

Sartocon[®] Slice Stainless Steel Holder

Low hold-up Volume
Tangential Flow Filtration
Holder For Sartocon[®] Slice
Cassettes



Product Information

The Sartocon[®] Slice stainless steel holder is optimized for the use with up to five Slice Tangential Flow Filtration (TFF) cassettes (0.1–0.5 m²). It is designed for low to medium volume applications from 1–100 L.

Target Use

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

Features and Benefits

- Cassettes in vertical position
- Unique optimized In- and Outlets
- Fits up to five cassettes
- Low hold up volume
- Adjustable feet
- Self draining | Self venting
- No air lock
- Provides process flexibility
- Minimized working volume
- Secure stand

Sartocon® Slice cassettes have a surface area of 0.1 m² each and are available with Polyethersulfone and Hydrosart® membrane types.

The filterholder can be thermally sterilized by autoclaving or inline steaming if the pressure compensation tools are used. Typical applications for Sartocon® Slice 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 system is suitable for cell harvesting and clarification processes.

The Sartocon® Slice filterholder features the latest advances in TFF design:

- Vertical orientation
- Self draining system
- Low minimal hold-up volume
- Monolithic design
- Large retentate channels
- cGMP construction and operation

The Sartocon® Slice cassette (filter area: 0.1 m²) has the same hydrodynamic flow path design and identical materials like the larger Sartocon® cassette (Hydrosart® 0.6 m², PESU 0.7 m²).

This unique design within the Sartocon® family provides the user with predictable performance for scaling studies.

The Sartocon® Slice holder completes the Sartocon® Slice cassette in an ideal way and gives the perfect tool for every low volume TFF application. The Sartocon® Slice holder has sanitary tri-clamp connections. This ensures a safe and reliable connection to other equipment. The feed and retentate ports as well as the two filtrate ports are located on one side. This allows fixed piping and a compact system with low minimum working volume. The adjustable feet guarantees a firm stand of the holder on the bench.

Port locations are situated to avoid air locks and to ensure proper cleaning and product recovery.

Technical Specifications

Number of cassettes	1–5
Surface area	0.1–0.5 m ²
Material	AISI 316 L stainless steel (German standard 1.4435), mirror polished
Surface finish Product contact	Ra ≤ 0.4 µm
Support plate	AISI 316 L stainless steel
Port connection	<ul style="list-style-type: none"> ▪ Outer diameter 25 mm (TC25) ▪ Inner diameter 10 mm (DN10) ▪ DIN11850, Reihe 2
Max. operating pressure at 20°C	4 bar (58 psi)

Ordering Information

Product Name	Order Number
Sartocon® Slice Holder	17521---002
Pressure compensation tools	17521---028
Sartocon® Slice Flush Plate PP-DWST	305030C-C
Sartocon® Slice Adapter Plates	17521---029-C 17521---920-C

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 more information, visit**
www.sartorius.com