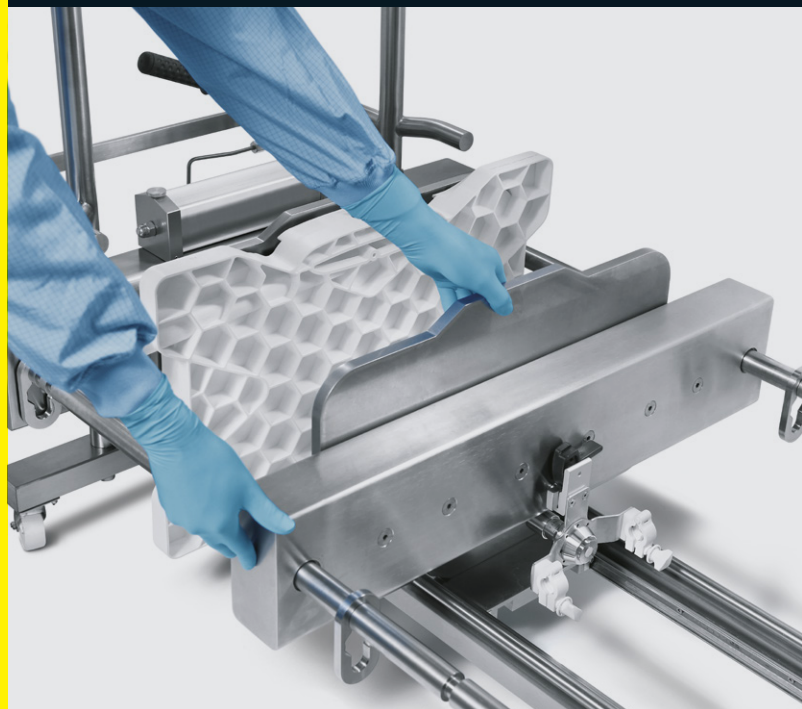


Filter Holders

For Sartoclear® Depth Filters
and Sartobind® Membrane
Adsorbers



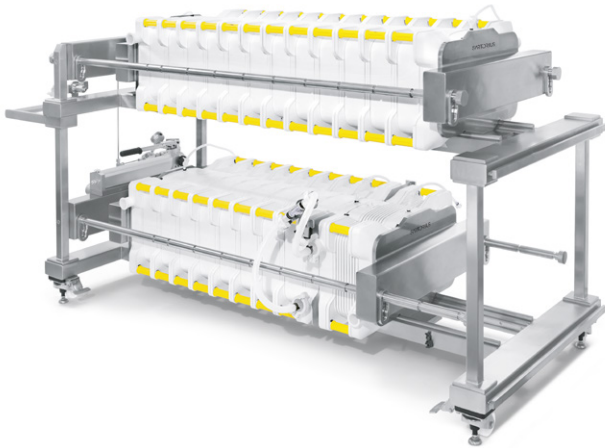
Product Information

Filter holders are hydraulic clamping systems especially designed for the use of modular Sartoclear® clarification filters and Sartobind® membrane chromatography cassettes. The holders are available in three standard formats for pilot-scale and commercial manufacturing.

One Platform for Two Technologies

Sartoclear® depth filters and Sartobind® cassettes are two complementing technologies using the same holder and manifold plates. Use Sartoclear® depth filters for applications with moderate solids content (see table for more application guidance).

If you need large scale membrane chromatography for virus | VLP large protein capture, use Sartobind® anion or cation exchanger cassettes. And to scale above 2000 L monoclonal antibody production for aggregate removal, Sartobind® phenyl cassettes would be the right choice.

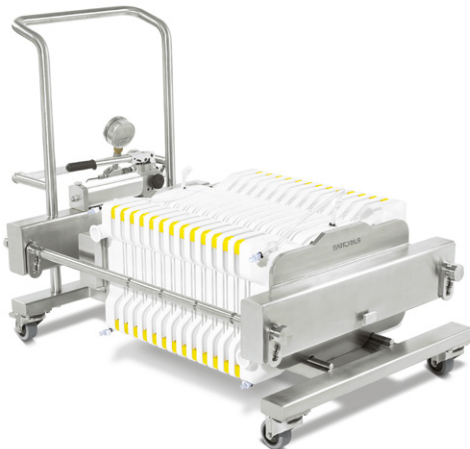


Double-process holder with Sartoclear® filters

Flexibility

The filter holders offer maximum flexibility for your process. Between one and 12 Sartoclear® depth filters can be installed in a process layer of the process or double-process filter holder. This enables bioreactors with volumes of between 50 and 2000 liters to be harvested with a double-process holder. In the pilot filter holder between one and 13 Sartobind® cassettes can be set up.

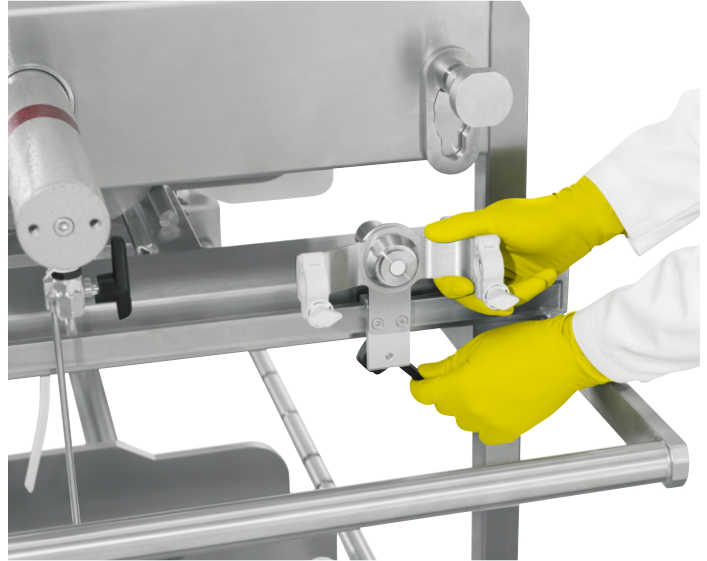
The separation plate allows you to install common filter, membrane chromatography (adsorber) cassette combinations in series or parallel on one layer.



Pilot filter holder with Sartobind® membrane chromatography cassettes

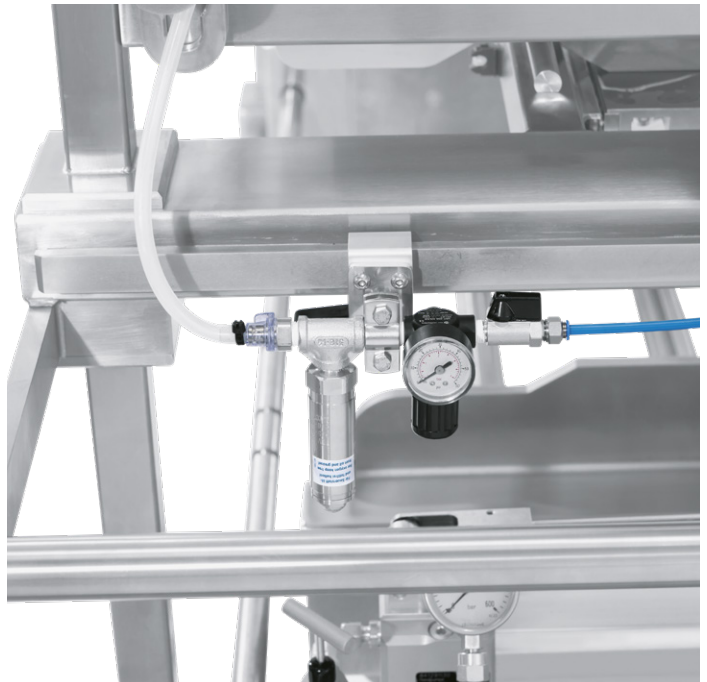
Easy Handling

The filter holders are made for use in a biopharmaceutical environment. The systems are designed for easy cleaning and ergonomic handling. The installation of varying numbers of filters is intuitive, fast and easy.

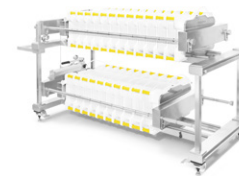
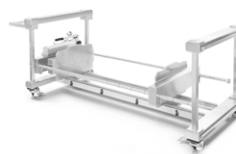


User Friendly

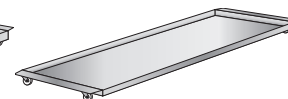
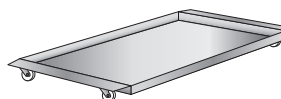
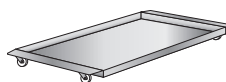
The filter holders are equipped with easily accessible mounting bars for quick and secure installation of commonly used accessories such as pressure sensors, pressure safety devices and bioburden reduction filters.



Technical Data



	Pilot Filter Holder	Process Filter Holder	Double Process Filter Holder
Layers	1	1	2
Capacity for 0.8 m ² Sartoclear® double-layer filters	1-6	1-12	1-24
Capacity for 1.0 m ² Sartoclear® single-layer filters	1-6	1-12	1-24
Capacity for 0.8 L or 1.6 L Sartobind® cassettes	1-13	Request	Request
Dimensions H x W x D [cm]	96 x 133 x 79	91 x 230 x 93	122 x 230 x 93
Weight	160 Kg	500 Kg	950 Kg
Max. total weight	290 Kg	780 Kg	1430 Kg
Materials of construction	Stainless steel	Stainless steel	Stainless steel
CE pressure equipment directive	Designed according to the requirements of the pressure directive 97/23/EC		



Drip pan	Pilot Filter Holder	Process Filter Holder	Double Process Filter Holder
Dimensions H x W x D [cm]	4 x 110 x 67	4 x 184 x 70	4 x 184 x 70
Materials of construction	Stainless Steel	Stainless Steel	Stainless Steel

Order Information

	Description	Order No.
Holders		
Pilot filter holder	For Sartoclear® depth filters or Sartobind® cassettes	2ZGL--0005
Process filter holder	For Sartoclear® depth filters or Sartobind® cassettes	2ZGL--0006
Double process filter holder	For Sartoclear® depth filters or Sartobind® cassettes	2ZGL--0007
Accessories		
Drip pan	For pilot filter holder For easy cleaning of spilled fluids	2ZGL--0008
	For process filter holder and double process filter holder For easy cleaning of spilled fluids	2ZGL--0015
Separation plate	For serial or parallel filtration on one layer	2ZGL--0013
Pressure safety device	Prevents excessive pressure during compressed air draining of the filter cassettes	2ZGL--0014




The clamping force is adjusted easily with the manually operated hydraulic pump

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com