

Agilent Captiva Filter Vials

Instructions for use

Filtrating samples before analysis can generate longer column lifetime, less downtime, and improved sample integrity. Agilent Captiva filter vials do the filtering for you, giving you one less step in your workflow.

Benefits include:

- **Greater convenience:** Filtration takes place inside the vial, reducing steps and tools.
- **Cleaner samples:** Fewer touchpoints in your sample journey mean fewer chances for contamination.
- **Higher-quality data:** Even small amounts of particulates can clog your column inlet, causing high backpressure, retention-time shift, and resolution loss.

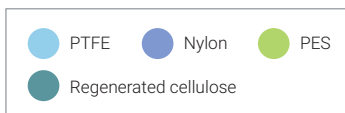
Reduce the steps in your GC or LC workflow with Agilent Captiva filter vials—now with preslit septa options



Color-coded for your convenience

Pore size identified by septum color

- 0.2 µm: white
- 0.45 µm: red



The Agilent Captiva filtration portfolio is designed to ensure you can choose the best product for your lab. Whether you prefer the convenience of a filter vial, the value of an econofilter, or the certified performance of a premium syringe filter, Captiva filtration has the solution for your particulate removal needs.

Learn more at www.agilent.com/chem/filtration

Agilent Captiva filter vials provide a convenient way to filter samples before LC or GC analysis. They replace the combination of syringe filters, syringes, autosampler vials, septa, and caps with a single unit.

Easy as 1-2-3



1. Fill:
Use a pipette to add sample to the fill line.



2. Cover:
Twist gently to ensure a secure seal.



3. Plunge:
Press the plunger slowly over ~three seconds to filter the sample

When using in an autosampler, adjust draw position to 10 mm, and uncheck bottom sensing.

Ordering Information

Description	Part Number (Nonslit Septa)	Part Number (Preslit Septa)
0.45 µm PTFE filter vial, 100/pack	5191-5933	5610-2122
0.20 µm PTFE filter vial, 100/pack	5191-5934	5610-2123
0.45 µm Nylon filter vial, 100/pack	5191-5935	5610-2118
0.20 µm Nylon filter vial, 100/pack	5191-5936	5610-2119
0.45 µm RC filter vial, 100/pack	5191-5939	5610-2124
0.20 µm RC filter vial, 100/pack	5191-5940	5610-2125
0.45 µm PES filter vial, 100/pack	5191-5941	5610-2120
0.20 µm PES filter vial, 100/pack	5191-5942	5610-2121
Vial closure tool	5191-5943	

Learn more:

www.agilent.com/chem/filtration

Buy online:

www.agilent.com/chem/store

Find an Agilent or authorized distributor:

www.agilent.com/chem/contactus

U.S. and Canada

1-800-227-9770

agilent_inquiries@agilent.com

Europe

info_agilent@agilent.com

Asia Pacific

inquiry_lsca@agilent.com

India

india_lsca_marketing@agilent.com

This information is subject to change without notice.

© Agilent Technologies, Inc. 2022
Published in the USA, July 15, 2022
Publication number 5994-0814EN



5191-5907

Filter	PES	PTFE	RC	NY
Solvents				
Acetone	-	**	**	**
Acetonitrile	-	**	**	N/A
Benzene	-	*	**	**
Benzyl Alcohol	-	**	**	**
<i>n</i> -Butyl Acetate	-	**	**	**
<i>n</i> -Butanol	*	**	**	**
Carbon Tetrachloride	-	*	**	**
Chloroform	-	*	**	**
Cyclohexane	-	**	**	**
Diethylacetamide	-	-	**	**
Diethyl Ether	-	**	**	**
Dimethyl Formamide	-	**	*	*
Dioxane	-	**	**	**
Ethanol, 98 %	**	**	**	**
Ethyl Acetate	-	**	**	**
Ethylene Glycol	**	**	**	**
Formamide	**	**	**	**
Gasoline	*	**	**	**
Glycerin	**	**	**	**
<i>n</i> -Heptane	**	**	**	**
<i>n</i> -Hexane	**	**	**	**
Isopropanol	**	**	**	**
Isopropyl Acetate	-	**	**	**
Methanol, 30 %	**	**	**	**
Methanol, 98 %	*	**	**	**
Methyl Acetate	-	**	**	**
Methylene Chloride	-	**	**	**
Methyl Ethyl Ketone	-	**	**	**
Methyl Isobutyl Ketone	-	**	**	**
Monochlorobenzene	-	**	**	**
Pyridine	-	**	**	**
Tetrahydrofuran	-	*	**	**
Toluene	-	*	**	**
Trichloroethane	N/A	**	N/A	N/A
Xylene	-	*	**	**
Acids				
Acetic Acid, 25%	*	**	**	-
Acetic Acid, 80%	N/A	**	**	-
Hydrochloric Acid, 20 %	**	**	-	-
Hydrofluoric Acid, 25 %	*	**	*	-
Nitric Acid, 25 %	*	*	-	-

- ** Compatible
- * Limited compatibility
- Not compatible
- N/A Not analyzed

** Chemical compatibilities can be influenced by various factors. Therefore, we recommend that you confirm compatibility with the liquid you want to filter by performing a trial filtration run before you start your actual filtration.

Filter	PES	PTFE	RC	NY
Phosphoric Acid, 1 %	**	**	-	-
Sulfuric Acid, 25 %	*	**	*	-
Trichloroacetic Acid, 10%	N/A	**	**	-
Bases				
Ammonium Hydroxide, 25 %	*	**	*	**
Sodium Hydroxide, 1 N	**	**	*	**
Aqueous Solutions				
Formalin, 30 %	*	**	*	**
Hydrogen Peroxide, 30 %	**	**	-	-
Sodium Hypochlorite, 5 %	**	**	-	-
pH Range				
pH 1 to 14	-	**	-	-
pH 1 to 13	**	**	-	-
pH 3 to 14	*	**	*	**
pH 3 to 12	**	**	**	**
pH 4 to 8	**	**	**	**