

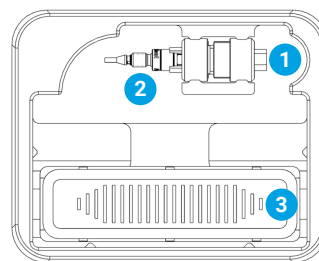
# Agilent InfinityLab Quick Change Inline Filter with Rigid Capillary

Engineered for efficiency and ease of use



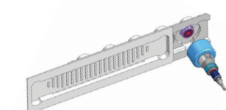
## Kit contents

InfinityLab Quick Change inline filter assembly

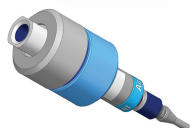


1. Filter housing
2. Quick Turn fitting
3. Filter disks in touchless packaging, 5/pk

## Procedure to connect the filter onto your system



1. Using the touchless packaging in which your filter disks are packed, insert a filter disk into the inline filter part with the blue screw nut.



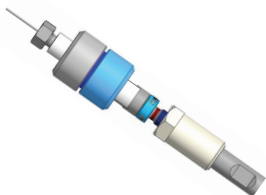
2. By hand, turn both parts of the filter housing at the same time until you hear a click. The click means that your inline filter is sealed up to 1,300 bar. **No tools are needed.**  
*Note: It is usual for the product to feel loose after the housing is closed, but this will not affect the seal.*



3. Disconnect the fitting connection at your column.



4. Connect the capillary onto the female end of the inline filter. Depending on the type of fitting you use, wrenches might be needed to tighten the fitting.



5. Turn on the flow until the eluent\* comes out of the inline filter at the male end of the filter. (This step is to remove the air out of the new filter in order to avoid air bubbles entering into the column.)
6. Turn the inline filter onto your column by hand tightening.

**Limits:** Depending on the users, different pressure limit between 800-1,300 bar can be reached by hand tightening. If higher pressure limits is required, please tighten the fitting with the wrench by 1/4-turn.

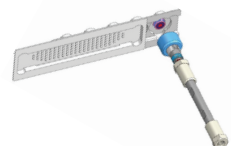
## Procedure to replace the filter disk



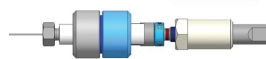
1. By hand, turn both nuts of the filter housing anticlockwise to open the filter. You can also hold one nut and turn the other.



2. Remove the used filter disk.



3. Insert a new filter disk into the filter housing part with the blue nut using the touchless packaging in which your filter disk is packed.



4. By hand, turn both parts of the filter housing at the same time until you hear a click.

5. If the filter disconnected from the column at step 1, turn the inline filter onto your column again by hand tightening.

**Warning:** Don't disconnect the Quick Turn fitting from the column when the system pressure is on because it might detach the fitting from the filter. Please check the spring-loaded function of the Quick Turn fitting before reconnecting.

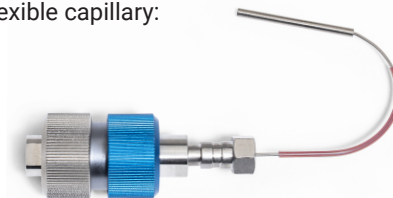
\* Don't use corrosive solvent such as strong halogenated solvents or strong acidic or basic solutions for step 5 because the vapor from the drop of the eluent might cause corrosion of the parts inside the fitting.

## Part numbers

Inline Filter Assemblies	p/n
InfinityLab Quick Change inline filter assembly with <b>rigid capillary</b> for UHPLC (incl. 5 filter disks 2.1 mm id, 0.2 µm porosity)	5067-1607
InfinityLab Quick Change inline filter assembly with <b>rigid capillary</b> for HPLC (incl. 5 filter disks 4.6 mm id, 0.5 µm porosity)	5067-1606
InfinityLab Quick Change inline filter assembly for UHPLC (incl. 5 filter disks 2.1 mm ID, 0.2 µm porosity), with 90 mm <b>flexible capillary</b>	5067-1603
InfinityLab Quick Change inline filter assembly for HPLC (incl. 5 filter disks 4.6 mm ID, 0.5 µm porosity), with 90 mm <b>flexible capillary</b>	5067-1602
Filter Disks	p/n
Filter disks 2.1 mm ID, 0.2 µm porosity, 5/pk, in touchless packaging	5067-1610
Filter disks 2.1 mm ID, 0.5 µm porosity, 5/pk, in touchless packaging	5067-1611
Filter disks 4.6 mm ID, 0.2 µm porosity, 5/pk, in touchless packaging	5067-1612
Filter disks 4.6 mm ID, 0.5 µm porosity, 5/pk, in touchless packaging	5067-1613
Filter disks 4.6 mm ID, 2.0 µm porosity, 5/pk, in touchless packaging	5067-1614
Replacement Capillaries for Flexible Capillary Version	p/n
Capillary stainless steel 0.12 mm id, 90 mm length, 2x extra-long fittings, preswaged on one end, nonswaged on the other end, for inline filter for UHPLC	5023-3344
Capillary stainless steel 0.17 mm id, 90 mm length, 2x extra-long fittings, preswaged on one end, nonswaged on the other end, for inline filter for HPLC	5023-3343
Quick Turn Fitting**	p/n
Quick Turn Fitting	5067-5966
Ferrule for Quick Turn fitting	5043-0924

\*\* go to package insert (publication number 01200-90115) to see instruction for replacement of the ferrule / Quick Turn fitting

For connections of the inline filter onto the Autosampler or pump head, you could use the InfinityLab Quick Change inline filter with flexible capillary:



## Maintenance

It is recommended to monitor the backpressure of your routine applications. Replace the filter disk if backpressure is 10% above the regular value. A clogged filter disk contributes to carryover and retention time shifts.

DE72893907

[www.agilent.com](http://www.agilent.com)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2023

Published in US, April 1, 2023

Part Number 5994-6005EN