

Sartolab® MultiStation

Hands-Free Parallel Filtration
of up to Six Samples



Benefits

- One single vacuum source to enable the filtration of up to 6 samples at the same time
- Hands-free filtration
- Time saving (no extra installation time to vacuum for each single filter unit)

Product Information

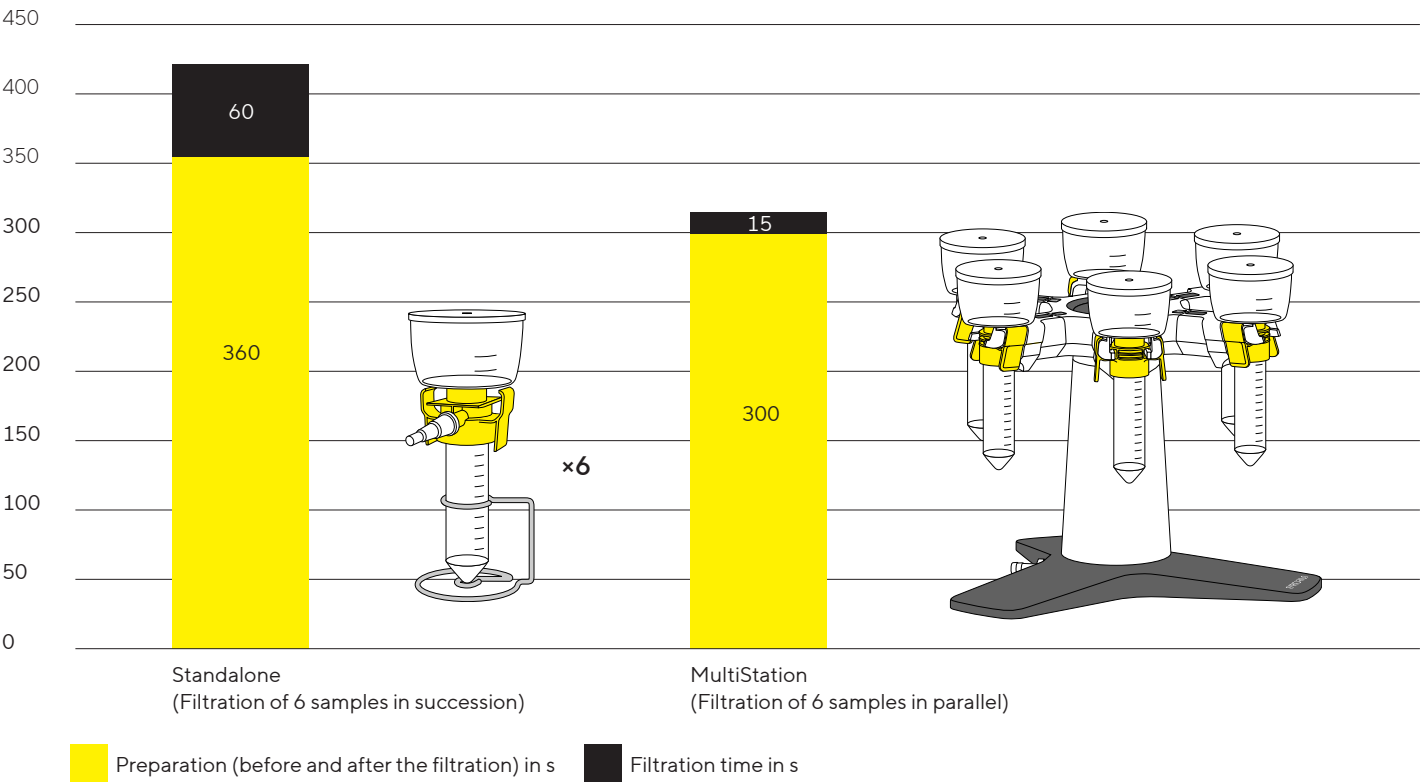
Sartolab® MultiStation is a stand especially designed to hold 1 to 6 vacuum filtration units. It allows a filtration of all samples at the same time. The MultiStation is permanently connected to a vacuum source; the vacuum filtration units can be easily installed in the MultiStation for quick and easy filtration of samples without extra time consuming connection.

Sartolab® MultiStation functions with the Sartolab® RF| BT vacuum filtration units. The adapter connecting the funnel to the sample tube or receiver flask perfectly fits into the bracket of the MultiStation.

One click and the filtration unit hangs in the device. That assures a perfect stability of the filter; with a second click, the vacuum is automatically applied and the filtration begins.

The rotating head is multi-directional and enables easy handling of your samples in front of you. Furthermore, all 6 brackets are numbered to avoid any confusion during the filtration of multiple samples.

Example of Filtration of 6 × 50 mL Samples with and without the Sartolab® MultiStation



Specifications

Material (visible parts)	ABS Aluminum Stainless steel
Dimensions (Length × Width × Height)	307 × 348 × 281 mm
Weight	4.6 kg
Tubing adaptor	Designed for tubing with an inner diameter between 4 and 10 mm and with a wall thickness of minimum 3 mm.

Ordering Information

Order Number	Description	Quantity
SDLC01	Sartolab® MultiStation for the filtration of up to 6 samples in parallel	1 unit

Consumables

Sartolab® RF (Vacuum Filtration Units with Receiver Flasks)

Order Number	Description	Membrane	Pore Size (µm)	Volume of the Funnel (mL)	Volume of the Bottle (mL)	Sterilization	Packing Unit
180E01-----E8	Sartolab® RF 50	PES	0.22	50	50	No	96
180F01-----E8	Sartolab® RF 50	PES	0.45	50	50	No	96
180E01-----2	Sartolab® RF 50	PES	0.22	50	50	Yes	24
180F01-----2	Sartolab® RF 50	PES	0.45	50	50	Yes	24
180E02-----E	Sartolab® RF 150	PES	0.22	150	150	Yes	12
180F02-----E	Sartolab® RF 150	PES	0.45	150	150	Yes	12
180D03-----E	Sartolab® RF 250	PES	0.1	250	250	Yes	12
180E03-----E	Sartolab® RF 250	PES	0.22	250	250	Yes	12
180F03-----E	Sartolab® RF 250	PES	0.45	250	250	Yes	12
180E04-----E	Sartolab® RF 500	PES	0.22	500	500	Yes	12
180F04-----E	Sartolab® RF 500	PES	0.45	500	500	Yes	12
180D05-----E	Sartolab® RF 1000	PES	0.1	1000	1000	Yes	12
180E05-----E	Sartolab® RF 1000	PES	0.22	1000	1000	Yes	12
180F05-----E	Sartolab® RF 1000	PES	0.45	1000	1000	Yes	12

Sartolab® BT (Bottle Top Filters without Receiver Flasks)

Order Number	Description	Membrane	Pore Size (µm)	Volume of the Funnel (mL)	Sterilization	Packing Unit
180E12-----2	Sartolab® BT 150	PES	0.22	150	Yes	24
180E13-----2	Sartolab® BT 250	PES	0.22	250	Yes	24
180E14-----2	Sartolab® BT 500	PES	0.22	500	Yes	24
180E15-----2	Sartolab® BT 1000	PES	0.22	1000	Yes	24
180F15-----2	Sartolab® BT 1000	PES	0.45	1000	Yes	24

Sartolab® Bottles

Order Number	Description	Volume (mL)	Sterilization	Packing Unit
180-22-----2	Sartolab® bottle	150	Yes	24
180-23-----2	Sartolab® bottle	250	Yes	24
180-24-----2	Sartolab® bottle	500	Yes	24
180-25-----2	Sartolab® bottle	1000	Yes	24

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
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