

HPLC Purity Certificate

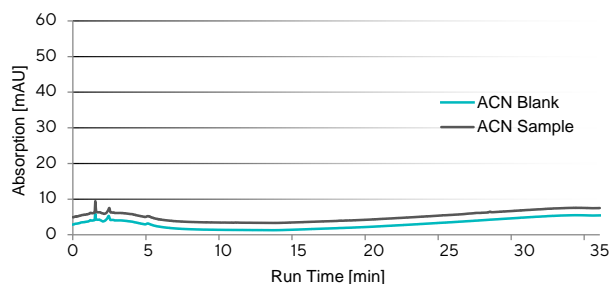
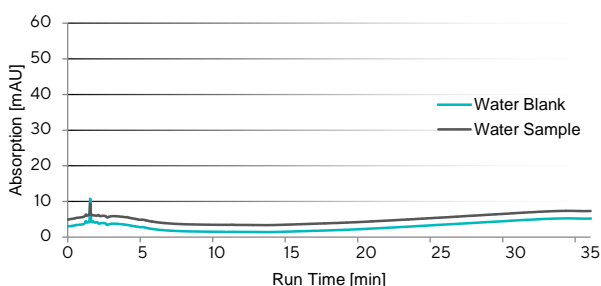
Minisart® RC

Order no.: 17764 ----- Q

Lot no.: 220876103

This is to certify that the product described above has been analyzed for extractables by testing the incorporated membrane by HPLC-UV|Vis.

Chromatograms of the HPLC-UV|Vis analyses of the extracts



HPLC Method

Column: C18, 250 x 4.6 mm, Column Temperature: 40°C, Injection Volume: 20 µL, Flow Rate: 1 mL/min, Detection Wavelength: $\lambda = 220$ nm, Analysis Time: 35 min

Mobile Phases: A) Acetonitrile, B) Water

Gradient: 60% A for 10 min, 60% to 100% A in 20 min, 100% A for 5 min

Filter device eluted with 0.5 mL extraction solvent per cm² membrane area (surface-area to volume ratio: 2 cm²/mL).

For regenerated cellulose (RC), the extraction strength of the solvent acetonitrile has been tested to be comparable to methanol.

2022-08-10

Date

Dr. Anna Vreemann
Site Quality Manager

Manufactured by Sartorius Stedim Biotech GmbH
37070 Goettingen, Germany
Phone: +49.551.308.0
Fax: +49.551.308.3289
www.sartorius.com