



Vivaclear Centrifugal Filters

For in vitro use only



Technical specifications

Rotor	40–45° Fixed angle rotor 500 µl
-------	---------------------------------

Pore size	0.8 µm
-----------	--------

Dimensions

Total length	43 mm
--------------	-------

Filtrate collection tube diameter	11 mm
-----------------------------------	-------

Active membrane area	0.34 cm ²
----------------------	----------------------

Hold-up volume, membrane plus support	< 5 µl
---------------------------------------	--------

Maximum RCF	14,000 × g
-------------	------------

Materials of construction

Body	Polypropylene
------	---------------

Membrane	Polyethersulphone
----------	-------------------

Filtrate collection tube	Polypropylene
--------------------------	---------------

Ordering information

	Pack size	Cat. No
Vivaclear Mini 0.8 µm PES	100	VK01P042

Description

Vivaclear centrifugal filters are disposable microfiltration devices for the fast and reliable clarification | filtration of biological samples in the range 100 µl to 500 µl. They can be used in fixed angle rotors accepting 2.2 ml centrifuge tubes.

Product Features

- High-flux Polyethersulphone membrane
- 0.8 µm pore size
- Low hold up volume (<5 µl)
- Fast and reproducible performance

Applications

- Clarification of samples before loading onto Vivapure protein purification spin columns
- Removal of particles and particulates
- Filtration of plasma and serum
- Filtration of cells or cell debris

Operation

1. Fill device with up to 500 µl sample to be clarified.
2. Centrifuge at up to 2,000 × g for 2–10 minutes, or until sample has passed through the filter.
3. Remove and discard the filter.
4. The clarified filtrate can be used immediately, or the tube can be closed for short-term storage.

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

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.91.3586102
Japan +81.3.3740.5407

Specifications subject to change without notice.

Printed and copyrighted by
Sartorius Stedim Biotech GmbH · W/sart-294 · G
Publication No.: SLU2010-e08035
Order No.: 85030-519-55