

Vivaclear

Centrifugal Filters for Rapid Clarification of Low Volume Samples



Benefits

- Convenient filtration of low volume samples
- Superior flow rates ensure rapid clarification
- High sample recoveries with minimal volume loss
- Reproducible performance in multiple applications

Product Overview

Vivaclear centrifugal filters are disposable microfiltration devices for low volume sample filtration or clarification. Specifically designed for low sample volumes in the range of 100 to 500 μL , they feature a very low hold up volume for minimal sample loss.

The high flux polyethersulfone membrane ensures fast and reliable clarification, pre-filtration or particle removal from multiple sample types.

Product information

Vivaclear extends the familiarity of centrifugal ultrafiltration to microfiltration applications, including clarification, pre-filtration and particle removal.

Simply fill the device with the sample to be filtered and centrifuge at up to 2,000 g until the entire sample has passed through the membrane. Filtration typically takes only 2 - 10 minutes and the sample is ready to use immediately, or can be stored by sealing the filtrate tube with the integrated cap.

Typical Applications

- Clarification of samples prior to IEX purification with Vivapure® Mini H
- Sample clarification prior to concentration with Vivaspin® 500
- Pre-filtration of plasma and serum
- Cell culture or lysate clarification
- Particle or precipitate removal

Technical Specifications

Operating conditions	
Rotor type	Fixed angle
Rotor angle	40 - 45°
Maximum RCF	14,000 g

Dimensions	
Total length	43 mm
Diameter, filtrate tube	11 mm
Active membrane area	0.34 cm ²
Pore size	0.8 µm
Hold-up volume, membrane and support	<5 µL

Materials of construction	
Body	Polypropylene
Membrane	Polyethersulfone
Filtrate tube	Polypropylene

Ordering Information


Vivaclear	Quantity	Order No.
0.8 µm PES	100	VK01P042

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 contacts, visit
www.sartorius.com