

Agilent Multidraw Infinity III Bio LC Upgrade Kit (G7137-68711)

Installation Note

This note describes how to install the Agilent Multidraw Infinity III Bio LC Upgrade Kit (G7137-68711) into an Agilent 1290 Infinity III Bio Multisampler (G7137A).

General Information 1

Delivery Checklist 2

Installing the Multidraw Infinity III Bio LC Upgrade Kit 3

Configuration of the Controller 4

General Information

The Multidraw Infinity III Bio LC Upgrade Kit can be installed in the Agilent 1290 Infinity III Bio Multisampler.

With the kit, you can add a maximum of 500 μL or 1500 μL to the injection volume (100 μL analytical head) of your injector.

The maximum injection volume will vary, depending on the analytical head and sample loop installed.

NOTE

The delay volume of your Multisampler is extended when using the extended seat capillaries from the multidraw kit. When calculating the delay volume of the Multisampler, double the volume of the extended capillaries (used in partial filling mode). The delay volume can be reduced by bypassing the Multisampler once the sample has reached the head of the column, see your *User Manual* for more information.

When using the Multisampler in multidraw mode, the syringe ejects an equivalent of the injected volume into waste. Therefore it is recommended to verify the correct installation of the waste tube to the waste outlet of the Multisampler.

Delivery Checklist

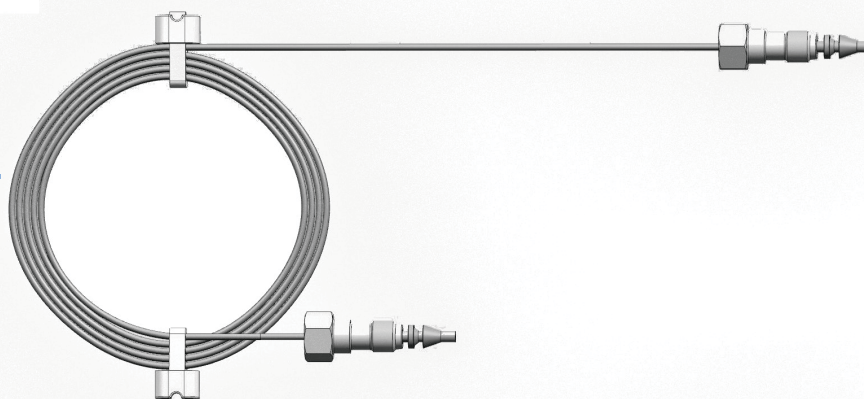
NOTE

Do not connect a waste outlet directly to the plastic waste tube (connected to port 4 of the injection valve). This can lead to siphoning effects that will influence the precision of the Multisampler.

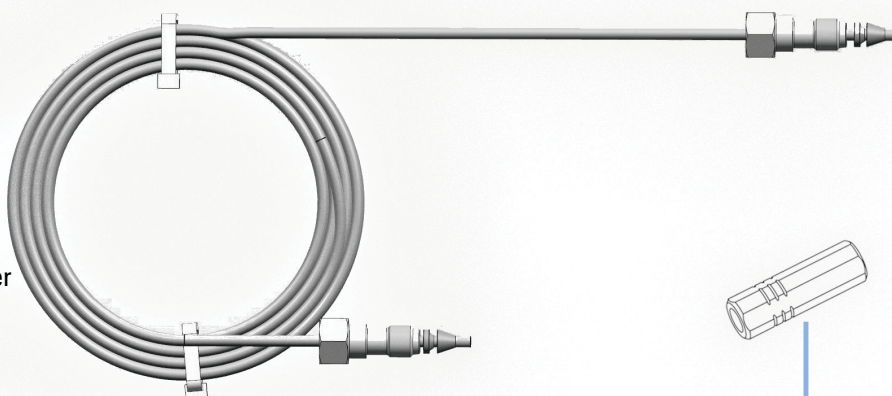
Make sure that all parts and materials have been delivered with the upgrade kit. Report missing or damaged parts to your local Agilent Technologies sales and service office.

p/n	Description
G7137-68307 	Extended seat capillary, MP35N, 500 µL, 0.5 mm id
G7137-68308 	Extended seat capillary, MP35N, 1500 µL, 0.9 mm id
5023-2625 	Union MP35N

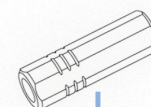
Extended seat capillary
500 µL



Extended seat capillary
1500 µL
(larger external diameter
than 500 µL loop)



Union



NOTE

Both extended seat capillaries have SL fittings, compatible with VICI and IDEX/Rheodyne Valves.

Installing the Multidraw Infinity III Bio LC Upgrade Kit

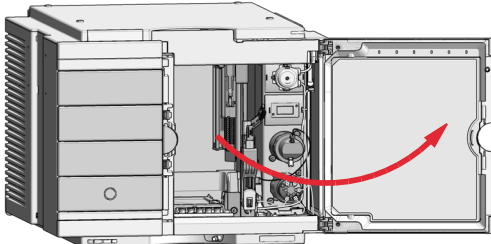
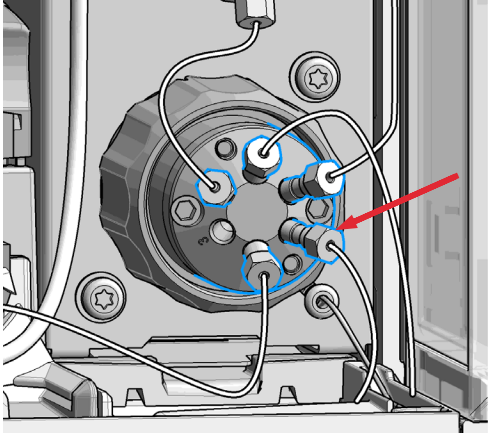
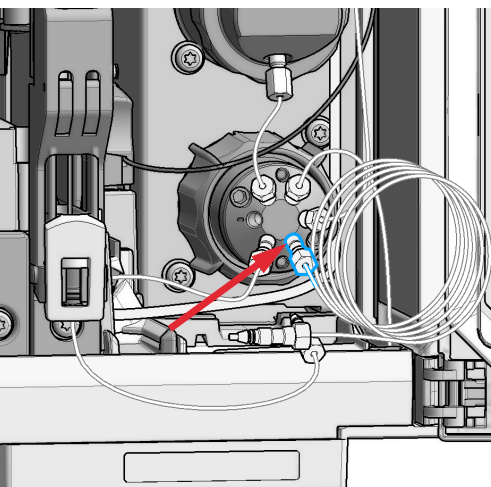
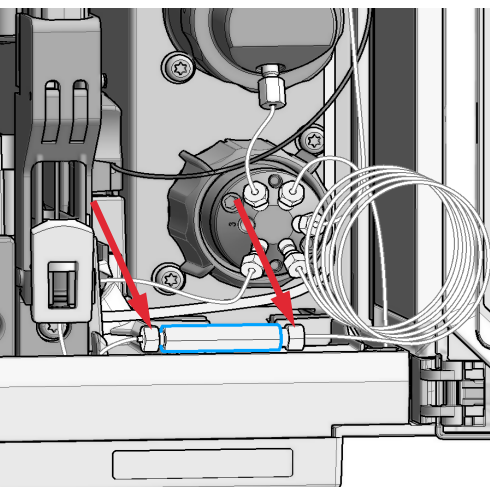
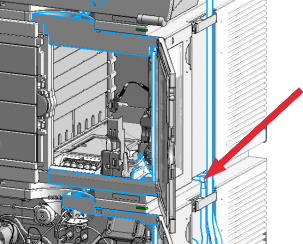
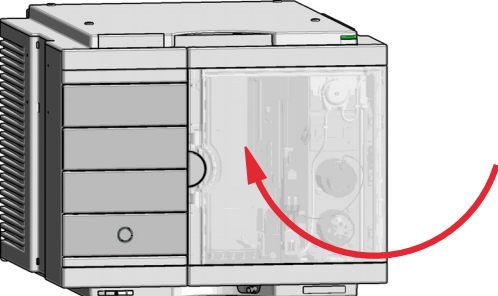
Tools required

p/n

Wrench, 1/4 inch 

Description

1/4 inch wrench (one supplied in the autosampler accessory kit)

1 Open the front door. 	2 Disconnect the seat-capillary fitting from the injection valve (port 5). 
3 Install the 500 or 1500 µL seat extension loop (longer side) to the injection valve (port 5). 	4 Install the union between the seat capillary and the other side of the seat extension loop. Store the extension loop in the leak tray. 
5 Connect the waste tube to the waste guidance. 	6 Close the front door. 
7 Connect the waste tube (shipped with the start up kit of your Multisampler).	

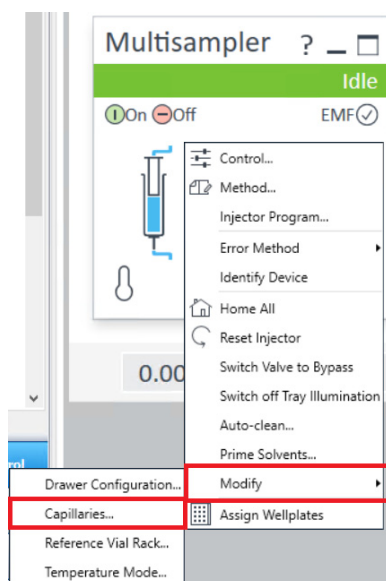
Configuration of the Controller

The configuration of your controller is necessary to enable the multidraw mode.

When setting an injection larger as the configured injection volume the multidraw mode is active. Multiples of the injection syringe volume are stored in the extended seat capillary prior to switching the injection valve.

OpenLab CDS ChemStation C.01.10 Update 3 and OpenLab 2.5 and above

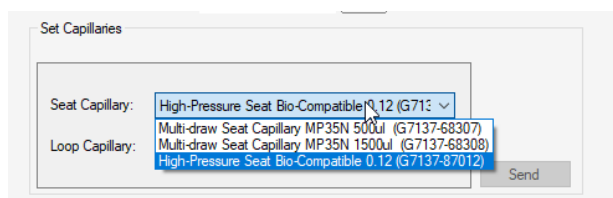
- 1 Select **Capillaries** in the instrument function.



- 2 In the **Configuration** menu, change seat capillary to the value of the installed extended seat capillary and click **OK**.

Lab Advisor B.02.16 and above

- 1 Connect to the stack in **System Overview**.
- 2 Select **Instrument Control** and open **Controls** of the Multisampler.
- 3 On the **Configuration** tab, select the volume for your installed extended seat capillary (500 µL or 1500 µL) from the **Set Capillary** drop-down menu.



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

www.agilent.com

© Agilent Technologies Inc. 2020-2024

Printed in Germany

Edition: 10/2024



Part No: G7137-90110 Rev. Rev. B
Document No: D0004814 Rev. Rev. B

