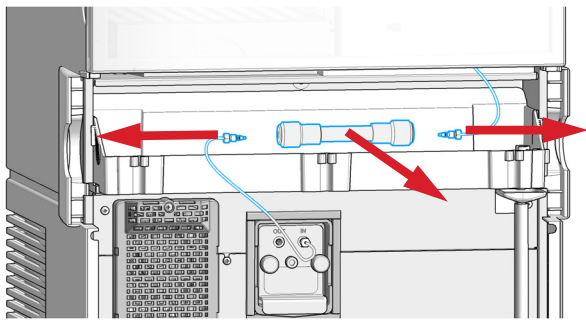
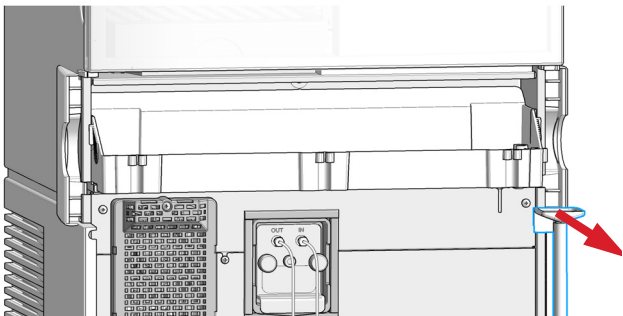
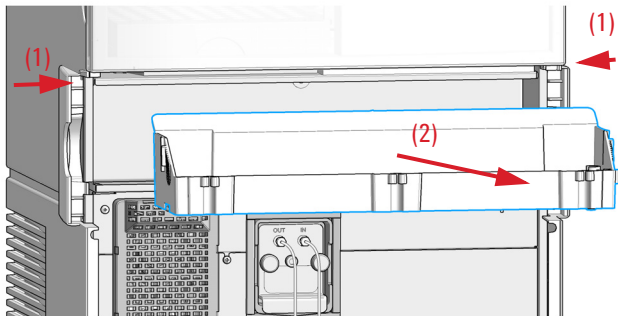
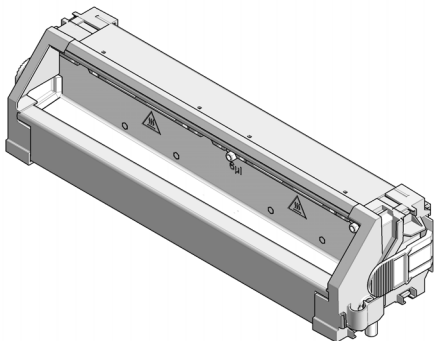
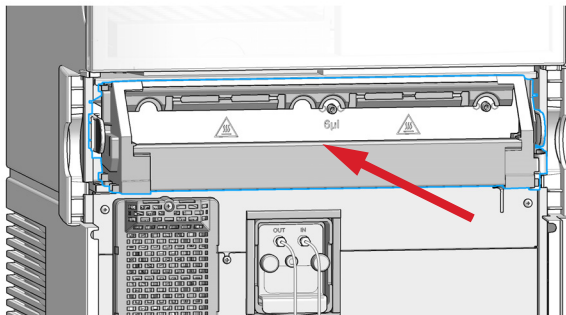


## Install the Column Compartment Upgrade Kit

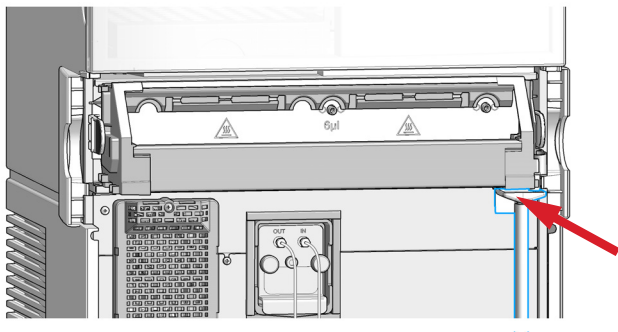
<b>Parts required</b>	<b>p/n</b>	<b>Description</b>
	G4297A	1220 Infinity II Column Compartment Kit

**Software required** Lab Advisor Software

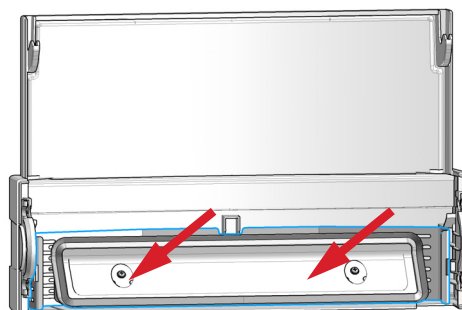
**Preparations** Switch off the instrument.

<p><b>1</b> Remove the lower front cover.</p> 	<p><b>2</b> Disconnect the column and remove it.</p> 
<p><b>3</b> Remove the leak drain tube.</p> 	<p><b>4</b> Press the knurled portion at either side of the column tray inwards (1) and remove the column tray (2).</p> 
<p><b>5</b> Unpack the Column Compartment upgrade kit and separate the two parts.</p> 	<p><b>6</b> Click the Column Compartment into position in place of the column tray.</p>  <p>The electrical connection to the Column Compartment is made automatically.</p>

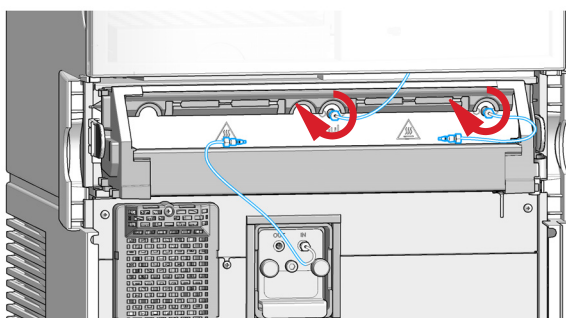
**7** Replace the leak drain tube.



**8** Click the Column Compartment insulation into place in the lower front cover, with the cutout in the Column Compartment insulation support at the bottom.

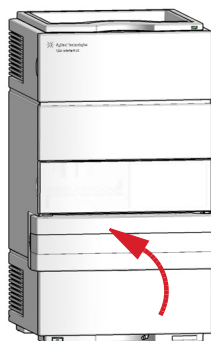


**9** Install and place the capillaries into the column compartment.



**10** Install the column and reconnect the capillaries.

**11** Close the lower front cover.



**Next Steps:**

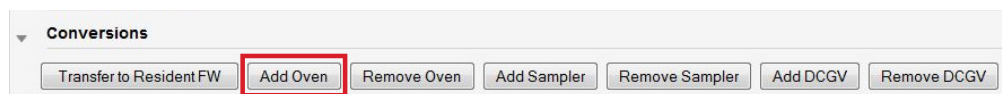
- 12** Configure the Column Compartment in the Lab Advisor software.
- 13** Power cycle the instrument and start your Chromatographic Data System.
- 14** Use Auto configure and the instrument now should be registered under its new main assembly number.

## Configure the Instrument in the Software

- 1 Start the LabAdvisor software.
- 2 Power on and connect your Instrument to LabAdvisor.
- 3 Open **Instrument Control**.



- 4 Open **Conversions** in any of the sub-modules.
- 5 Click **Add Oven** button.



The hardware configuration change will be finalized with the next power-cycle.

- 6 Power cycle the instrument.
- 7 Start your Chromatographic Data System and use **Auto configure**.

The instrument now should be registered under its new main assembly number.



G4297-90001

Part Number: G4297-90001

Edition: 01/2017  
Printed in Germany

© Agilent Technologies, Inc 2016, 2017

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