

# Sartoflow<sup>®</sup> 10 Stainless Steel Holder

Hydraulic Filter Holder for  
Sartocon<sup>®</sup> Cassettes



## Description

### Sartoflow<sup>®</sup> 10 Holder

The Sartoflow<sup>®</sup> 10 stainless steel holder is optimized for use with up to 10 Sartocon<sup>®</sup> production scale Tangential Flow Filtration (TFF) cassettes or two Sartocubes<sup>®</sup> (max. 7 m<sup>2</sup>). It is designed for applications from 30 L to 1,000 L volume.

Target use:

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

Sartocon<sup>®</sup> cassettes with a filter area of up to 0.7 m<sup>2</sup> each, are available with Hydrosart<sup>®</sup> and Polyethersulfone membrane types.

Sartoflow® 10 holders cover the whole range of pharmaceutical and biotechnological TFF applications like concentration and diafiltration of proteins, vaccines, viruses, antibodies, oligo nucleotides, endotoxin removal etc. The holder is also suitable for cell harvesting and clarification processes. The family of Sartorius TFF holders feature the latest advances in TFF design:

- Vertical orientation of cassettes
- Self-draining systems
- In-line steamable
- No welds
- Minimized hold-up volume

The Sartoflow® 10 filter holder is part of a modular micro- and ultrafiltration system and fits up to 10 Sartococon® filter cassettes. It is designed for process development and pilot scale production in biopharmaceutical applications.

The Sartococon® cassettes are placed between the filter and clamping plate, without a need for additional gaskets. It is not necessary to remove the clamping plate for installation of the cassettes. Place the cassettes on the guide rods and close the holder with the manual driven oil hydraulic pump. The TFF holder and the hydraulic pump are optionally mounted on a stainless steel skid for easy handling.

The clamping pressure can be controlled by an integrated oil pressure gauge. This allows very accurate and reproducible clamping conditions. Defined clamping conditions are essential, especially for steaming in place applications, but it is also very valuable when cassettes have to be changed frequently. The Sartoflow® 10 holder offers high performance production scale technology for every process development and small scale production facility.

The vertical positioning of the cassettes allows complete draining of retentate and permeate channels by gravity during harvesting and steaming procedures.

The Sartoflow® 10 flow distribution plate is machined from a single piece of stainless steel. Its sanitary tri-clamp connections are not welded but machined from the same plate. This eliminates potential problems associated with welds and ensures a long service life and safe and reliable connection to other equipment. The feed and retentate connectors are located on one side of the plate. This allows easy integration into any skid design and ensures a compact system with low minimum working volume.

Connectors are situated to eliminate potential air locks and to ensure thorough cleaning and total product recovery.

Features	Benefits
Vertical position	Self-draining Self-venting
Tri-clamp connectors	Safe fit of accessories
No welds	No corrosion
Steamable	No cross contamination
Fits up to 10 cassettes or two Sartocubes®	Up to 1,000 L working volume
Hydraulic closure	Precise and reproducible clamping conditions

## Technical Data

### Specifications

Materials of Construction	
Filter and clamping plate	316 L (1.4435) stainless steel, grinded
Connectors	Sanitary Tri Clamp flange (DN25)
Other components	316 L stainless steel

Surface finish, grinded	Ra ≤ 0.4 µm
Closure	Lukas hydraulic ram ACM 10/150
Oil pressure gauge	0-400 bar   0-6000 psi
Pressure limit	430 bar
Maximum number of cassettes	10 Sartococon® cassettes
Dimensions (L×W×H)	680 × 420 × 508 mm
Weight	Approx. 70 kg

## Accessories and Order Information

Handhydraulic set for SF 10 or SF 20 Set: Handhydraulic pump manometer, pressure hose, needle valve	1ZW---0004	Sartocon® dummy modules PP-DWST 20 mm 40 mm 100 mm	302020D-C 302040D-C 302100D-C
Pressure gauge, 0-6 bar, 1-1/2"	17546---003	Sartoflow® Wedge	U1-SCUBE-WEDGE
Diaphragm valve, 1-1/2"	17546---005	Sartoflow® Heatshield	1ZJO--0001
Silicone gasket (FDA), 1-1/2"	17546---012	Threaded spacer nominal length 100mm	17000---309
Clamp for 1-1/2" Tri-clamp	SB-18-00-0194	Threaded spacer nominal length 65 mm	17000---314
Sartocon® manifold for permeate outlet	17546---016		
Hose connector to 1-1/2" Tri-clamp flange	17546---018		
Sartocon® adapter plate (round to oval adapter)	302020A		
Sartocon® single-use adapter plates	17546---113-C		
Sartocon® flush plate PP-DWST	302030C		
		<b>Included with Filter Holder</b>	<b>Product Number</b>
		Sartoflow® Wedge	U1-SCUBE-WEDGE
		Sartoflow® Heatshield	1ZJO--0001
		Threaded spacer nominal length 100mm	17000---309
		Threaded spacer nominal length 65 mm	17000---314
		Dummy module (100 mm)	302100D-C
			<b>Units</b>
			2
			1
			2
			1
			1

## Sartoflow® 10, 316 L (1.4435)


Order Number	Norm	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	x
179-6DZ221157EM	Zoll	22.1	15.7	x
179-6DI297181EM	ISO	29.7	18.1	x
179-6DJ230230EM	JIS-G	23	23	x

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