

Minisart® RC Syringe Filters for HPLC Sample Preparation

RC versus PVDF used with Methanol and Acetonitrile

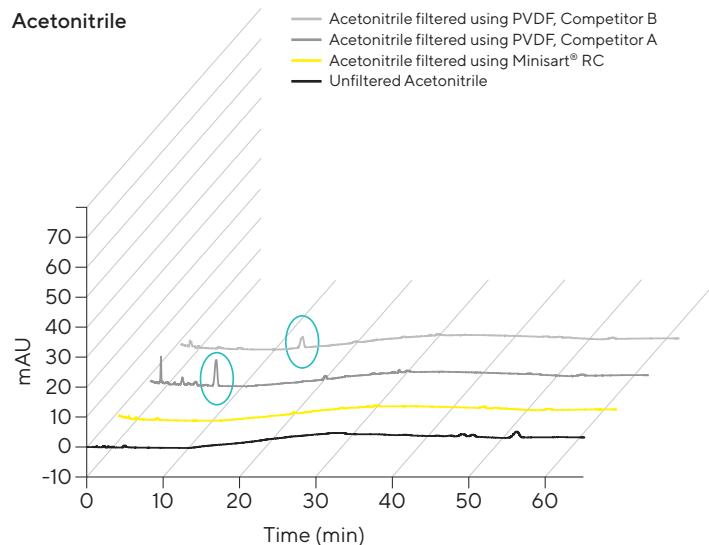
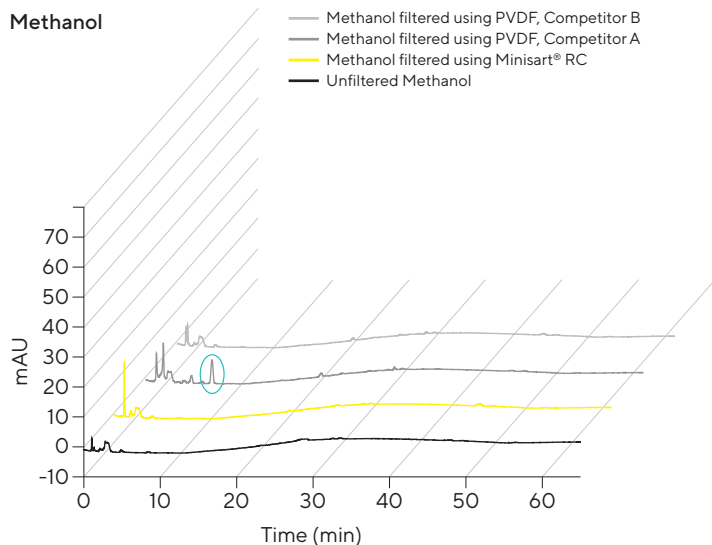
Comparison with PVDF

PVDF is a membrane material often used for HPLC sample preparation. The fluorochemical process needed to produce this raw material can be harmful to the environment. Moreover, PVDF requires an additional coating step with methacrylates to make its membrane surface hydrophilic. In addition, unlike RC syringe filters, those with PVDF membranes are not suitable for DMSO and other amides, or for ketones, esters, and ethers.

Conclusion

Minisart® RC is just as suitable, and often even better, for HPLC sample preparation than PVDF. The broader chemical compatibility of Minisart® RC enables usage with DMSO and for an even wider range of applications.





HPLC Procedure

Column C18: 250 × 4.6 mm, Flow Rate: 1 mL/min, Wavelength: 210 nm

HPLC Injection Volume: 20 µl, Analysis Time: 65 min, Temperature: 40 °C, Mobile Phases: A) Acetonitrile | B) Water, Gradient: Hold 60% A for 10 min | 60% to 95% A in 20 min | 95% to 100% A in 35 min

Reliable and Ultrapure

Clean & Green

Are you using high-purity filter paper in your lab? Our regenerated cellulose (RC) membrane incorporated in Minisart® RC syringe filters is just as pure, but allpurpose and more convenient.

Our RC polymer is naturally hydrophilic so no additional coating step is necessary. Try our Clean and Green solution to prepare your samples.

RC is produced using renewable raw materials in an eco-friendly process, which uses biodegradable components and recycles all solvents employed during manufacture.

Ordering Information

Ø mm	Pore Size	Sterile*	Qty Pk	Order No.
Minisart® RC (Regenerated Cellulose + PP)				
25 mm	0.2 µm	Yes	50	17764-----ACK
25 mm	0.2 µm	No	50	17764-----K
25 mm	0.2 µm	No	200	17764-----S
25 mm	0.2 µm	No	500	17764-----Q
25 mm	0.45 µm	No	50	17765-----K
25 mm	0.45 µm	No	200	17765-----S
25 mm	0.45 µm	No	500	17765-----Q

Ø mm	Pore Size	Sterile*	Qty Pk	Order No.
15 mm	0.2 µm	Yes	50	17761-----ACK
15 mm	0.2 µm	No	50	17761-----K
15 mm	0.2 µm	No	500	17761-----Q
15 mm	0.45 µm	No	50	17762-----K
15 mm	0.45 µm	No	500	17762-----Q
4 mm	0.2 µm	No	50	17821-----K
4 mm	0.2 µm	No	500	17821-----Q
4 mm	0.45 µm	No	50	17822-----K
4 mm	0.45 µm	No	500	17822-----Q

* Sterile Minisart® are individually packaged and have been sterilized by ethylene oxide (EO).
 Not presterilized Minisart® can be sterilized by autoclaving at 121 °C for 30 min | or by using EO.

Would You Like to Use Other Membranes?

Please refer to Minisart® SRP (hydrophobic PTFE for highest chemical compatibility).

Do You Need Minisart® with Pre-filters for Filtration of Highly Particle Laden Samples?

Please refer to Minisart® NY Plus which incorporate a glass fiber pre-filter.

Are You Looking for CE-certified Minisart®?

Please request CE-certified article numbers for Minisart® NML, Minisart® HY and Minisart® SRP. Sartorius offers a full range of syringe filters dedicated for various filtration applications.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com