

Agilent 210/218 - Isocratic Pump

Installation of the Stainless Steel Tubing Kit

Introduction

This technical note shows you how to install the Stainless Steel Tubing Kit for the isocratic pump of the Agilent 218 Solvent Delivery Module.

The tubing is pre-cut and bent to fit between components when placed on the mast in the correct location. Follow the directions in this note carefully when assembling and installing the mast kit, injection valve, prime/purge valve and 3-way pump head prime valves.



Delivery Checklist

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

The follow items are supplied with the Isocratic Pump Stainless Steel Tubing Kit, 1/16 in x 0.02 in i.d. (p/n 9910138000):

p/n	Description
1610141800	Nut, stainless steel, 10 – 32 x 1 /16 in, pack of 2. Use part number AL206085 when ordering replacements.
1610143200	Ferrule, stainless steel, 1 /16 in, pack of 2. Use part number AL286075 when ordering replacements.
3710064300	Tubing, pump head to pressure module, 1/16 in x 0.02 in i.d.
3710064700	Tubing, purge valve to manual injection valve, 1/16 in x 0.02 in i.d.
3710065000	Tubing, pressure module to purge valve, 1/16 in x 0.02 in i.d.

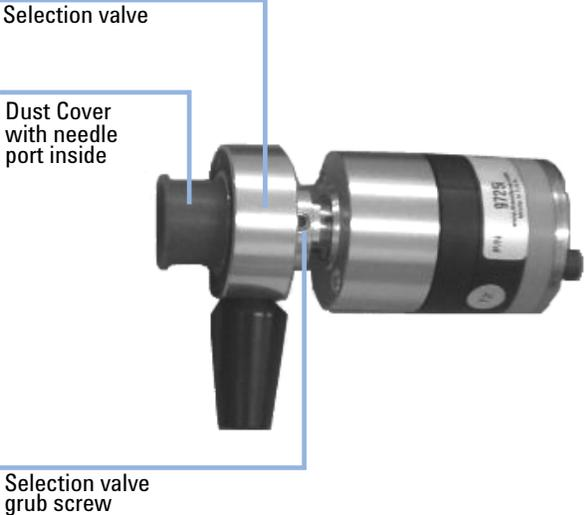
Installing the Mast Kit

The mast kit includes installation of the:

- Mast
- Manual injection valve
- Prime/purge valve
- 3-way pump head primve valve (If included with the pumphead that was purchased.)

Assembling the Manual Injection Valve and Bracket

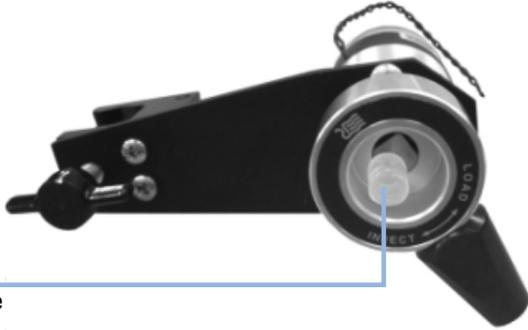
Tools required	Description
	Hex key, 2 mm

<p>1 Unpack the injection valve and attached black cable.</p>  <p>Selection valve</p> <p>Dust Cover with needle port inside</p> <p>Selection valve grub screw</p>	<p>2 Remove the red dust cover and gently pull out the needle port from the valve.</p>
<p>3 Remove the selection handle from the valve by loosening the two grub screws securing it to the injection valve body using the provided 2 mm hex key.</p>  <p>Flat side of metal post</p>	<p>4 Remove the two screws and injection valve bracket from the accessories bag that came with the injection valve.</p> <p>5 Attach the bracket to the injection valve with the two screws.</p>

Installing the Mast Kit

Assembling the Manual Injection Valve and Bracket

- 6** Replace the selection handle and tighten the two grub screws securing the handle using the 2 mm hex key. The grub screws should be flush with the flat side of the metal post.

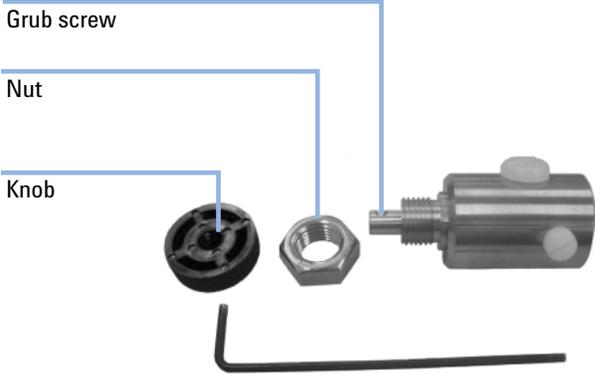


Needle
port

- 7** Replace the needle port in the valve.

Assembling the Prime/Purge Valve and Bracket

Tools required	Description
	Hex key, 2 mm

<p>1 Unpack the stainless steel purge valve and purge valve bracket.</p> 	<p>2 Remove the knob and nut from the purge valve. The knob is held in place with a grub screw which is loosened using a 2 mm hex key.</p> 
<p>3 Secure the valve to the valve bracket using the nut.</p> 	<p>4 Replace the knob on the valve and tighten the grub screw using the provided 2 mm hex key.</p> 

Assembling the 3-way Pump Head Prime Valve

Tools required

Description

Hex key, 2 mm

CAUTION

Overtightening the fittings may cause damage to the fittings.

→ Do not over-tighten the fittings.

→ Finger-tighten the nut only and use only the provided hex key to tighten the selection knob.

1 Unpack the stainless steel 3-way valve and bracket.



2 Using the 2 mm hex key, loosen the grub screw and slide off the selection knob.

Ridged washer

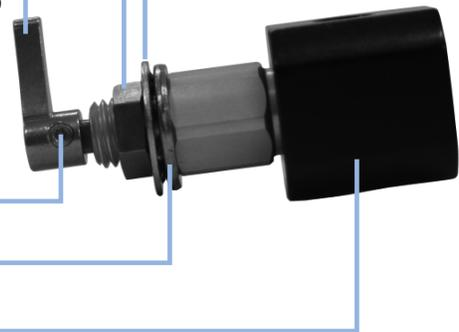
Nut

Selection knob

Grub screw

Flat washer

Valve body



3 Unscrew the nut and then remove the two washers from the 3-way valve.

4 Attach the bracket so that the three holes on the 3-way valve face away from the thumbscrew.

Flat side of metal post



Holes in 3-way valve body

5 Place the flat washer and then ridged washer on the 3-way valve. Screw the nut on to secure the two washers.

6 Using the clear tubing supplied in the pump head kit, install the shorter piece.

7 Slide the selection knob onto the 3-way valve.

NOTE

The engraved lines on the top of the selection knob should point to the holes in the black body of the 3-way valve. The grub screw should be flush with the flat side of the metal post on the 3-way valve.



8 Secure the selection knob to the 3-way valve using only the provided 2 mm hex key. Do not over-tighten.

Installing the Mast

Make sure the manual injection valve, prime/purge valve and 3-way pump head prime valve have a bracket attached before installing them onto the mast. See the corresponding procedures in this note for more information about the installation of the brackets for each valve.

NOTE

Perform the procedure only if a mast kit is not already installed.

- 1 Remove the mast, brackets and two pan-head screws from the Mast Kit Assembly.
- 2 Attach a mast clamp to the fittings at the front corner of the 210/218 pump or one at the lower front corner of the pump and one at the top front corner of the Agilent detector, if fitted, see [Figure 1](#) on page 8.

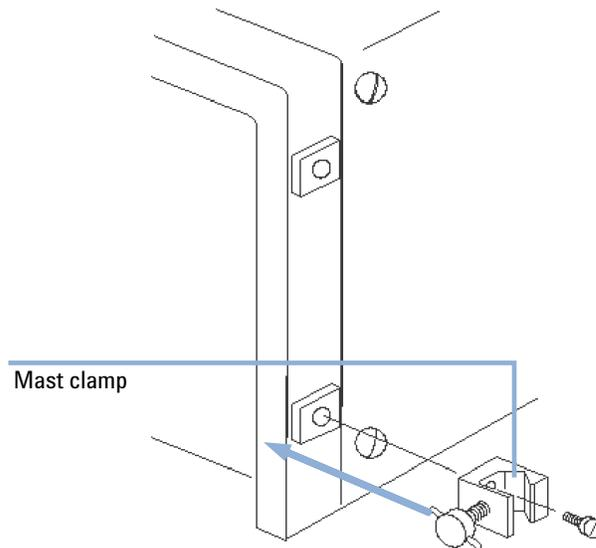


Figure 1 Mast clamp installation

- 3 Secure the stainless-steel mast within the clamps so the lower end of the mast is level with the bench.

- 4 Attach each valve to the mast in the heights described in [Table 1](#) on page 9 and shown [Figure 2](#) on page 9.

Table 1 Valve heights

Valve	Height
Injection Valve	285 mm
Prime/purge Valve	185 mm
3-way Valve	100 mm

NOTE

Valve heights are measured from the top of the bench to the bottom of the valve brackets. Tubing length and bends may vary slightly. Use the heights listed as a starting point and move the valves up or down to accommodate the tubing, if required. Avoid bending the tubing when installing it.

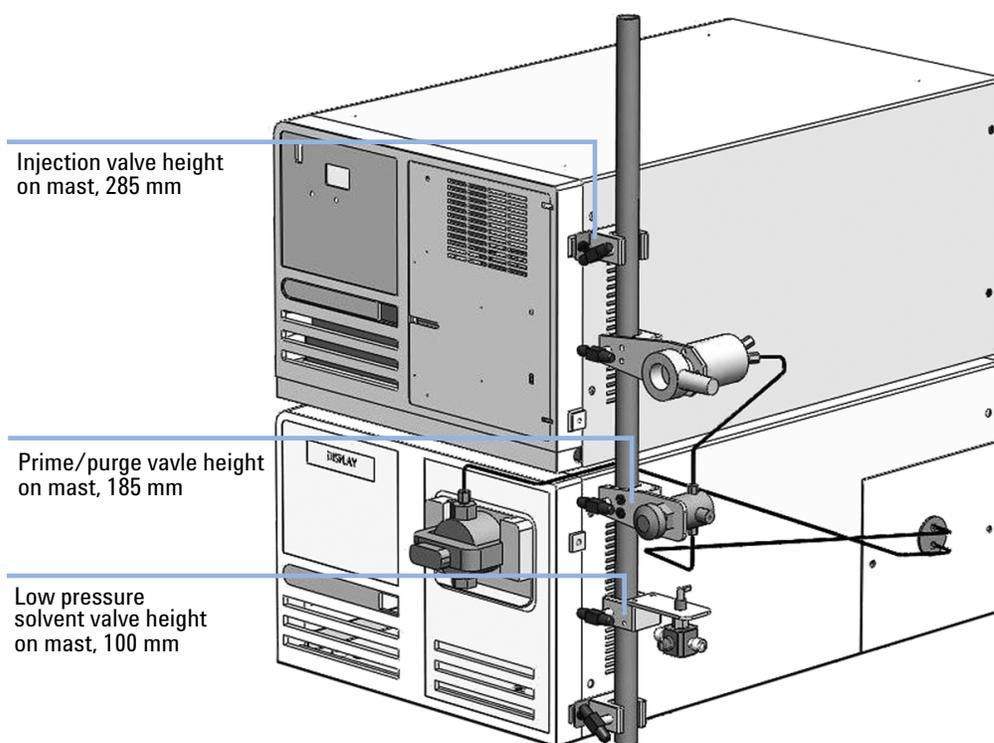


Figure 2 Component locations on the mast

Installing the Stainless Steel Tubing

Installing the Tubing

Tools required

Description

Wrench

Preparations

Install the mast kit on your modular HPLC system, see chapter *Installing the Mast Kit* for more information.

WARNING

When opening capillary or tube fittings, solvents may leak out.

The handling of toxic and hazardous solvents and reagents can carry health risks.

- Observe appropriate safety procedures (for example, wear goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the solvent vendor, especially when toxic or hazardous solvents are used.

CAUTION

Overtightening the fittings may cause damage to the fittings.

- Do not over-tighten the fittings.
- Finger-tighten the nut only and use only the provided hex key to tighten the selection knob.

NOTE

The stainless steel tubing is supplied pre-cut and bent to the appropriate size when used with an Agilent pump and mast kit. Use the nuts and ferrules supplied with each component when installing tubing. The two nuts and ferrules provided in this kit are spares.

- 1 If you are replacing tubing on an existing system, remove the old tubing first.

- Carefully route the tubings according to [Figure 3](#) on page 11. No additional bending of the tubings should be required.

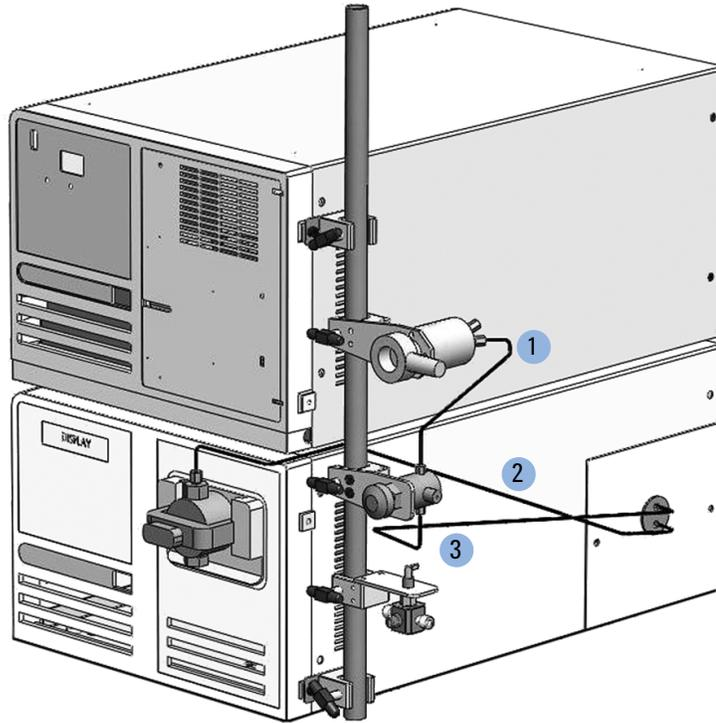
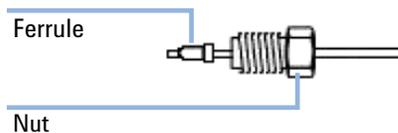


Figure 3 Installed tubing Kit

1	Tubing, purge valve to manual injection valve, 1/16 in x 0.02 in i.d. (p/n 3710064700)
2	Tubing, pump head to pressure module, 1/16 in x 0.02 in i.d. (p/n 3710064300)
3	Tubing, pressure module to purge valve, 1/16 in x 0.02 in i.d. (p/n 3710065000)

- Place the nut and then the ferrule on the tubing.



- Place the loosely assembled fittings into the valve ports and make sure the tubing is bottomed out inside the port.

NOTE

Ensure the appropriate fittings are used and insert the tubings fully into the correct components before tightening the fittings.

Installing the Stainless Steel Tubing

Installing the Tubing

- 5 Tighten the nut finger-tight and afterwards an additional $\frac{3}{4}$ of a turn with a wrench.

The manual injection valve is typically set up so that the:

- sample loop connects to ports 1 and 4
- tubing from the purge valve connects to port 2
- tubing to the column connects to port 3
- ports 5 and 6 are waste lines



8510249900

Part Number: 8510249900

Edition: 11/12
Printed in Germany

© Agilent Technologies, Inc 2012

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