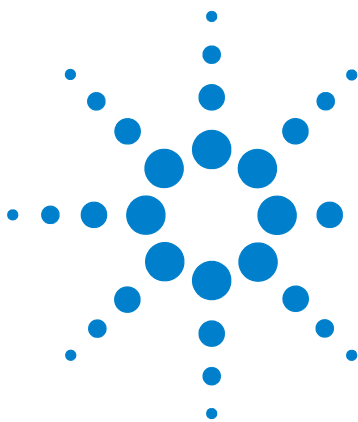




Installation Note for Multidraw Upgrade Kit G1313-68711

General Information	1
Delivery Checklist	2
Installing the Multidraw Upgrade Kit	3
Configuration of the Controller	4
ChemStation Configuration	4
Control Module Configuration	4
Instant Pilot	4

In this note we describe how to install the multidraw upgrade kit into an Agilent 1200 Series autosampler.

General Information

The multidraw upgrade kit can be installed in any Agilent 1200 Series autosampler. With the kit you can add a maximum of 400 μl or 1400 μl to the injection volume (100 μl analytical head) of your injector. The total injection volume is then 500 μl or 1500 μl for the standard autosampler with 100 μl analytical head. If the optional 900 μl analytical head is installed in the autosampler you can add only the 1400 μl injection volume capillary. This allows to expand the injection volume of your autosampler to a maximum of 1800 μl (multi-injection mode only allows storage of integer multiples of the syringe volume in the seat capillary).

NOTE

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

When using the autosampler in multidraw mode the syringe ejects an equivalent of the injected volume into waste. Therefore it is recommended to install a separate waste tube (shipped with the start-up kit of your autosampler) to the waste outlet of the autosampler.



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 autosampler.

Make sure all parts and materials have been delivered with the upgrade kit. The delivery checklist is shown in Table 1. Please report missing or damaged parts to your local Agilent Technologies sales and service office.

Table 1 Multi Draw Upgrade Kit

Description	Quantity	Part Number
Extended seat capillary, 400 μ l	1	G1313-87307
Extended seat capillary, 1400 μ l	1	G1313-87308
Union	1	0100-0900

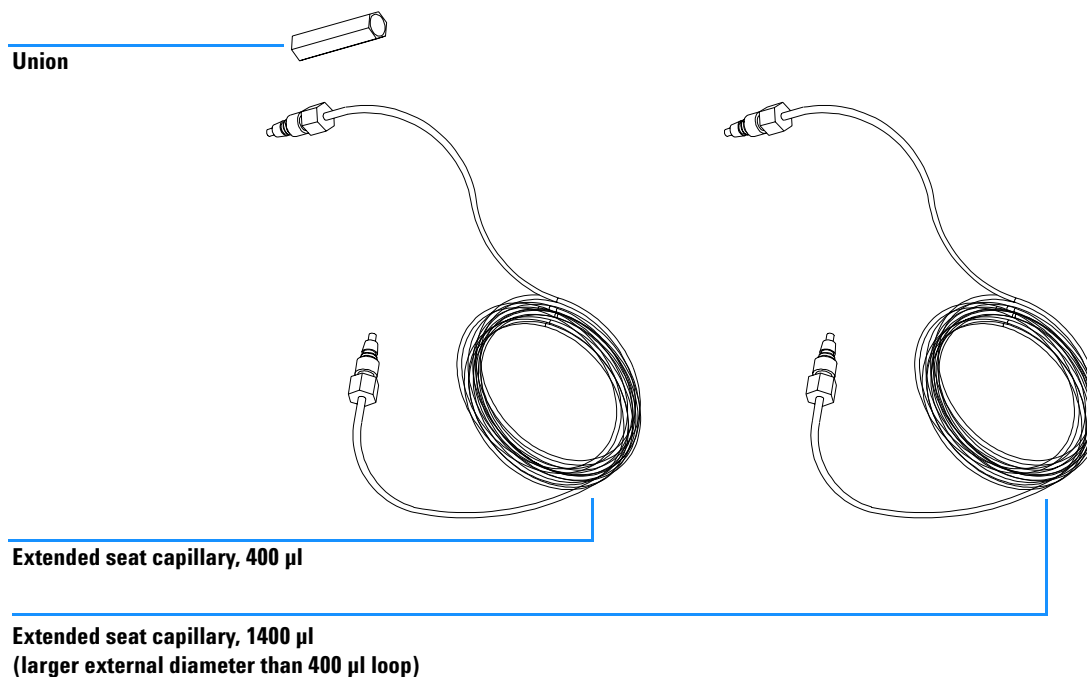


Figure 1 Multidraw Upgrade Kit

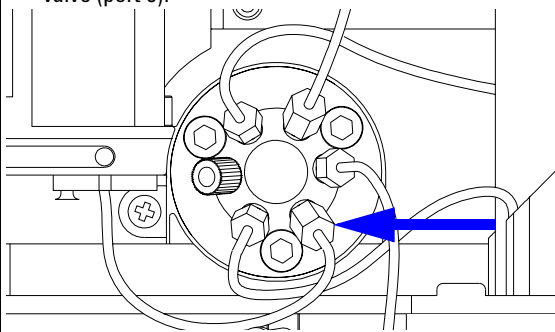
Installing the Multidraw Upgrade Kit

Tools required	1/4 inch wrench 2 × (one supplied in the autosampler accessory kit)
-----------------------	---

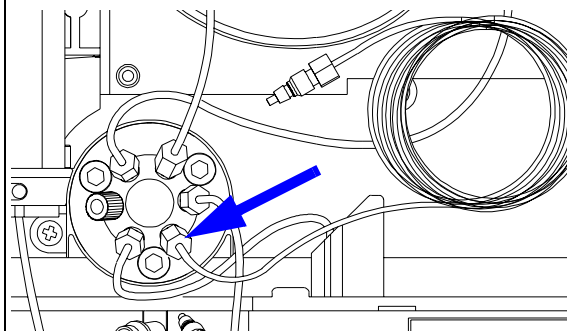
Before beginning this procedure:

- 1 Remove the front cover.

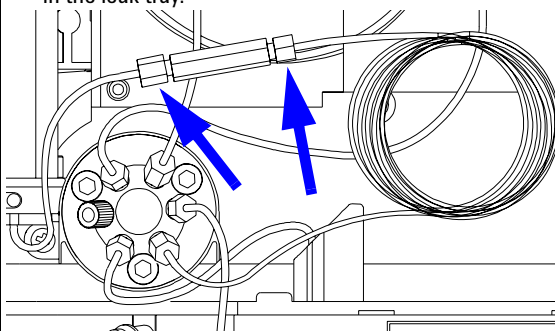
- 1 Disconnect the seat-capillary fitting from the injection valve (port 5).



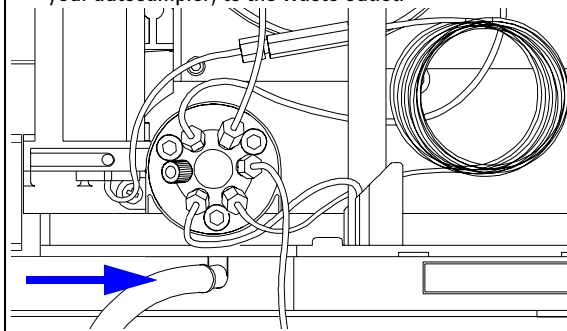
- 2 Install the 400 or 1400 µl seat extension loop (longer side) to the injection valve (port 5)



- 3 Install the union between the seat capillary and the other side of the seat extension loop. Store the extension loop in the leak tray.



- 4 Connect the waste tube (shipped with the start up kit of your autosampler) to the waste outlet.



On completion of this procedure:

- 1 Install the front cover.

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.

ChemStation Configuration

- 1 Select **More Injector** in the instrument function.
- 2 In the configuration menu change seat capillary to the value of the installed extended seat capillary (400 µl or 1400 µl) and press **OK**.

Control Module Configuration

- 1 Press **System > Configure** and select the Autosampler from the pull down menu.
- 2 Select the **volume** for your installed extended seat capillary (400 µl or 1400 µl) from the Seat Capillary pull down menu and press **Done**.

Instant Pilot

- 1 In the startup screen select **More > 1 Configure**.
- 2 Then select **Autosampler** and specify the **Volume** of the installed seat capillary and the **Multiple Draw Wait** time.

