

Tangential Flow Filtration Filter Holders

Predictable Performance from Scaling Studies up to Commercial Manufacturing Simplifying Progress

SARTURIUS

Sartorius Stedim Biotech at a Glance

At Sartorius Stedim Biotech, we work together to achieve one primary goal: to provide our customers with the best solutions for their process needs. That's why we deliver across the board. Our products and services cover the biopharmaceutical industry's entire process chain – from development and quality assurance to production. We are your total solution provider and experts in single-use technologies and added-value services.

With Sartorius Stedim Biotech, you have gained a partner with extensive engineering capabilities in biotechnology and pharmaceutical processes. We will provide you with highperformance membranes that are integral elements in your biopharmaceutical manufacturing process.

On top of our cutting-edge skills in developing membranes and designing cassettes, our core area of expertise is in engineering the highest quality Tangential Flow Filtration (TFF) filters and systems.

These integrated capabilities extend from laboratory to pilot scale, right on up to clinical manufacturing. We offer automated production systems that enable bioprocess manufacturers to integrate upstream and downstream processes. In any case, you can be sure that you will be benefiting from high-quality products and services.





Sartorius Stedim Biotech TFF - Simply Convincing

Your requirement is to efficiently process high-value products with the maximum possible yield. High-performance membranes are an integral part of downstream processing applications, such as concentration and diafiltration of proteins, vaccines, viruses, antibodies, oligonucleotides, endotoxin removal, etc. TFF is also an efficient method for cell harvesting and clarification processes.

Every single step in a biopharmaceutical process requires high-performance filtration units to ensure maximum product yield in all manufacturing phases. Highvalue biological compounds can be produced only within precisely defined environments for consistent and robust, downstream processing. This is where Sartorius Stedim Biotech TFF products and technologies excel.

The unique design within the TFF cassette family provides end users with predictable performance for scaling studies. Each cassette has the same hydrodynamic flow path geometry and identical materials of construction. The cassettes are self-sealing and feature low protein binding. The vertical positioning of the cassettes in the holder enables complete drainage of retentate and permeate channels by gravity during recovery procedures. With our products and services, we offer precisely defined quality standards, a wide choice of performance specifications, and incomparably high yield rates.

The Sartorius Stedim Biotech TFF cassette family is scalable from lab to process.

TFF Process-Scale Filter Holders for Maximum Reliability

The family of Sartorius TFF filter holders features the latest advances in TFF design for full flexibility.

- Self-draining and self-venting thanks to vertical positioning
- Large retentate channels for low pressure drop
- Easy installation
- Low hold-up volume
- Monolithic design
- All ports located on the product plate

This unique design within the Sartocon® family provides you with predictable performance from process volumes down to scaling studies or in daily lab work.

Sartocon® Slice 200 Holder

The Sartocon® Slice 200 stainless steel holder is optimized for use with up to two Slice 200 TFF cassettes (max. 400 cm²). It is designed for low-volume applications from 20 mL to 5 L. The Sartocon® Slice 200 holder uses female stainless steel Luer lock connectors. This ensures safe and reliable connections to ancillary equipment. The Slice 200 holder is the perfect tool for any low-volume TFF application.

The Sartocon® Slice 200 cassettes have the same flow channel dimensions as do Sartocon® process filtration cassettes. Therefore, it is ideal for your scaling studies.

Target Applications

- Scaling studies
- Membrane screening studies
- Product development
- Preclinical trials



Sartocon® Slice Holder

The Sartocon® Slice stainless steel holder has been designed for low-to medium-volume applications from 300 mL to 100 L and optimized for use with up to five Sartocon® Slice TFF cassettes (0.1–0.5 m²). The filter holder can be sterilized by autoclaving or inline steaming if pressure compensation tools are used.

Target Applications

- Process development
- Preclinical trials
- Clinical trials
- Scaling studies
- Pilot-scale and small-volume production

This unique design within the Sartocon® family provides you with predictable performance all the way from process volumes down to scaling studies or daily labscale procedures.



Outstanding Design for Full Flexibility

- Self-draining and self-venting due to vertical positioning
- Large, oval retentate channels for low pressure drop
- Easy installation
- Low hold-up volume
- Monolithic design
- All ports located on the product plate
- Hard-piped installation with low hold-up volume

Target Applications

- Process development
- Preclinical trials
- Clinical trials
- Pilot lots
- Production

Our process filter holders are available in a choice of versions compliant with DIN, ISO or the Japanese Industrial Standard. All holders are manufactured using high-grade 316 L stainless steel. For maximum flexibility, you can order any process filter holder with a mirror finish.

Sartocon® 2 Plus

The Sartocon® 2 plus AISI 316 L stainless steel holder (1.4435) has been optimized for use with Sartocon® production-scale TFF cassettes of up to 7 m². It is suitable for applications from 30 L to >2,000 L process volumes. Its easy-to use design and high quality materials make this holder the perfect tool.



Sartoflow® 10

The Sartoflow® 10 holder (1.4435 stainless steel) has been optimized for use with Sartocon® production-scale TFF cassettes featuring up to 7 m² of membrane area. This holder is suitable for applications from 30 L to >2,000 L volumes.



This holder is designed for thermal sterilization in an autoclave.



An easy-to-use design and high-quality materials make this holder the perfect tool.

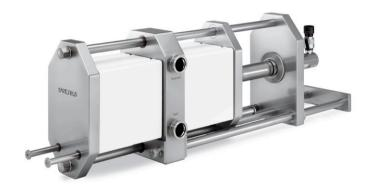


Sartoflow® 20

The Sartoflow® 20 AISI 316 L stainless steel holder (1.4435) has been optimized for use with up to 14 m² Sartocon® TFF cassettes. It is designed for volumes ranging from 200 L to >2,500 L.



The clamping pressure can be controlled by a manual pump or an automatic hydraulic pump.



Sartoflow® 40 | 60 | 80

To match the advanced level of our membrane development and cassette designing capabilities, our core area of expertise is in engineering the highest quality TFF filters, holders and systems. This extends from lab to pilot scale, to clinical manufacturing, all the way to fully automated production systems.

The stacked holders have been designed following the proven principle of the Sartoflow® 20 filter holder with an enlarged filter plate for the integration of feed, retentate and permeate piping. Depending on specific process requirements, the piping can be mirror inverted. Sartoflow® 40 | 60 | 80 holding devices offer a combination of the smallest footprint possible, safety and operator convenience.

Like all other Sartocon® and Sartoflow® filter holders from Sartorius Stedim Biotech, the stacked filter holder can be drained backwards by gravity. Each level of a Sartoflow® 40 | 60 | 80 holder can hold up to 14 m² of Sartocon® or Sartocube® cassettes, adding to a total filter area of 28 m², 42 m² or 56 m². The family of Sartorius TFF filter holders features the latest advances in TFF design for full flexibility. Thanks to vertical positioning of our cassettes inside the filter holders, the feed flow easily evacuates air from the system. Sartorius Stedim Biotech offers the perfect equipment solution for single-use or reusable TFF applications, such as concentration and diafiltration of proteins, vaccines, viruses, antibodies and oligonucleotides.

Sartoflow® 40



Sartoflow® 60



Sartoflow® 80



Ordering Information

TFF Accessories

Product	Order number	Description			
Adapter plates	17546113-C	Sartocon® Single-use Adapter (PP-DWST)			
	17521114-C	Sartocon® Slice Single-use Adapter (PP-DWST)			
	17546116-C	Sartocon® Single Use Adapter-Set (PP-DWST)			
	175Z40GDB-0001	Sartocon® Adapter (316 L) - Round to oval ports			
	302020A-C	Sartocon® Adapter (PP-DWST) – Round to oval ports			
	17521029-C	Sartocon® Slice Adapter (PP-DWST) – Round to oval ports			
Silicone gaskets	17546114	Gasket for Sartocon® Adapter plate (316 L)			
	17521030C	Sartocon® Slice oval hole gasket			
	17521031C	Sartocon® Slice round hole gasket			
	17546112C	Sartocon® oval hole gasket			
	17546115	Sartocon® round hole gasket			
Flush plates	302030C-C	Sartocon® Flush plate (PP-DWST & for cleaning EPDM O-ring) for feed and permeate flushing			
	305030C-C	Sartocon® Slice Flush plate (PP-DWST & EPDM O-ring) for feed and permeate flushing			
Dummy module plates,	302020D-C	Sartocon® Dummy-Module (PP-DWST) 20 mm – equivalent of 1 cassette			
without flow channels	302040D-C	Sartocon® Dummy-Module (PP-DWST) 40 mm – equivalent of 2 cassettes			
	302100D-C	Sartocon® Dummy-Module (PP-DWST) 100 mm – equivalent of 5 cassettes			
Supporting materials	1ZJO0001	Sartoflow® Heat Shield			
	U1-SCUBE-WEDGE	Sartoflow® Wedge for standard Sartoflow® 10 and 20 holders			
Threaded spacers for	17000309	Nominal length = 100 mm Effective length = 85 mm			
Sartoflow® 10 20	17000314	Nominal length = 65 mm Effective length = 50 mm			
Trolley	179SFG-18	Trolley for Sartoflow® 20 with drip tray for storage of components (e.g., pump, hoses); with two wheels and two swivel castors with brake			
Hand hydraulic set	1ZW0004	Hand hydraulic set for Sartoflow® 10 for SF 10 or SF 20 Sartoflow® 20: Hand hydraulic pump, pressure gauge, pressure hose, needle valve			

Filter Holders and Filtration Sets

Product	Description	Order number	
Slice 200 Sartocon® Slice 200 filter holder (316 L) – 200 – 400 cm²		1752501	
3.5 m² SU Filter Holder	3.5 m² SU Filter Holder with Handhydraulic	BB-34164022	
Slice microfiltration set	Sartocon® Slice Microfiltration set (316 L) – 0.1 – 0.5 m², filter holder, pressure gauges, valves, clamps, permeate manifold, torque wrench	17521105	
Slice ultrafiltration set	Sartocon® Slice Ultrafiltration set (316 L) – 0.1 – 0.5 m² with filter holder, pressure gauges, valves, clamps, permeate manifold, torque wrench	17521106	
Microfiltration set Sartocon® 2 Plus	Sartocon® Microfiltration set (316 L) – 7 m², filter holder, pressure gauges, valves, clamps, permeate manifold, torque wrench	17546201	
Ultrafiltration set Sartocon® 2 Plus Sartocon® Ultrafiltration set (316 L) – 7 m², filter holder, pressure gauge, valves, clamps, permeate manifold, torque wrench		17546202	

Sartocon® 2 Plus, AISI 316 L (1.4435)

Order number	Standard	F R (mm)	P (mm)	Mirror-inverted
17546DD260160E	DIN	26	16	_
17546DZ221157E	Zoll	22.1	15.7	
17546DI297181E	ISO	29.7	18.1	
17546DJ230230E	JIS-G	23	23	
17546DD260160EM	DIN	26	16	
17546DZ221157EM	Zoll	22.1	15.7	
17546DI297181EM	ISO	29.7	18.1	
17546DJ230230EM	JIS-G	23	23	

Sartoflow® 10, AISI 316 L (1.4435)

Order number	Standard	F R (mm)	P (mm)	Mirror-inverted
179-6DD260160E	DIN	26	16	
179-6DZ221157E	Zoll	22.1	15.7	
179-6DI297181E	ISO	29.7	18.1	
179-6DJ230230E	JIS-G	23	23	
179-6DD260160EM	DIN	26	16	
179-6DZ221157EM	Zoll	22.1	15.7	
179-6DI297181EM	ISO	29.7	18.1	
179-6DJ230230EM	JIS-G	23	23	

Sartoflow® 20, AISI 316 L (1.4435)

Order number	Standard	F R (mm)	P (mm)	Mirror-inverted
179-2DD380260E	DIN	38	26	
179-2DZ348221E	Zoll	34.8	22.1	
179-2DI384237E	ISO	38.4	23.7	
179-2DJ357230E	JIS-G	35.7	23	
179-2DD380260EM	DIN	38	26	
179-2DZ348221EM	Zoll	34.8	22.1	
179-2DI384237EM	ISO	38.4	23.7	
179-2DJ357230EM	JIS-G	35.7	23	
179-6DJ230230EM	JIS-G	23	23	

F = Feed R = Retentate P = Permeate

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



Representation Formation, visit

www.sartorius.com