

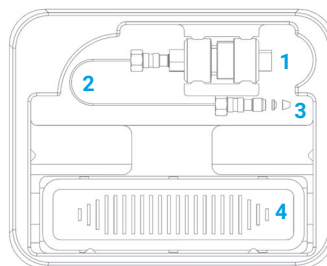
Agilent InfinityLab Quick Change Inline Filter

Engineered for efficiency and ease of use



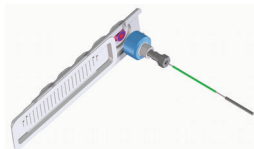
Kit contents

InfinityLab Quick Change inline filter assembly

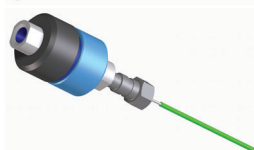


1. Filter housing (two parts)
2. Capillary, SST, 90 mm length
3. Loose fitting for the nonswaged end of the capillary
4. Filter discs in touchless packaging, 5/pk

Procedure to connect the inline filter onto your system



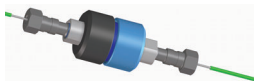
1. Using the touchless packaging in which your filter discs are packed, insert a filter disc into the inline filter part with the blue screw nut.



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. Using the loose fitting that is included with your product, screw your inline filter into port 6* of the autosampler injection valve.
4. Recommended: Flush the inline filter with your mobile phase shortly before the next step.

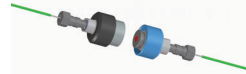


5. Use your system capillary to connect the other end of the inline filter to your system.

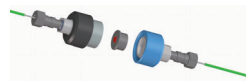
* Position 6 of the injection valve of autosampler is recommended, but you can connect the inline filter into another position if you prefer. To minimize sample dispersion, connect the inline filter to a position outside of your sample flow path, such as onto the outlet valve at the pump head.

Note: If the dimension of the capillary that is included in the assembly does not fit into your system configuration, you can replace it with other capillaries. You can also use other fittings, such as InfinityLab Quick Connect or Quick Turn fittings, for your capillary connections.

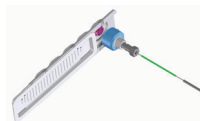
Procedure to replace the filter disc



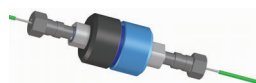
1. By hand, turn both nuts anticlockwise to open the filter housing. You can also hold one nut and turn the other. There is **no need** to disconnect the filter from your system.



2. Remove the used filter disc.



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



4. 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.

Part numbers

Inline Filter Assemblies	p/n
InfinityLab Quick Change inline filter assembly for UHPLC (incl. 5 filter discs 2.1 mm id, 0.2 µm porosity), with 0.12 mm × 90 mm flexible capillary	5067-1603
InfinityLab Quick Change inline filter assembly for HPLC (incl. 5 filter discs 4.6 mm id, 0.5 µm porosity), with 0.17 mm × 90 mm flexible capillary	5067-1602
Filter Discs	
Filter discs 2.1 mm id, 0.2 µm porosity, 5/pk	5067-1610
Filter discs 2.1 mm id, 0.5 µm porosity, 5/pk	5067-1611
Filter discs 4.6 mm id, 0.2 µm porosity, 5/pk	5067-1612
Filter discs 4.6 mm id, 0.5 µm porosity, 5/pk	5067-1613
Filter discs 4.6 mm id, 2.0 µm porosity, 5/pk	5067-1614
Replacement Capillaries	
Capillary stainless steel 0.12 mm id, 90 mm length, 2x extralong 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 extralong fittings, preswaged on one end, nonswaged on the other end, for inline filter for HPLC	5023-3343

Maintenance

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

DE.5105208333

www.agilent.com

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020
Published in US, December 1, 2020
Part Number 5994-2779EN