



Sartolab® P20 Pressure Filter Units

Compact Design for the Filtration of Large Volumes

Simplifying Progress

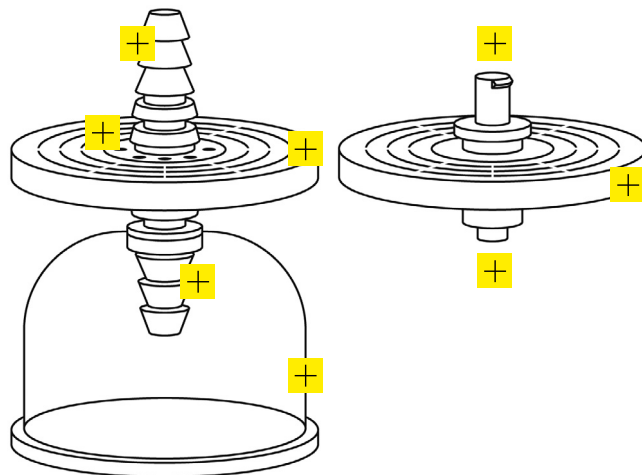
SARTORIUS

Sartolab® P20

Ready-to-use pressure filter units for the clarification and sterile filtration of media and aqueous solutions in batches from 100 mL to 5 L.

- Highest flow rates with a large surface of filtration (20 cm²)
- No loss of protein with a low binding 0.2 µm polyethersulfone (PES) membrane
- Low hold-up volume due to an optimized membrane support
- Several inlet & outlet connectors combinations (hose barb or Luer-Lock)
- With or without automatic PTFE venting holes
- With or without filling bell

Types available (click to expand): 



E-shop
Sartolab® P20



Datasheet
Sartolab® P20



User Manual
Sartolab® P20

Sartolab® P20

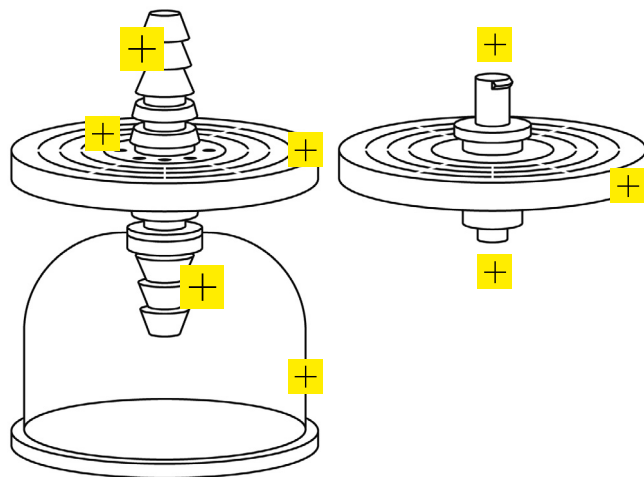
Types Available

Item 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

Hose Barb

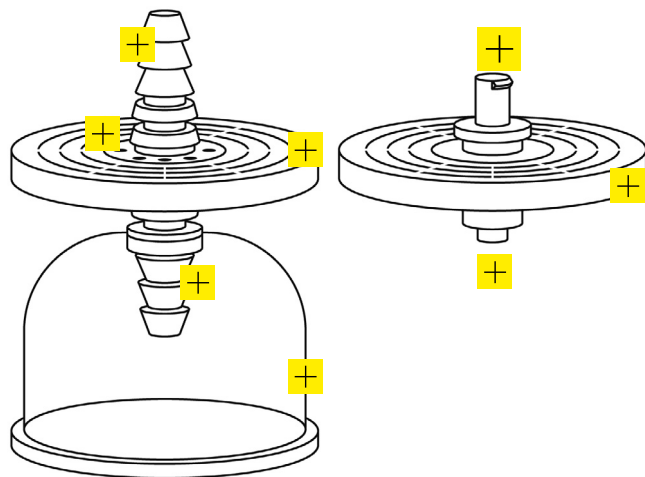
Compatible with tubing
with an internal diameter
of 6–12 mm



Sartolab® P20

Inlet: Female Luer-Lock

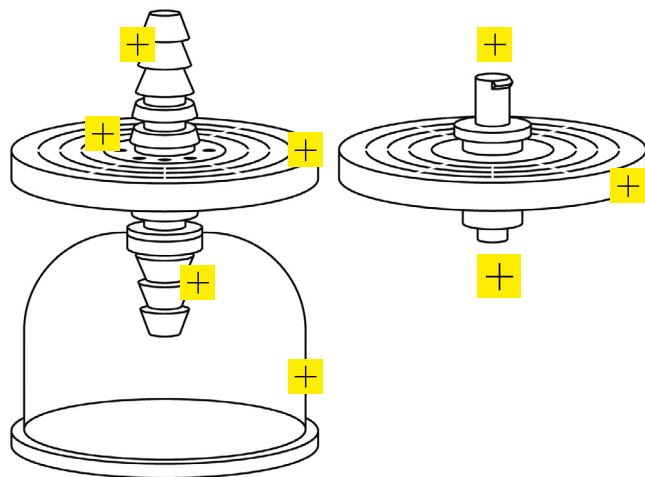
Compatible with Luer-Lock syringes and tubing with a male Luer-Lock connector



Sartolab® P20

Outlet: Male Luer-Lock

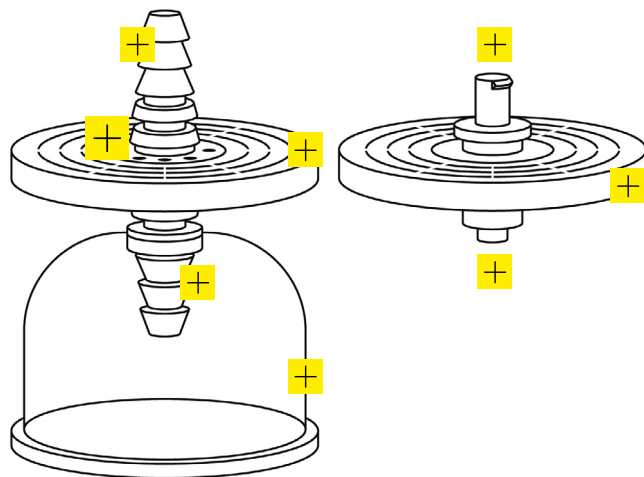
Compatible with tubing
with a female Luer-Lock
connector



Sartolab[®] P20

Automatic Venting Holes

with a PTFE membrane



Sartolab® P20

Membrane & Housing

Membrane

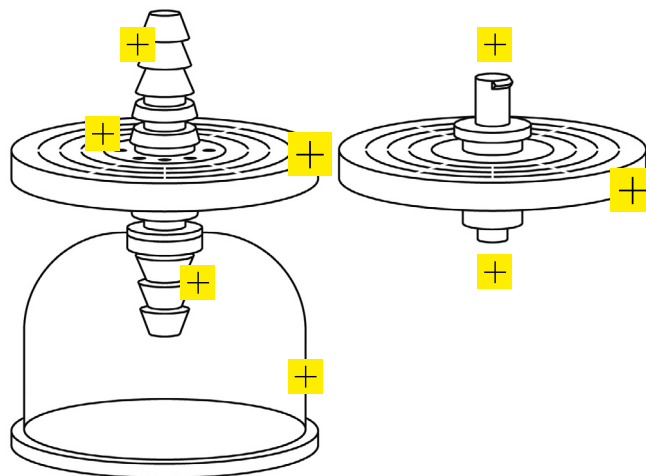
- Material: Polyethersulfone (PES)
- Pore size: 0.2 µm
- Effective filtration surface: 20 cm²
- Filter diameter: 56 mm

Housing

- Material: Polycarbonate

Dimensions

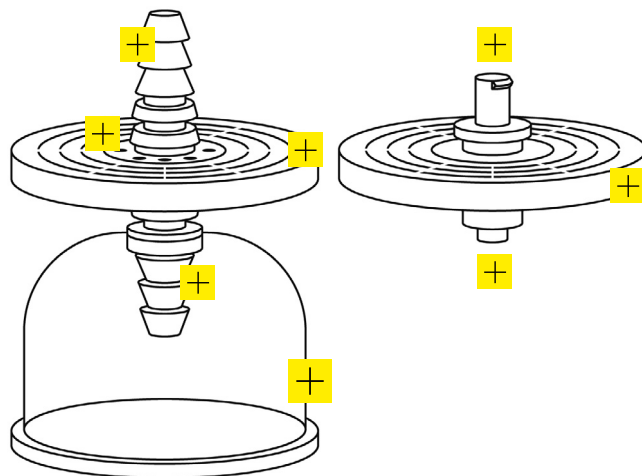
- Hose barb inlet and outlet: 63 x 61 mm
- Female Luer-Lock inlet and hose barb outlet: 50 x 61 mm
- Female Luer-Lock inlet and male Luer-Lock outlet: 31 x 61 mm



Sartolab® P20

Filling Bell with Lid

Used to protect the filtrate against external contamination

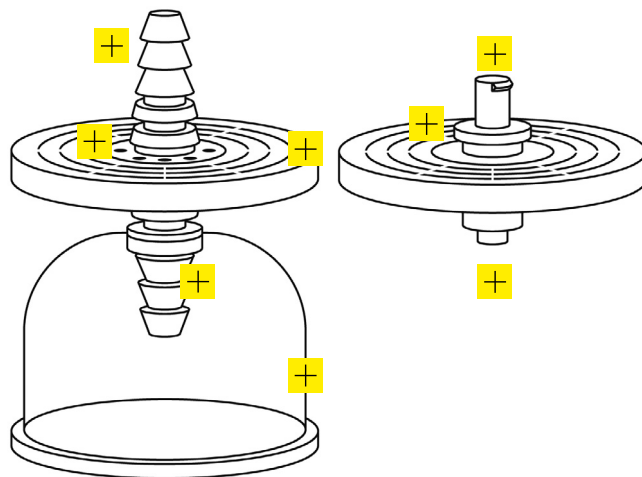


Sartolab® P20 Plus

Ready-to-use pressure filter units for the clarification and sterile filtration of media and aqueous solutions in batches from 100 mL to 10 L.

- Highest flow rates with a large surface of filtration (20 cm²)
- No loss of protein with a low binding 0.2 µm or 0.45 µm polyethersulfone (PES) membrane
- Depth filter as prefilter for highest throughput
- Low hold-up volume due to an optimized membrane support
- Several inlet & outlet connectors combinations (hose barb or Luer-Lock)
- With or without automatic PTFE venting holes
- With or without filling bell

Types available (click to expand): [+](#)



E-shop
Sartolab® P20



Datasheet
Sartolab® P20



User Manual
Sartolab® P20

Sartolab® P20 Plus

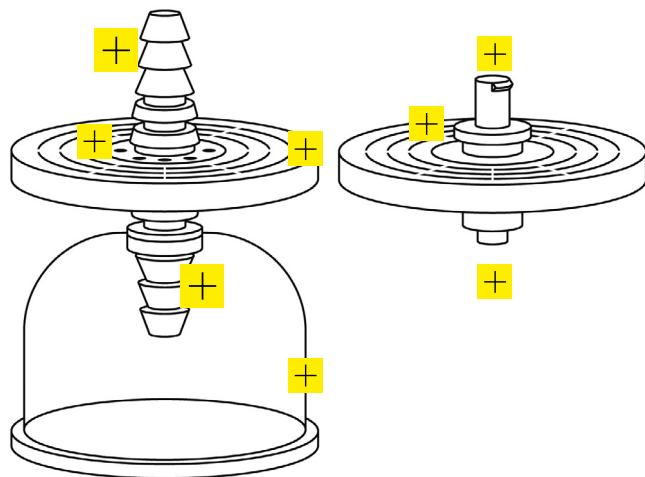
Types Available

Item 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 Plus

Hose Barb

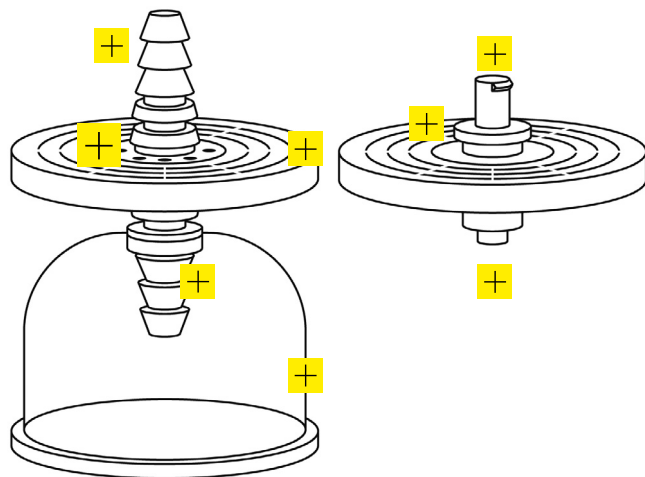
Compatible with tubing
with an internal diameter
of 6-12 mm



Sartolab® P20 Plus

Automatic Venting Holes

with a PTFE membrane

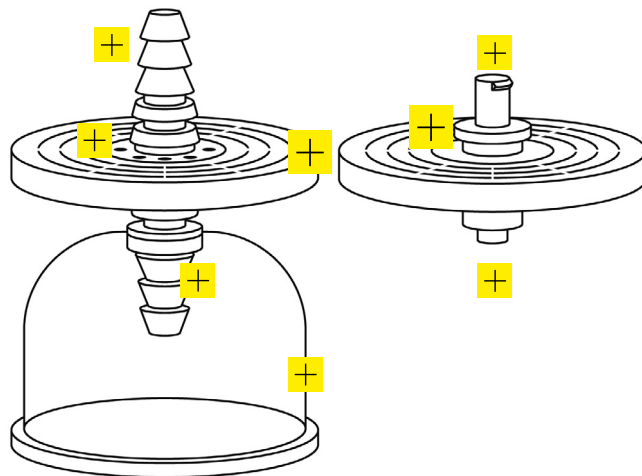


Sartolab® P20 Plus

- Prefilter made of 100 % binder-free high-purity quartz microfibers in combination with either a 0.2 µm PES membrane for sterile filtration or 0.45 µm PES membrane for clarification
- Effective filtration surface: 20 cm²
- Filter diameter: 56 mm
- Housing diameter: 61 mm
- Housing material: Polycarbonate

Housing Dimensions

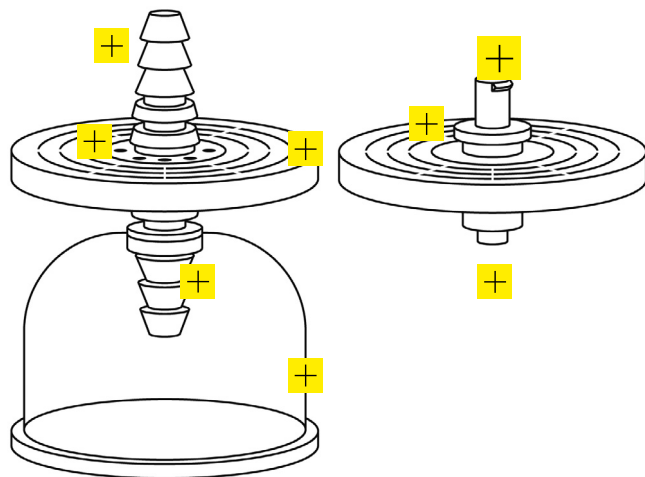
- Hose barb inlet and outlet: 63 x 61 mm
- Female Luer-Lock inlet and hose barb outlet: 50 x 61 mm
- Female Luer-Lock inlet and male Luer-Lock outlet: 31 x 61 mm



Sartolab® P20 Plus

Inlet: Female Luer-Lock

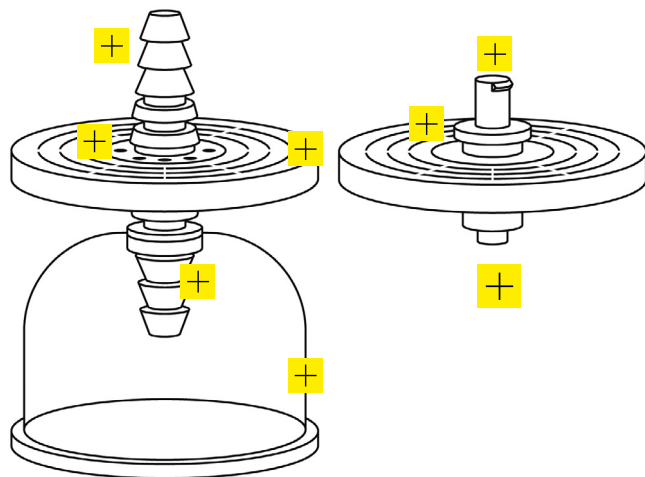
Compatible with Luer-Lock syringes and tubing with a male Luer-Lock connector



Sartolab® P20 Plus

Outlet: Male Luer-Lock

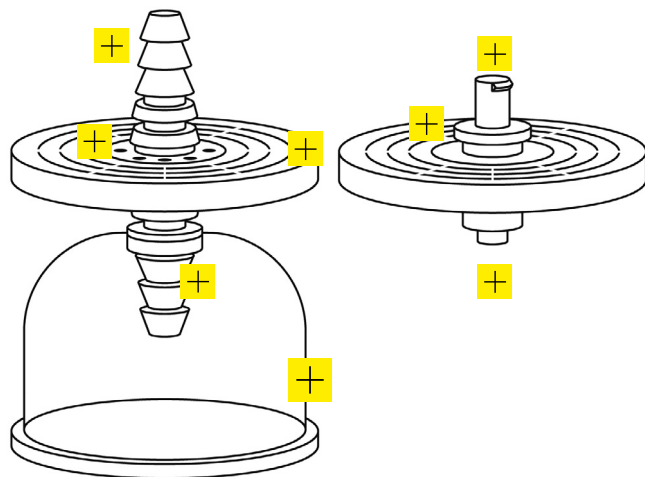
Compatible with tubing
with a female Luer-Lock
connector



Sartolab® P20 Plus

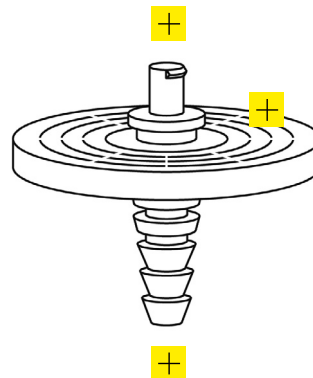
Filling Bell with Lid

Used to protect the filtrate against external contamination



Sartolab® P20 Prefilter

- Prefilter made of 100 % binder-free high-purity glass microfibers
 - Highest flow rates with a large surface of filtration (20 cm²)
 - Low hold-up volume due to an optimized membrane support
 - 10 pieces per pack, non-sterile
- **Item Number: 18072-----D**



E-shop
Sartolab® P20



Datasheet
Sartolab® P20

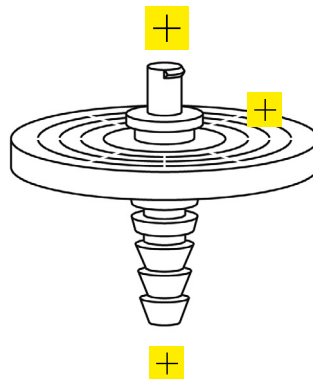


User Manual
Sartolab® P20

Sartolab® P20 Prefilter

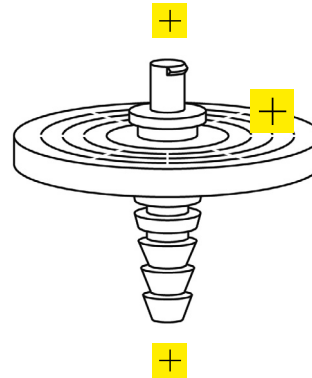
Inlet: Female Luer-Lock

Compatible with Luer-Lock syringes and tubing with a male Luer-Lock connector



Sartolab® P20 Prefilter

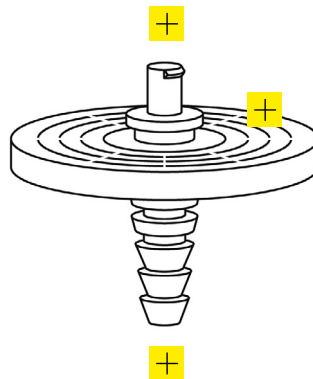
- Prefilter made of 100 % binder-free high-purity glass microfibers
- Effective filtration surface: 20 cm²
- Filter diameter: 56 mm
- Housing diameter: 61 mm
- Housing material: Polycarbonate
- Housing dimensions: 50 x 61 mm



Sartolab® P20 Prefilter

Outlet: Hose Barb

Compatible with tubing
with an internal diameter
of 6-12 mm



Accessories

Item Number	Description	
18099	Sartolab® Integrity Test Holder	
6988094	Tubing Set for Luer-Lock Connectors	
180MP-1	Diaphragm Pump	
180MP-1B	Pump Puls Footswitch	
180MP-1F	Pump Tripod Holder	
180MP-1G	Pump Mounting Plate	



E-shop
Sartolab® P20



Datasheet
Sartolab® P20

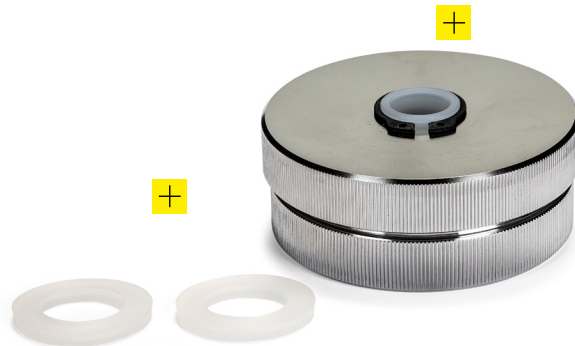


User Manual
Sartolab® P20

Sartolab® Integrity Test Holder

For integrity testing, a bubble point test can be performed on the Sartolab® P20 and Sartolab® P20 Plus after use

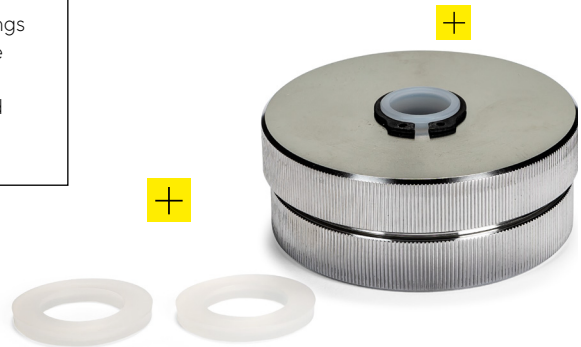
- If the filter has PTFE venting holes, use the integrity test holder to close the holes and enable a bubble point.
 - The holder is not necessary for Sartolab® P20 and Sartolab® P20 Plus without automatic venting.
- **Item Number: 18099**



Sartolab® Integrity Test Holder

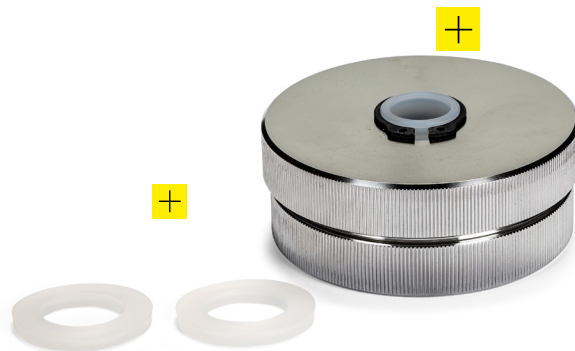
The silicone gasket seals the openings above the PTFE venting membrane

The integrity test holder is delivered with 2 extra silicone gaskets.



Sartolab® Integrity Test Holder

Material of the device: stainless steel



Tubing Set for Luer-Lock Connectors

Item number: 6988094

Set contains:

- Silicone tubing (Length: 2.5 m, internal diameter: 4 mm, thickness: 1.5 mm)
- Suction foot: a weight to keep the tubing into the liquid
- Male Luer-Lock connectors with a blue union nut (5 pieces in a reclosable bag)



Diaphragm Pump

Item number: 180MP-1

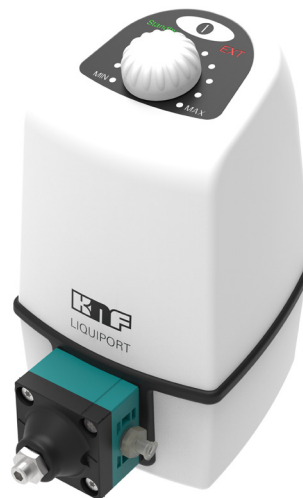
The Liquiport PML 16980 pump by KNF is optimized for use with Sartolab® P20 pressure filtration devices and features:

- An integrated pressure relief valve set at 3.5 bar for secure operation
- Dry running capability
- Integrated overheating protection
- Splashproof construction with an IP65 rating
- Compact design with a small footprint

Material

- Pump head: PP
- Valve plate: / seals FFKM
- Diaphragm: PTFE-coated
- Housing: PA, TPE, PC

Net weight: 1 kg



Datasheet
Diaphragm Pump & Accessories



User Manual
Sartolab® P20

Pump Puls Footswitch

Item number: 180MP-1B

The pulse footswitch for Liqueport PML 16980 provides some advantages such as:

- Hands-free operation: Enhances ergonomics
- Increased safety: Reduces the risk of contamination in sensitive environment
- Improved efficiency: Provides quick and easy start/stop control

Material

- Steel housing painted in black

Net weight: 0.22 g



Datasheet
Diaphragm Pump & Accessories



User Manual
Sartolab® P20

Pump Tripod Holder

Item number: 180MP-1F

The pulse footswitch for Liquiport PML 16980 provides some advantages such as:

- Space Efficiency: Elevates the pump off the work surface, freeing up valuable bench space
- Improved Safety: By securely holding the pump in place, a tripod holder minimizes the risk of accidental spills or damage, contributing to a safer working environment.

Material

- Stainless steel 1.4301

Net weight: 0.12 g



Datasheet
Diaphragm Pump & Accessories



User Manual
Sartolab® P20

Pump Mounting Plate

Item number: 180MP-1G

The mounting plate for Liquiport PML 16980 provides several advantages such as:

- Space Efficiency: Allows the pump to be mounted on vertical surfaces, freeing up valuable workspace.
- Improved Safety: By securely fixing the pump in place, a mounting plate reduces the risk of movement, contributing to a safer working environment.

Material

- Stainless steel 1.4301

Net weight: 0.2 g



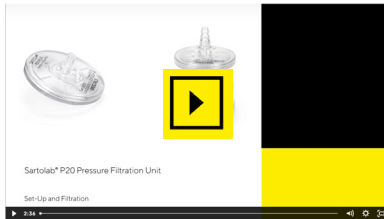
Datasheet
Diaphragm Pump & Accessories



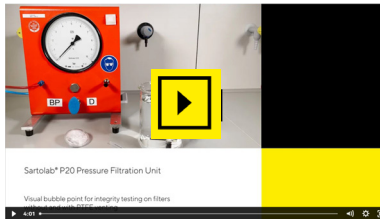
User Manual
Sartolab® P20



Videos



Setup and Filtration of Sartolab®
P20 Pressure Devices



How to Perform a Visual Bubble
Point with a Sartolab P20
Pressure Filtration Device



Cleaning & Maintenance of the
KNF Liquoport 100 Diaphragm
Pump

Germany

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

USA

Sartorius Corporation
3874 Research Park Dr.
Ann Arbor, MI 48108
Phone +1 734 769 1600

 For further contacts, visit
www.sartorius.com