



# Sartocon® 2 plus Stainless Steel Holder

Crossflow holder for Sartocon® cassettes

Sartorius Stedim Biotech Crossflow Holder



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.

The Sartocon® 2 plus device is available in an electro and a mechanical polished version.

Feature	Benefits
Cassettes in vertical position	self draining
Tri-clamp connectors	safe fit of accessories
No welds	no corrosion
Steamable	no cross contamination

Fits up to ten Sartocon® cassettes	high flexibility for up to 1000 l working volume
------------------------------------	--

## Specifications

### Materials of Construction

Filter and Clamping plate	316L stainless steel
Connectors	1"–1 ½" sanitary Tri Clamp flange (DN25)
Nuts for tie rods	nickel coated bronze
Other components	316L Stainless steel

## Description

### Sartocon® 2 plus

The Sartocon® 2 plus stainless steel holder is optimized for the use with up to ten Sartocon® production scale Crossflow cassettes or two Sartocubes® (max. 7 m<sup>2</sup>). It is designed for applications from 30 l to 1000 l volume.

Target use:

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

Sartocon® cassettes with a filter area of up to 0.7 m<sup>2</sup> each, are available with Polyether-sulfone and Hydrosart membrans.

The Sartocon® 2 plus system covers 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
- autoclavable and inline steamable
- no welds
- minimized hold up volume

The Sartocon® 2 plus holding device is part of a modular micro- and ultrafiltration system and fits up to ten Sartocon® filter cassettes. It is designed for process development and pilot scale production in pharmaceutical and biotechnology applications. Easy handling design and high quality materials make this holder the perfect tool.

The Sartocon® 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 a torque wrench.

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

Thermal sterilisation in an autoclave or steaming in place even with cassettes installed is possible.

The Sartocon® 2 plus's flow distribution plate is machined from a single piece of stainless steel. It's sanitary 1 ½" 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.

**Technical Data**

Surface finish	Ra ≤ 0.6 µm
Maximum number of cassettes	10 Sartococon® cassettes
Dimensions (L×W×H)	371×270×348 mm
Weight	approx. 50 kg

**Ordering Informations and Accessories**

Sartococon® 2 plus microfiltration set	17546---201
Sartococon® 2 plus ultrafiltration set	17546---202
Torque wrench for Sartococon® 2 plus holder	17128
Socket, 27 mm	6986135
Pressure gauge, 0-6 bar, 1-1 1/2" autoclavable	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
Sartococon® 2 plus manifold for permeate outlet	17546---016
Hose connector to 1-1 1/2" Tri Clamp flange	17546---018
Sartococon® Adapterplate (round to oval adapter)	302020A-C
Sartococon® Single Use Adapter Plates	17546---113-C
Sartococon® Flush Plate PP-DWST	302030C-C
Sartococon® Dummy Modules PP-DWST	
20 mm	302020D-C
40 mm	302040D-C
100 mm	302100D-C

**Ultrafiltration**

Membrane Material	MWCO	Order Numbers Sartococon®	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	302146AL07K--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 Sartococon®	Order Numbers Sartocube®
Hydrosart®	0.2 µm	3021860706W--SW	3021860730W-BSW
	0.45 µm	3021860606W--SW	3021860630W-BSW
Polyethersulfone (PESU)	0.1 µm	3021545806W--SW	3021545830W-BSW

**Sartocon® 2 plus, 1.4404**

SAP	Norm	F   R		P	e-polished	mirror-inverted
		[mm]	[mm]			
17546ID260160	DIN	26	16			
17546ID260160E	DIN	26	16	×		
17546IZ221157	Zoll	22.1	15.7			
17546IZ221157E	Zoll	22.1	15.7	×		
17546II297181	ISO	29.7	18.1			
17546II297181E	ISO	29.7	18.1	×		
17546IJ230230	JIS-G	23	23			
17546IJ230230E	JIS-G	23	23	×		
17546ID260160-M	DIN	26	16		×	
17546ID260160EM	DIN	26	16	×	×	
17546IZ221157-M	Zoll	22.1	15.7		×	
17546IZ221157EM	Zoll	22.1	15.7	×	×	
17546II297181-M	ISO	29.7	18.1		×	
17546II297181EM	ISO	29.7	18.1	×	×	
17546IJ230230-M	JIS-G	23	23		×	
17546IJ230230EM	JIS-G	23	23	×	×	

**Sartocon® 2 plus, 1.4435**

SAP	Norm	F   R		P	e-polished	mirror-inverted
		[mm]	[mm]			
17546DD260160	DIN	26	16			
17546DD260160E	DIN	26	16	×		
17546DZ221157	Zoll	22.1	15.7			
17546DZ221157E	Zoll	22.1	15.7	×		
17546DI297181	ISO	29.7	18.1			
17546DI297181E	ISO	29.7	18.1	×		
17546DJ230230	JIS-G	23	23			
17546DJ230230E	JIS-G	23	23	×		
17546DD260160-M	DIN	26	16		×	
17546DD260160EM	DIN	26	16	×	×	
17546DZ221157-M	Zoll	22.1	15.7		×	
17546DZ221157EM	Zoll	22.1	15.7	×	×	
17546DI297181-M	ISO	29.7	18.1		×	
17546DI297181EM	ISO	29.7	18.1	×	×	
17546DJ230230-M	JIS-G	23	23		×	
17546DJ230230EM	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](http://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.: SPC2052-e12034  
Order No.: 85030-530-84  
Ver. 03 | 2012