

# Sartolab<sup>®</sup> RF 50

Vacuum Filtration Unit  
for Volumes up to 50 mL

## User Benefits

- Highest flow rates with a large filtration surface
- No loss of protein with a low protein affinity membrane
- Low hold-up volume due to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit (no additional transfer of liquid necessary for storage)
- Can be used as a stand-alone system or with the Sartolab<sup>®</sup> Multistation, which allows parallel filtration of up to 6 samples

## Product Information

Sartolab<sup>®</sup> RF 50 vacuum filtration units are single-use units for the filtration of sample volumes up to 50 mL. They are available with either a 0.22 or 0.45 µm polyethersulfone membrane.

Both membranes provide fast flow rates and ensure low protein binding and extractables. The 0.22 µm version is ideal for sterile filtration of cell cultures, buffers, and aqueous solutions, while the 0.45 µm version is ideal for clarification.

Sartolab<sup>®</sup> RF 50 are available either individually packed and sterile or in bulk non-sterile.



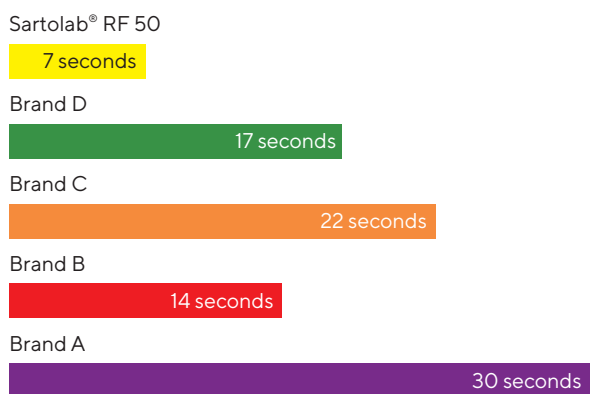
The Sartolab® RF 50 vacuum filtration unit comprises the following:

- a funnel, with dust cover
- a 50 mL conical tube, with graduation and writing field
- a tubing connector for vacuum connection (sterile versions only)
- a screw cap for storage of your filtrate (individually wrapped for the sterile versions and bulk wrapped for the non-sterile versions)

The design of the yellow adapter connecting the funnel to the conical tube ensures a vacuum-tight seal and enables the filtration unit to be used on the Sartolab® Multistation for filtration of up to 6 samples in parallel with one vacuum source. However, Sartolab® RF 50 can also be used alone by directly connecting the unit's tubing connector to your vacuum source.

Sartolab® RF 50 is the fastest 50 mL vacuum filtration unit compared to 4 of the leading competitor's units.

Comparison of filtration times for 50 mL cell media + 10% FBS in five 0.22 µm 50 mL vacuum filtration units



## Materials

Funnel with dust cover	Polystyrene (PS)
Membrane filter	0.22 µm asymmetric polyethersulfone (Cat. No. 180E01) 0.45 µm polyethersulfone (Cat. No. 180F01)
Funnel adapter	High Density Polyethylene (HDPE)
Tubing connector for vacuum connection	High Density Polyethylene (HDPE)
Conical tube	Polypropylene
Conical tube cap	High Density Polyethylene (HDPE)

## Specifications

Membrane diameter	58.5 mm
Effective filtration area	21 cm <sup>2</sup>
Hold-up volume	1.2 mL
Filtration capacity	50 mL
Size of the 50 mL conical tube	Height: 115.5 mm External/internal diameter: 29.5 mm/27.48 mm
Autoclavable	No
Sterilization method	E-Beam (beta) irradiation
Storage temperature of the conical tube	-80° C to max 100° C
Packaging	Single-packaged, sterile
Operating pressure	-350 to -750 mbar

## Ordering Information


Cat. No.	Description	Quantity
180E01-----2	Sartolab® RF 50, 0.22 µm, PES	24 units
180F01-----2	Sartolab® RF 50, 0.45 µm, PES	24 units
180E01-----E8	Sartolab® RF 50, 0.22 µm, PES, non-sterile	96 units
180F01-----E8	Sartolab® RF 50, 0.45 µm, PES, non-sterile	96 units

**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 information, visit  
[www.sartorius.com](http://www.sartorius.com)