



**New!**  
 Minisart® RC now with  
**lot-specific\* HPLC-Certificates!**  
[www.sartorius.com/lab-filtration](http://www.sartorius.com/lab-filtration)  
**Download now!**  
 \*for lot numbers > 70194103

# HPLC-Certificate

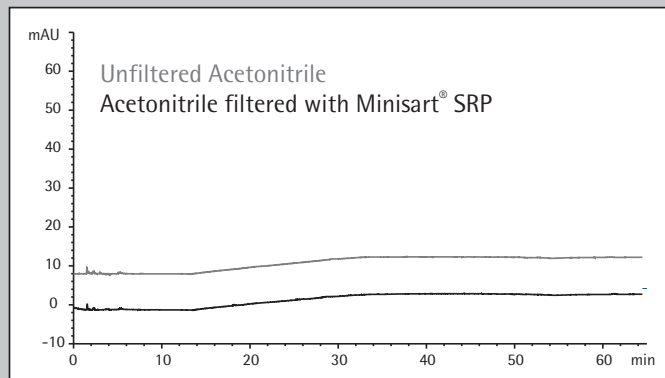
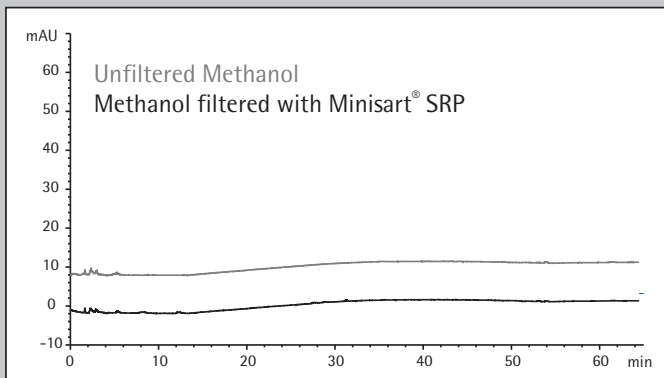
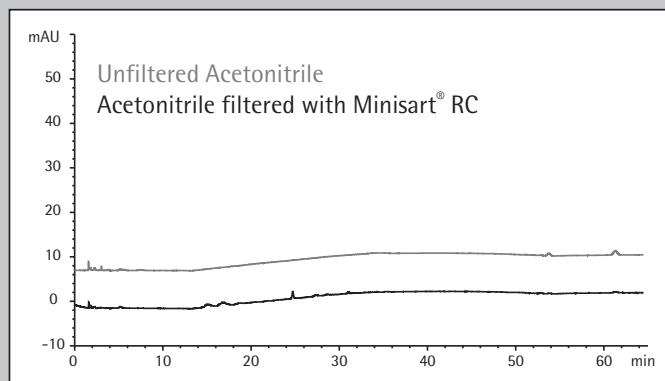
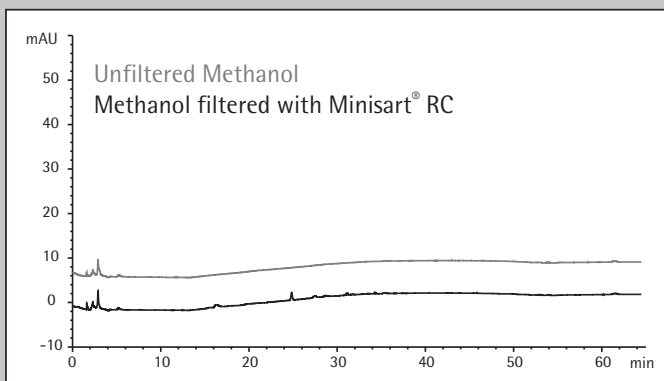
Minisart® Syringe Filters for Sample Preparation – reliable and ultrapure

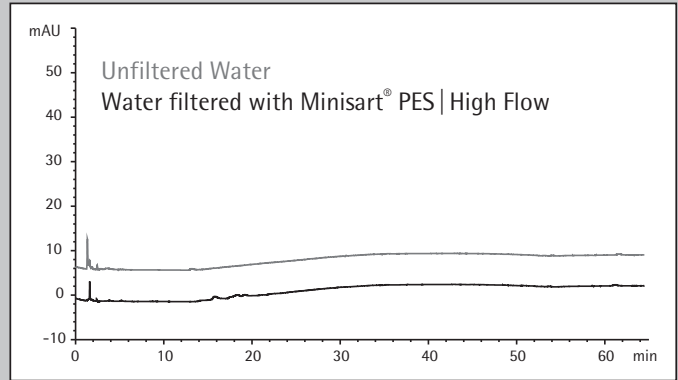
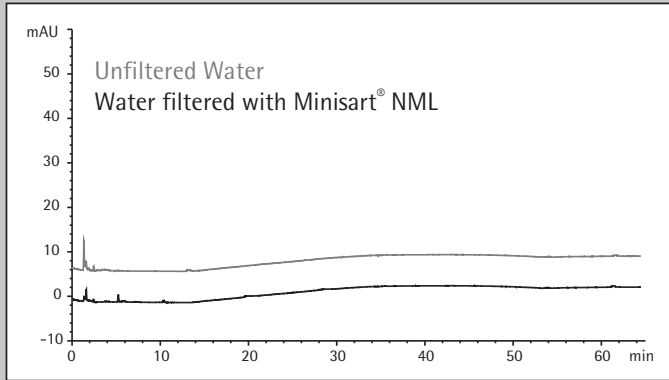
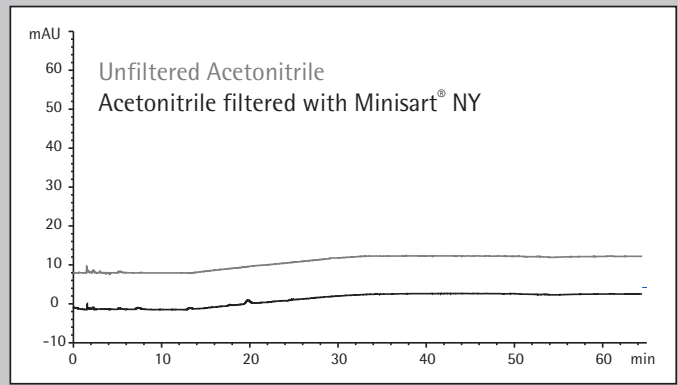
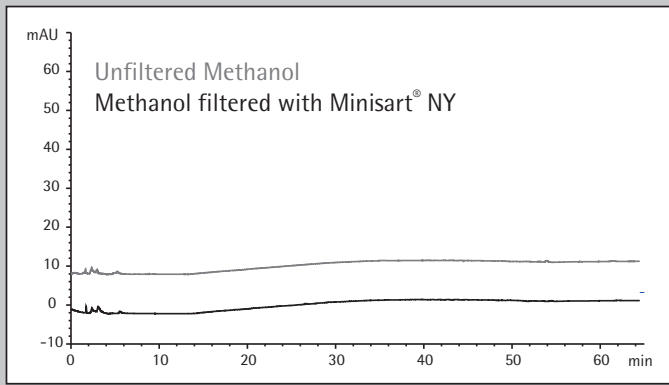
## Sartorius Manufacturer Statement

The following article numbers are tested and released for HPLC.  
 Rinsed with methanol, acetonitrile, or water they show extremely low leachable and extractable levels.

Minisart® RC with <b>Methanol &amp; Acetonitrile</b>	17761	17762	17764	17765	17821	17822		
Minisart® SRP (PTFE) with <b>Methanol &amp; Acetonitrile</b>	17558	17559	17573	17574	17575	17576	17820	17844
Minisart® NY & NY Plus with <b>Methanol &amp; Acetonitrile</b>	17845	17846	1776B	1776C	1784B	1784C		
Minisart® PES   High Flow   NML (CA) with <b>Water*</b>	1776D	16532	16533	16537	16541	16553	16534	16555
	16592	17593	17594	17597	17598			

\*Minisart® PES (1776D) is suitable for up to 30% methanol





**HPLC Procedure**

Column:

C18: 250 × 4.6 mm

Flow Rate: 1 ml/min

Wavelength: 220 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

Certified by

Dr. Hartmut Hennig,  
Site Manager QA Filtration Products

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen

Phone +49.551.308.0

Fax +49.551.308.3289

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

Specifications subject to change  
without notice.

Printed and copyrighted by  
Sartorius Stedim Biotech GmbH  
Publication No.: SL-6192-e170204  
Order No.: 85037-547-44  
Ver. 02 | 2017