

# Sartolab® P20 Pressure Filter Units

Compact Design for the  
Filtration of Large Volumes



## Benefits

- Highest flow rates with a large surface of filtration (20 cm<sup>2</sup>)
- No loss of protein with a low binding membrane
- Low dead volume due to an optimized membrane support
- Versions available with a prefilter for high particle load solutions

## Product Information

Sartolab® P20 devices are ready-to-use pressure filter units for general laboratory use like the clarification and sterile filtration of media and aqueous solutions in batches from 100 mL to 10 L. Sartolab® P20 Plus with an incorporated prefilter is recommended for difficult-to-filter solutions like, for example, media that contains serum. The Sartolab® P20 Prefilter is designed to clarify solutions with high particle loads.

### Membrane of Choice

Polyethersulfone (PES) is the membrane of choice for the Sartolab® P20 pressure filter units, as it combines very low protein binding properties with the highest flow rates. The Sartolab® P20 pressure filter unit is available either with 0.2 µm or 0.45 µm PES membranes, with or without a prefilter made of high purity quartz microfibers. An additional version containing glass microfiber only is also available for clarification purposes.

### Compact Design

Sartolab® P20 pressure filter units have been designed to filter batches from 100 mL to 10 L, either using a syringe or in-line with a peristaltic pump, or a pressure vessel. Sartolab® P20 pressure filter units are available in different configurations, with or without PTFE automatic venting, with or without a filling bell (including cover) on the outlet and with a combination of different inlet and outlet connectors to meet the needs of most applications.

## Technical Specifications

Different Filter Materials	0.2 µm polyethersulfone 0.45 µm polyethersulfone High purity binder-free quartz microfibers High purity binder-free glass microfibers
Housing Material	Transparent polycarbonate
Filter Housing Diameter	61 mm
Filtration Area	20 cm <sup>2</sup>
Holdup Volume	Sartolab® P20: 1 mL Sartolab® P20 Plus: 1.2 mL Sartolab® P20 Prefilter: 1 mL
Filtration Range	Sartolab® P20: 100 mL to 5 L Sartolab® P20 Plus: 100 mL to 10 L Sartolab® P20 Prefilter: 100 mL to 10 L
Recommended Max. Inlet Pressure	Sartolab® P20 and Sartolab® P20 Plus: 4 bar Sartolab® P20 Prefilter: 0.8 bar
pH Range	1–10
Housing Burst Pressure	≥ 5 bar
Autoclavable	121°C
Packaging	Sterile versions: single-packaged in a tyvek pouch Non sterile versions: bulk-packaged

# Ordering Information

## Sartolab® P20

Order Number	Filter Material	Inlet	Outlet	PTFE Venting	Filling Bell	Sterilization (EO)	Qty./ Pack
18075-----D	0.2 µm PES	Female Luer-Lock	Hose barb	no	no	yes	10
18075-----UPN	0.2 µm PES	Female Luer-Lock	Hose barb	no	no	no	100
18089-----D	0.2 µm PES	Hose barb	Hose barb	yes	yes	yes	10
18090-----D	0.2 µm PES	Female Luer-Lock	Male Luer-Lock	no	no	yes	10

## Sartolab® P20 Plus

Order Number	Filter Material	Inlet	Outlet	PTFE Venting	Filling Bell	Sterilization (EO)	Qty./ Pack
18068-----D	Quartz microfibers & 0.2 µm PES	Female Luer-Lock	Hose barb	yes	yes	yes	10
18076-----N	Quartz microfibers & 0.45 µm PES	Hose barb	Hose barb	no	no	no	100
18091-----D	Quartz microfibers & 0.2 µm PES	Hose barb	Hose barb	yes	yes	yes	10
18092-----D	Quartz microfibers & 0.2 µm PES	Female Luer-Lock	Male Luer-Lock	no	no	yes	10

## Sartolab® P20 Prefilter


Order Number	Filter Material	Inlet	Outlet	PTFE Venting	Filling Bell	Sterilization (EO)	Qty./ Pack
18072-----D	Glass microfibers	Female Luer-Lock	Hose barb	no	no	no	10

## 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](http://www.sartorius.com)