



SARTOFLOW[®] 10 Stainless Steel Holder

Hydraulic Crossflow holder for Sartoco[®] cassettes

Sartorius Stedim Biotech Crossflow Holder



Description

SARTOFLOW[®] 10 Holder

The SARTOFLOW[®] 10 stainless steel holder is optimized for the use with up to ten Sartoco[®] production scale Crossflow cassettes or two Sartocubes[®] (max. 7 m²). It is designed for applications from 30 l to 1000 l volume.

Target use:

- process development
- preclinical trials
- clinical trials
- pilot lots
- production

Sartoco[®] cassettes with a filter area of up to 0.7 m² each, are available with Hydrosart[®], and Polyethersulfone, membrane types.

SARTOFLOW[®] 10 systems cover the whole range of pharmaceutical and biotechnological crossflow applications like concentration and diafiltration of proteins, vaccines, viruses, antibodies, oligo nucleotides, endotoxin removal etc. The system is also suitable for cell harvesting and clarification processes. The family of Sartorius Stedim Biotech Crossflow holders feature the latest advances in crossflow design:

- vertical orientation of cassettes
- self draining systems
- inline 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 ten Sartoco[®] filter cassettes. It is designed for process development and pilot scale production in biopharmaceutical applications.

The Sartoco[®] cassettes are placed between filter and clamping plate without a need for additional gaskets. It's not even necessary to remove the clamping plate for installation of the cassettes. Just place the cassettes on the guide rods and close the holder with the manual driven oil hydraulic pump. The crossflow 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. It's sanitary tri-clamp connections are not welded but machined from the same plate. This eliminates potentially problems associated with welds and ensures 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 assures a compact system with low minimum working volume.

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

Feature	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 2 Sartocubes [®]	up to 1000 l working volume
Hydraulic closure	precise and repro- ducible clamping conditions

Specifications

Materials of Construction

Filter and Clamping plate	316L (1.4435) stainless steel, grinded
Connectors	sanitary Tri Clamp flange (DN25)
Other components	316L stainless steel

Technical Data

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 Sartoco [®] cassettes
Dimensions (L×W×H)	680×420×508 mm
Weight	approx. 70 kg

Accessories and Ordering Informations

Handhydraulic Set for SF 10 or SF 20 Set: Handhydraulic pump manometer, pressure hose, needle valve	1ZW---0004
Pressure gauge, 0–6 bar, 1–1 1/2"	17546---003
Diaphragm valve, 1–1 1/2"	17546---005
Silicone gasket (FDA), 1–1 1/2"	17546---012
Clamp for 1–1 1/2" Tri-clamp	17033
Sartoco [®] manifold for permeate outlet	17546---016
Hose connector to 1–1 1/2" Tri-clamp flange	17546---018
Sartoco [®] Adapterplate (round to oval adapter)	302020A-C
Sartoco [®] Single Use Adapter Plates	17546---113-C
Sartoco [®] Flush Plate PP-DWST	302030C-C
Sartoco [®] Dummy Modules PP-DWST	
20 mm	302020D-C
40 mm	302040D-C
100 mm	302100D-C

Ultrafiltration

Membrane Material	MWCO	Order Numbers Sartoco [®]	Order Numbers Sartocube [®]
Hydrosart [®]	2 kD	3021441906E--SW	3021441930E-BSW
	5 kD	3021442906E--SW	3021442930E-BSW
	10 kD	3021443906E--SW	3021443930E-BSW
	30 kD	3021445906E--SW	3021445930E-BSW
	100 kD	3021446806E--SW	3021446830E-BSW
PESU	1 kD	3021460907E--SW	n.a.
	5 kD	3021462907E--SW	n.a.
	8 kD	3021463407E--SW	n.a.
	10 kD	3021463907E--SW	3021463935E-BSW
	Albumin	30214AL07K--SW	n.a.
	30 kD	3021465907E--SW	3021463935E-BSW
	50 kD	3021465007E--SW	3021463035E-BSW
	100 kD	3021466807E--SW	3021466835E-BSW
	300 kD	3021467907E--SW	3021467935E-BSW

Microfiltration

Membrane Material	Pore size	Order Numbers Sartoco [®]	Order Numbers Sartocube [®]
Hydrosart [®]	0.2 µm	3021860706W--SW	3021860730W-BSW
	0.45 µm	3021860606W--SW	3021860630W-BSW
PESU	0.1 µm	3021545806W--SW	3021545830W-BSW

SARTOFLOW[®] 10, 316L (1.4435)

Order number	Norm	F/R [mm]	P [mm]	e-polished	mirror - inverted
179-6DD260160	DIN	26	16		
179-6DD260160E	DIN	26	16	×	
179-6DZ221157	Zoll	22.1	15.7		
179-6DZ221157E	Zoll	22.1	15.7	×	
179-6DI297181	ISO	29.7	18.1		
179-6DI297181E	ISO	29.7	18.1	×	
179-6DJ230230	JIS-G	23	23		
179-6DJ230230E	JIS-G	23	23	×	
179-6DD260160-M	DIN	26	16		×
179-6DD260160EM	DIN	26	16	×	×
179-6DZ221157-M	Zoll	22.1	15.7		×
179-6DZ221157EM	Zoll	22.1	15.7	×	×
179-6DI297181-M	ISO	29.7	18.1		×
179-6DI297181EM	ISO	29.7	18.1	×	×
179-6DJ230230-M	JIS-G	23	23		×
179-6DJ230230EM	JIS-G	23	23	×	×

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH. | W
Publication No.: SPC2050-e12034
Order No.: 85030-530-82
Ver. 03 | 2012