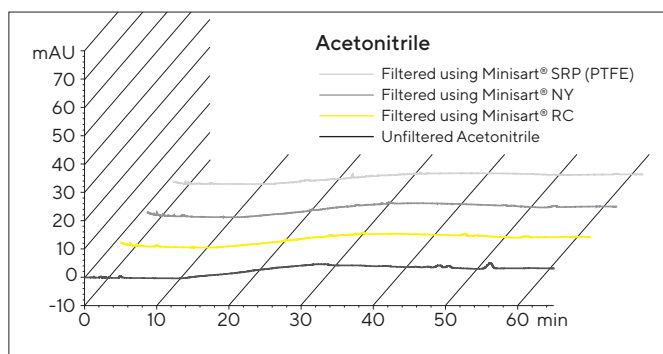
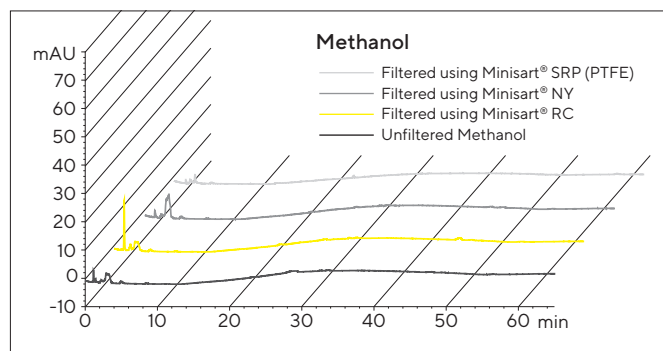


## Minisart® Syringe Filters for HPLC Sample Preparation

Sartorius Minisart® with a Polypropylene (PP) Housing are Available in a Choice of Membranes:  
Hydrophilic Regenerated Cellulose (RC), Nylon (PA) or Hydrophobic PTFE

The purpose of using syringe filters for sample preparation prior to analysis is to protect analytical equipment and columns. Sartorius Minisart® PP filters reliably remove particles from your samples without adding unwanted extractables or leachables.



### 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

Our regenerated cellulose (RC) membrane used in Minisart® RC, our nylon (polyamide) membrane incorporated in Minisart® NY and our coating-free hydrophobic PTFE membrane used in Minisart® SRP are exceptionally pure compared with other common filter membranes.

## Ordering Information

Ø mm	Pore Size	Sterile*	Qty/Pk	Order No.
<b>Minisart® RC (Regenerated Cellulose + PP)</b>				
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
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
<b>Minisart® NY (Nylon   Polyamide + PP)</b>				
25 mm	0.2 µm	Yes	50	17845-----ACK
25 mm	0.2 µm	No	500	17845-----Q
25 mm	0.45 µm	Yes	50	17846-----ACK
25 mm	0.45 µm	No	500	17846-----Q
15 mm	0.2 µm	No	50	1776B-----K
15 mm	0.2 µm	No	500	1776B-----Q
15 mm	0.45 µm	No	50	1776C-----K
15 mm	0.45 µm	No	500	1776C-----Q
25 mm	GF1 prefilter   0.2 µm	No	50	1784B-----K
25 mm	GF1 prefilter   0.2 µm	No	500	1784B-----Q
25 mm	GF1 prefilter   0.45 µm	No	50	1784C-----K
25 mm	GF1 prefilter   0.45 µm	No	500	1784C-----Q

Ø mm	Pore Size	Sterile*	Qty/Pk	Order No.
<b>Minisart® SRP (PTFE + PP)</b>				
25 mm	0.2 µm	Yes	50	S7575-----FXOSK
25 mm	0.2 µm	No	50	17575-----K
25 mm	0.2 µm	No	200	17575-----S
25 mm	0.2 µm	No	500	17575-----Q
25 mm	0.45 µm	No	50	17576-----K
25 mm	0.45 µm	No	200	17576-----S
25 mm	0.45 µm	No	500	17576-----Q
15 mm	0.2 µm	No	50	17558-----K
15 mm	0.2 µm	No	500	17558-----Q
15 mm	0.2 µm	Yes	50	17573-----ACK
15 mm	0.2 µm	No	50	17573-----K
15 mm	0.2 µm	No	500	17573-----Q
15 mm	0.45 µm	No	50	17559-----K
15 mm	0.45 µm	No	500	17559-----Q
15 mm	0.45 µm	No	50	17574-----K
15 mm	0.45 µm	No	500	17574-----Q
4 mm	0.2	No	500	17844-----Q
4 mm	0.45 µm	No	50	17820-----K
4 mm	0.45 µm	No	500	17820-----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. Non-sterile RC Minisart® types are optimized for sample preparation and are not suitable for sterile filtration according to the bacteria challenge test.

<sup>1</sup> GF prefilter: Ultrapure glass fiber with 0.7 µm particle retention on top of the nylon membrane.

## Would You Like to Use Other Membranes?

- Minisart® High Flow with polyethersulfone (PES) is optimal for filtration of additives and cell culture buffers. Due to the asymmetric membrane structure, the PES surface almost behaves like a pre-filter and enables high flow rates
- Minisart® NML (SFCA - Surfactant-free Cellulose Acetate) for sterile filtration and clarification of aqueous solutions

Sartorius offers a full range of syringe filters dedicated for various filtration applications.