



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

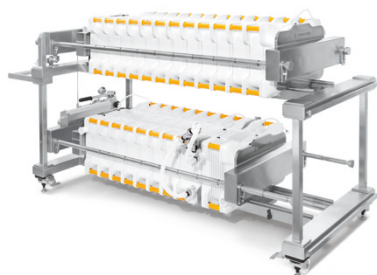
Sartoclear® Depth Filters and Sartobind® Cassettes are two complementing technologies using the same holder and manifold plates. Use Sartoclear® Depth Filters for applications with moderate solids content (see table for more application guidance).

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

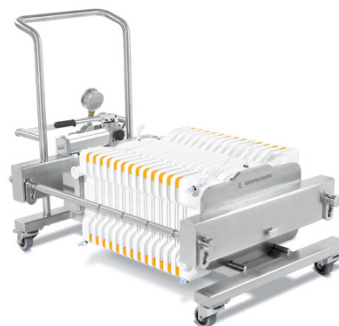
Flexibility

The filter holders offer maximum flexibility for your process. Between 1 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 1 and 13 Sartobind® cassettes can be set up.

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



Double Process Holder with Sartoclear® filters



Pilot filter holder with Sartobind® membrane chromatography cassettes

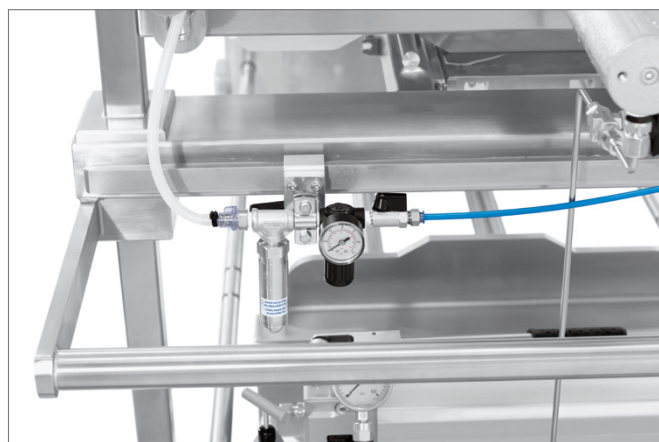
Easy handling

The filter holders are made for the 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.

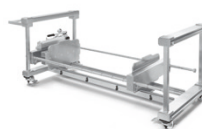


User friendly

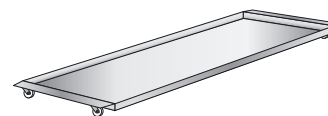
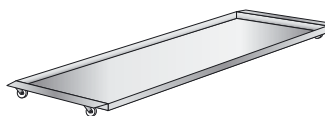
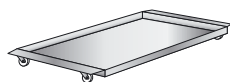
The filter holders are equipped with easy accessible mounting bars for the quick and secure installation of commonly used accessories such as pressure sensors, pressure safety devices and bioburden reduction filters.



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 × W × D (cm)	96 × 133 × 79	91 × 230 × 93	122 × 230 × 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 the requirements of the Pressure Directive 97/23/EC		



	Pilot Filter Holder	Process Filter Holder	Double Process Filter Holder
Drip pan			
Dimensions H × W × D (cm)	4 × 110 × 67	4 × 184 × 70	4 × 184 × 70
Materials of construction	Stainless Steel	Stainless Steel	Stainless Steel

Ordering 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

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49.551.308.0

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5 – 7
34302 Guxhagen
Phone +49.5665.407.0

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques – CS 91051
13781 Aubagne Cedex
Phone +33.442.845600

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex
Phone +33.442.845600

Austria

Sartorius Stedim Austria GmbH
Modectcenterstrasse 22
1030 Vienna
Phone +43.1.7965763.18

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.80

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi
Phone +36.23.457.227

Italy

Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41

Netherlands

Sartorius Stedim Netherlands B.V.
Phone +31.30.60.25.080
filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.647.38.40

Russian Federation

LLC "Sartorius Stedim RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone +34.913.586.098

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen
Phone +41.52.354.36.36

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ
Phone +44.1372.737159

Ukraine

LLC "Sartorius Stedim RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Africa

Republic of South Africa

Sartorius South Africa (Pty) Ltd
Unit 4, Alphen Square South
853 16th Road
Midrand 1685, RSA
Phone +27.11.315.5444

Americas

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1.800.368.7178

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Beijing Branch Office
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Guangzhou Branch Office
Room 1105
Xing Guang Ying Jing Building
No. 119, Shui Yin Road
Yue Xiu District, Guangzhou 510075
Phone +86.20.3836.4193

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.4331.4300

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700



► www.sartorius-stedim.com