



Sartolab[®] Multistation

Hands-Free Filtration of Multiple Samples



Watch video
Sartolab[®] Multistation



Watch animation
Sartolab[®] Multistation

Simplifying Progress

SARTORIUS

Sartolab® Multistation

Hands-Free Filtration of Multiple Samples

- Hands-free filtration
- One single vacuum source enables simultaneous filtration of up to 6 samples
- Time-saving (no installation time for each filter unit before use)



E-Shop
Sartolab® Multistation



See attached file
Datasheet Sartolab® Multistation



Watch animation
Sartolab® Multistation

Sartolab® Multistation

Hands-Free Filtration of Multiple Samples

- Hands-free filtration
- One single vacuum source enables simultaneous filtration of up to 6 samples
- Time-saving (no installation time for each filter unit before use)

One Vacuum Source Connection

- The Multistation is permanently connected to your vacuum source.



E-Shop
Sartolab® Multistation



See attached file
Datasheet Sartolab® Multistation



Watch animation
Sartolab® Multistation

Sartolab® Multistation

Hands-Free Filtration of Multiple Samples

- Hands-free filtration
- One single vacuum source enables simultaneous filtration of up to 6 samples
- Time-saving (no installation time for each filter unit before use)

Bracket for Hands-Free Filtration

- With one click, connect the filtration unit to the device, assuring perfect filter stability.
- With a second click, engage the vacuum automatically and begin filtering.



E-Shop
Sartolab® Multistation



See attached file
Datasheet Sartolab® Multistation



Watch animation
Sartolab® Multistation

Sartolab® Multistation

Hands-Free Filtration of Multiple Samples

- Hands-free filtration
- One single vacuum source enables simultaneous filtration of up to 6 samples
- Time-saving (no installation time for each filter unit before use)

360° Rotating Head

- Easily manipulate your samples with the rotating, multi-directional head.



E-Shop
Sartolab® Multistation



See attached file
Datasheet Sartolab® Multistation



Watch animation
Sartolab® Multistation

Sartolab® Multistation

Hands-Free Filtration of Multiple Samples

- Hands-free filtration
- One single vacuum source enables simultaneous filtration of up to 6 samples
- Time-saving (no installation time for each filter unit before use)

Numbered Brackets

- Easily keep track of your samples during filtration with the numbered brackets.



E-Shop
Sartolab® Multistation



See attached file
Datasheet Sartolab® Multistation



Watch animation
Sartolab® Multistation

Sartolab® Multistation

Hands-Free Filtration of Multiple Samples

- Hands-free filtration
- One single vacuum source enables simultaneous filtration of up to 6 samples
- Time-saving (no installation time for each filter unit before use)

Sartolab® Multistation

- **Order Number:** SDLC01



E-Shop
Sartolab® Multistation



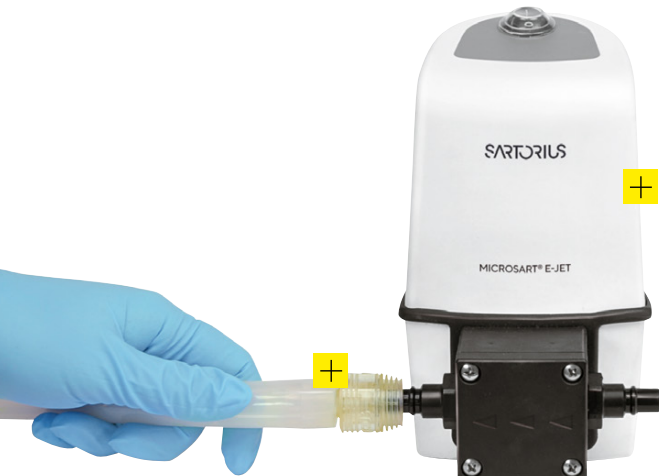
See attached file
Datasheet Sartolab® Multistation



Watch animation
Sartolab® Multistation

Microsart® E-jet Pump

- Small working surface due to compact design
- Easy-to-use
- Maintenance-free



E-Shop
Microsart® E-jet Pump



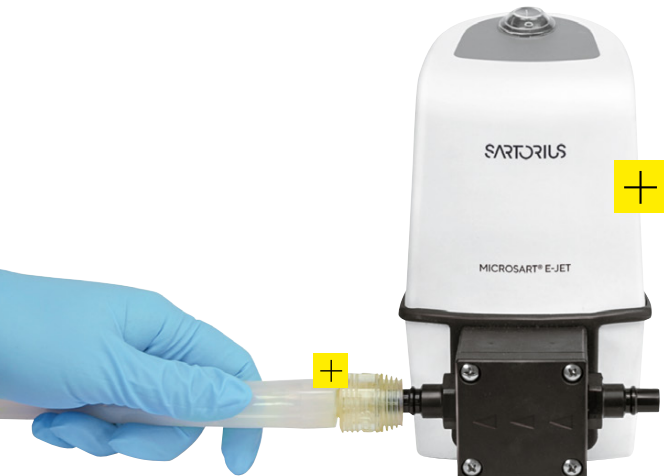
E-Shop
Vacuum Tube



See attached file
Microsart® E-jet Pump Datasheet

Microsart® E-jet Pump

- Small working surface due to compact design
- Easy-to-use
- Maintenance-free



Microsart® E-jet Pump

- Dimensions (W × L × H): 12 × 17 × 19 cm
- **Order Number:** 166MP-4



E-Shop
Microsart® E-jet Pump



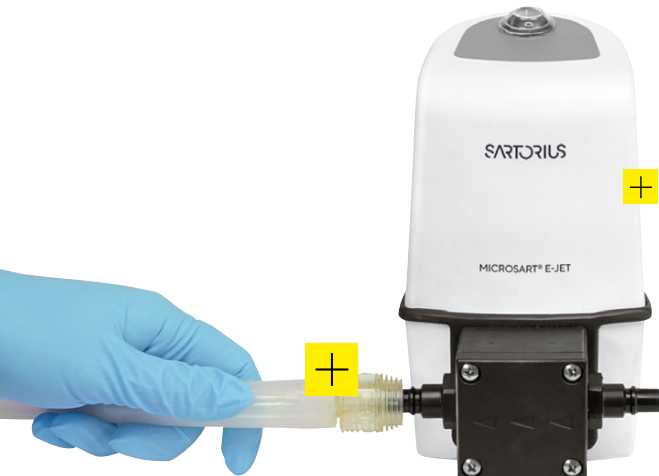
E-Shop
Vacuum Tube



See attached file
Microsart® E-jet Pump Datasheet

Microsart® E-jet Pump

- Small working surface due to compact design
- Easy-to-use
- Maintenance-free



Vacuum Tube

- Material: silicone
- Outer diameter: 15 mm
- Inner diameter: 7 mm
- Length: 200 mm
- Order Number: 1ZA---0007



E-Shop
Microsart® E-jet Pump



E-Shop
Vacuum Tube



See attached file
Microsart® E-jet Pump Datasheet

Consumables

Description

Sartolab® RF 50

Sartolab® RF 150-1000

Sartolab® BT 150-1000

Sartolab® Bottles

Sartolab® RF 50

Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF 50 Datasheet

Sartolab® RF 50

Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Yellow Screw Cap

- For storage of the filtrate
- Material: Polyethylene
- Packaging: Single-packaged, sterile



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF 50 Datasheet

Sartolab[®] RF 50

Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Sartolab[®] RF 50

Vacuum Filtration Units



Order Number	Description	Qty. per box
180E01-----2	Sartolab [®] RF 50, 0.22 µm, PES	24
180F01-----2	Sartolab [®] RF 50, 0.45 µm, PES	24
180E01-----E8	Sartolab [®] RF 50, 0.22 µm, PES, non-sterile	96
180F01-----E8	Sartolab [®] RF 50, 0.45 µm, PES, non-sterile	96



Sartolab[®] RF 50

Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Sartoclear Dynamics[®] Lab V50 Kits

Order Number	Description	Qty. per box
Sartolab[®] RF 50, 0.22 µm, PES		
SDLV-0050-01E0-2	Includes 1 pouch of 1 g DE filter aid per filtration unit (24 pouches)	24
SDLV-0050-02E0-2	Includes 2 pouches of 1 g DE filter aid per filtration unit (48 pouches)	24
Sartolab[®] RF 50, 0.45 µm, PES		
SDLV-0050-01F0-2	Includes 1 pouch of 1 g DE filter aid per filtration unit (24 pouches)	24
SDLV-0050-02F0-2	Includes 2 pouches of 1 g DE filter aid per filtration unit (48 pouches)	24



Sartolab® RF 50

Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Funnel With Dust Cover

- Material: Polystyrene (PS)
- Membrane filter:
 - 0.22 µm polyethersulfone (order no. 180E01)
 - 0.45 µm polyethersulfone (order no. 180F01)
- Membrane diameter: 58.5 mm
- Effective filtration area: 21 cm²
- Dead volume: 1.2 mL



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF 50 Datasheet

Sartolab® RF 50

Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Funnel Adapter

- 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 without an additional vacuum connector.
- Material: Polyethylene



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF 50 Datasheet

Sartolab® RF 50

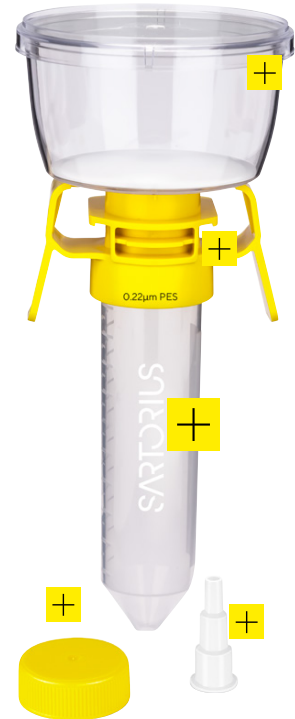
Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

50 mL Conical Tube

- Material: Polypropylene
- Volume: 50 mL
- Height: 115.5 mm
- External | internal diameter: 29.5 mm | 27.48 mm



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF 50 Datasheet

Sartolab® RF 50

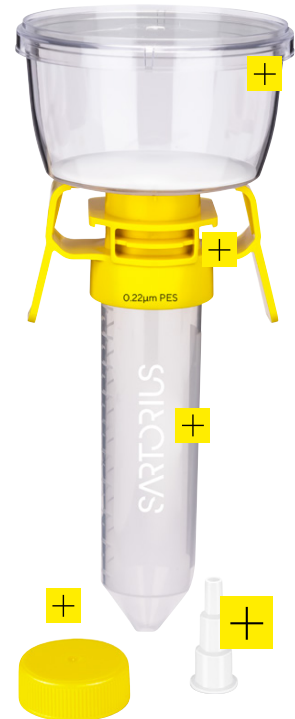
Vacuum Filtration Unit for Volumes up to 50 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Vacuum Connector ✕

- The vacuum connector is provided with each filtration unit for a direct connection to the vacuum tube (without the Multistation).
- Packaging: Single-packaged, sterile
- Not delivered with non-sterile versions of Sartolab® RF 50



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF 50 Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit



Sartolab® RF 150 - 1000 Vacuum Filtration Units



Order Number	Description	Membrane Type	Pore Size	Funnel Volume (mL)	Bottle Volume (mL)	Qty. per Box
180E02-----E	Sartolab® RF 150	Asymmetric PES	0.22 µm	150	150	12
180F02-----E	Sartolab® RF 150	PES	0.45 µm	150	150	12
180D03-----E	Sartolab® RF 250	PES	0.1 µm	250	250	12
180E03-----E	Sartolab® RF 250	Asymmetric PES	0.22 µm	250	250	12
180F03-----E	Sartolab® RF 250	PES	0.45 µm	250	250	12
180E04-----E	Sartolab® RF 500	Asymmetric PES	0.22 µm	500	500	12
180F04-----E	Sartolab® RF 500	PES	0.45 µm	500	500	12
180D05-----E	Sartolab® RF 1,000	PES	0.1 µm	1,000	1,000	12
180E05-----E	Sartolab® RF 1,000	Asymmetric PES	0.22 µm	1,000	1,000	12
180F05-----E	Sartolab® RF 1,000	PES	0.45 µm	1,000	1,000	12

Sartolab® RF 150-1000

Sartoclear Dynamics® Lab V Kits



Order Number	Description	Kit Contents	Qty. per Box
150 mL			
SDLV-0150-02E0-E	Sartoclear Dynamics® Lab V150, RF, 0.22 µm, 2 g	2 pouches of 1 g DE filter aid per filtration unit	12
SDLV-0150-05E0-2	Sartoclear Dynamics® Lab V150, RF, 0.22 µm, 5 g	1 pouch of 5 g DE filter aid per filtration unit	24
250 mL			
SDLV-0250-05E0-2	Sartoclear Dynamics® Lab V250, RF, 0.22 µm, 5 g	1 pouch of 5 g DE filter aid per filtration unit	24
SDLV-0250-10E0-2	Sartoclear Dynamics® Lab V250, RF, 0.22 µm, 10 g	1 pouch of 10 g DE filter aid per filtration unit	24
500 mL			
SDLV-0500-05E0-2	Sartoclear Dynamics® Lab V500, RF, 0.22 µm, 5 g	1 pouch of 5 g DE filter aid per filtration unit	24
SDLV-0500-10E0-2	Sartoclear Dynamics® Lab V500, RF, 0.22 µm, 10 g	1 pouch of 10 g DE filter aid per filtration unit	24
SDLV-0500-20E0-E	Sartoclear Dynamics® Lab V500, RF, 0.22 µm, 20 g	1 pouch of 20 g DE filter aid per filtration unit	12
1000 mL			
SDLV-1000-10E0-2	Sartoclear Dynamics® Lab V1000, RF, 0.22 µm, 10 g	1 pouch of 10 g DE filter aid per filtration unit	24
SDLV-1000-20E0-E	Sartoclear Dynamics® Lab V1000, RF, 0.22 µm, 20 g	1 pouch of 20 g DE filter aid per filtration unit	12
SDLV-1000-40E0-E	Sartoclear Dynamics® Lab V1000, RF, 0.22 µm, 40 g	2 pouches of 20 g DE filter aid per filtration unit	12

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Graduated Funnel with Dust Cover [×](#)

- Material: Polystyrene (PS)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Specifications

Membrane Filter:

0.1 µm, 0.22 µm, 0.45 µm PES

Membrane Diameter:

80 mm for 150 mL and 250 mL volumes
100 mm for 500 mL and 1,000 mL volumes

Effective Filtration Area:

43 cm² for 150 mL and 250 mL volumes
69 cm² for 500 mL and 1,000 mL volumes

Dead Volume:

2.7 mL for 150 mL and 250 mL versions
4.1 mL for 500 mL and 1,000 mL versions



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Funnel Adapter

The design of the yellow adapter connecting the funnel to the flask ensures a vacuum-tight seal and enables the filtration unit to be used on the Sartolab® Multistation without an additional vacuum connector.

Material: High Density Polyethylene (HDPE)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Vacuum Connector ✕

The vacuum connector is provided with each filtration unit for a direct connection to the vacuum tube (without the Multistation).

Material: High Density Polyethylene (HDPE)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Yellow Screw Cap ✕

- For storage of the filtrate
- Material: Polyethylene
- Individually wrapped to maintain sterility



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® RF 150-1000

Vacuum Filtration Unit for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Minimized risk of contamination with a complete ready-to-use unit

Ordering Information (click to expand): [+](#)

Receiver Flask

- Graduation and writing field
- Bottle neck size: 45 mm
- One-hand handling
- Indented bottom for stackability
- Material: Polystyrene (PS)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® BT 150-1000

Bottle Top Filters for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Flexibility to use any bottles suitable for vacuum with a neck size of 45 mm

Ordering Information (click to expand): [+](#)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® BT 150-1000

Bottle Top Filters for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Flexibility to use any bottles suitable for vacuum with a neck size of 45 mm

Sartolab® BT 150 - 1000 Bottle Top Filtration Units

Order Number	Description	Membrane Type	Pore Size	Funnel Volume (mL)	Qty. per Box
180E12-----E	Sartolab® BT 150	Asymmetric PES	0.22 µm	150	12
180E13-----E	Sartolab® BT 250	Asymmetric PES	0.22 µm	250	12
180E14-----E	Sartolab® BT 500	Asymmetric PES	0.22 µm	500	12
180E15-----E	Sartolab® BT 1,000	Asymmetric PES	0.22 µm	1,000	12
180F15-----E	Sartolab® BT 1,000	PES	0.45 µm	1,000	12



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet



Sartolab® BT 150-1000

Bottle Top Filters for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Flexibility to use any bottles suitable for vacuum with a neck size of 45 mm

Ordering Information (click to expand): [+](#)

Graduated Funnel with Dust Cover

- Material: Polystyrene (PS)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab[®] BT 150-1000

Bottle Top Filters for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Flexibility to use any bottles suitable for vacuum with a neck size of 45 mm

Specifications

Membrane Filter:

0.22µm, 0.45 µm PES

Membrane Diameter:

80 mm for 150 mL and 250 mL volumes
100 mm for 500 mL and 1,000 mL volumes

Effective Filtration Area:

43 cm² for 150 mL and 250 mL volumes
69 cm² for 500 mL and 1,000 mL volumes

Dead Volume:

2.7 mL for 150 mL and 250 mL versions
4.1 mL for 500 mL and 1,000 mL versions



Sartolab® BT 150-1000

Bottle Top Filters for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Flexibility to use any bottles suitable for vacuum with a neck size of 45 mm

Ordering Information (click to expand): [+](#)

Funnel Adapter

The design of the yellow adapter connecting the funnel to the flask ensures a vacuum-tight seal and enables the filtration unit to be used on the Sartolab® Multistation without an additional vacuum connector.

Material: High Density Polyethylene (HDPE)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® BT 150-1000

Bottle Top Filters for Volumes from > 50 mL up to 1000 mL

- Highest flow rates with an asymmetric membrane and a large surface of filtration
- No loss of proteins with a very low protein binding membrane
- Low dead volume thanks to an optimized membrane support
- Flexibility to use any bottles suitable for vacuum with a neck size of 45 mm

Ordering Information (click to expand): [+](#)

Vacuum Connector ✕

The vacuum connector is provided with each filtration unit for a direct connection to the vacuum tube (without the Multistation)

Material: High Density Polyethylene (HDPE)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® RF | BT Datasheet

Sartolab® Bottles

Single-Use Polystyrene Storage Bottles

- Ergonomically designed
- Available in 4 sizes (150 mL, 250 mL, 500 mL and 1,000 mL)
- Compatible with all Sartolab® RF | BT150-1000 vacuum filtration units

Ordering Information (click to expand): [+](#)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® Bottles Datasheet

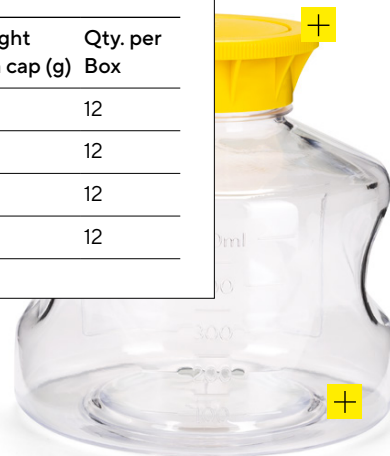
Sartolab® Bottles

Single-Use Polystyrene Storage Bottles

- Ergonomically designed
- Available in 4 sizes (150 mL, 250 mL, 500 mL and 1,000 mL)
- Compatible with all Sartolab® RF | BT150-1000 vacuum filtration units

Sartolab® Bottles

Order Number	Description	Nominal Capacity (mL)	Diameter (mm)	Height (mm)	Weight with cap (g)	Qty. per Box
180-22-----E	Sartolab® Bottle, 150 mL	150	87	77	49	12
180-23-----E	Sartolab® Bottle, 250 mL	250	87	99	57	12
180-24-----E	Sartolab® Bottle, 500 mL	500	112	114	87	12
180-25-----E	Sartolab® Bottle, 1,000 mL	1000	116	169	133	12



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® Bottles Datasheet

Sartolab® Bottles

Single-Use Polystyrene Storage Bottles

- Ergonomically designed
- Available in 4 sizes (150 mL, 250 mL, 500 mL and 1,000 mL)
- Compatible with all Sartolab® RF | BT150-1000 vacuum filtration units

Receiver Flask

- Graduation and writing field
- Bottle neck size: 45 mm
- One-hand handling
- Material: Polystyrene (PS)



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® Bottles Datasheet

Sartolab® Bottles

Single-Use Polystyrene Storage Bottles

- Ergonomically designed
- Available in 4 sizes (150 mL, 250 mL, 500 mL and 1,000 mL)
- Compatible with all Sartolab® RF | BT150-1000 vacuum filtration units

Yellow Screw Cap

- For storage of the filtrate
- Material: Polyethylene



E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® Bottles Datasheet

Sartolab® Bottles

Single-Use Polystyrene Storage Bottles

- Ergonomically designed
- Available in 4 sizes (150 mL, 250 mL, 500 mL and 1,000 mL)
- Compatible with all Sartolab® RF | BT150-1000 vacuum filtration units

Bottom

Intended for stackability




E-Shop
Sartolab® Vacuum Filtration Units



See attached file
Sartolab® Bottles Datasheet

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

 For further contacts, visit
www.sartorius.com

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906