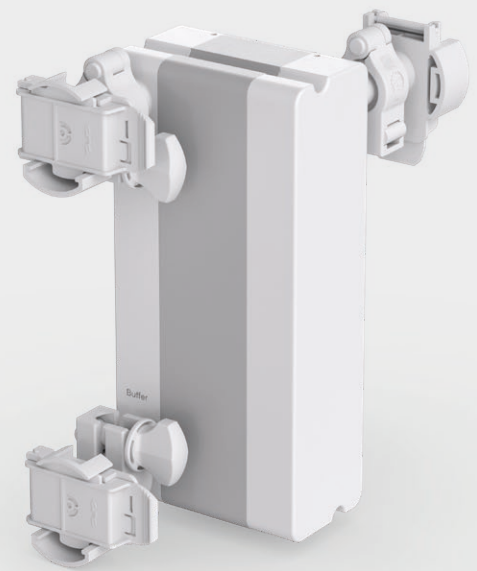


Sartocon® SP

Self-contained Tangential
Flow Filtration Cassettes for
Diafiltration in a Single Pass



Product Information

Sartocon® SP (Single Pass) cassettes are designed for continuous diafiltration with low buffer consumption that matches conventional diafiltration performance in a single pass.

Target Use

- Ideally suited for continuous diafiltration of monoclonal antibodies
- Process development for continuous and intensified processing
- Preclinical | clinical trials

Sartocon® SP cassettes are available with Hydrosart® membrane format in 30 kDa molecular weight cut-off and feature a unique three chamber design to facilitate highly effective, continuous diafiltration.

In combination with the Sartoflow® SP system, it's possible to achieve a more productive and cost effective mAb concentration and buffer exchange processing tool.

Hydrosart® Membrane

Hydrosart® stabilized cellulose is a high performance membrane that has been optimized for biopharmaceutical process applications. For long term continuous diafiltration processing the key benefits of Hydrosart® are low protein binding, non-fouling and easy and reproducible regeneration are critical parameters in particular when joined with unparalleled diafiltration performance.

Hydrosart® is stable across a broad pH range, is extremely hydrophilic, making it non-protein binding, and virtually non-fouling. Membrane regeneration, and depyrogenation are accomplished by using (1N) NaOH at elevated temperatures as required.

Product Profile

Sartocon® SP membrane cassettes exhibit minimal adsorption properties of proteins, viruses and other biological materials

Feature	Benefits
Single pass diafiltration flow path	Perfectly suited to continuous diafiltration
Three internal flow channels	Ultra efficient use of diafiltration buffer
Single buffer inlet	Reduced complexity
Available with two connection types: <ul style="list-style-type: none">▪ TC 25 - ½"▪ AseptiQuik®* S	Flexibility and ease of use. TC for process development AseptiQuik® for aseptic connectivity and cGMP commercial manufacturing
Pre-sterilized and flushed with WFI	Ready to use
Low protein-binding	Extended operating lifetime and high product yield

*AseptiQuik® is a registered trademark of the Colder Products Company.



Technical Data

Specifications

Materials of Construction

Membrane	Hydrosart® Stabilized Cellulose
Gaskets	Polypropylene PVDF
Spacer	Polypropylene
Sealing compound	Grey Silicone
Plates with connector	Polypropylene

Retention Rates Hydrosart®

Substance	Approx. Mol. Wt.	30 kDa [%]
Albumin	67,000	≥ 97.5

Hydrosart® Single Pass Diafiltration Cassettes

Molecular Weight Cut-off: 30 kDa
Effective Filtration Area: 0.2 m²
Connectors: TC | AseptiQuik®

Operating Parameters

Maximum Feed pressure, P_{in} 43.5 psi | 3 bar

Operating temperature 40 °C maximum
pH stability 2 – 14 (Hydrosart®)

Air Diffusion Rate

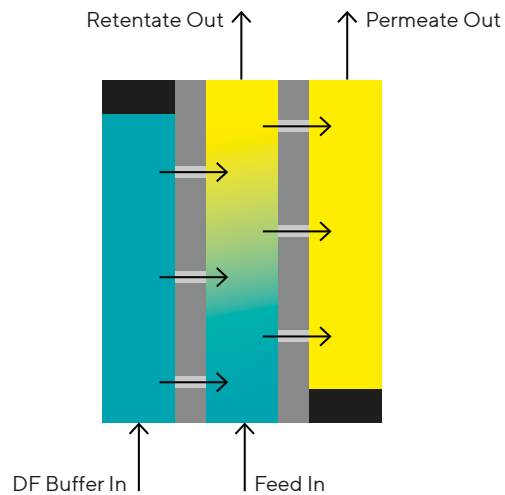
30 mL/min at P_{in} = 15 psi | 1 bar

Preconditioning

Pre-flushed to a low level of contaminants and gamma irradiated with 25 kGy.

Regulatory Compliance

All materials have passed the USP Biological Test.
All materials are BSE and TSE free. The filtrate meets or exceeds the current USP and EP for sterile Water for Injection, with respect to extractable substances, oxidizable substances and Bacteria Endotoxins.



Quality Control

Each Sartocoon® SP is individually assigned a serial number, integrity tested and certified.

Validation information is available upon request.
For further assistance, please contact your local Sartorius StedimBiotech Technical Service or our Goettingen-based Applications Department in Germany.

Technical References

Validation Guide
Publication no.: DIR: 3093174-000-01

Instructions for Use
Publication no.: 1000075275

Patent Rights

Sartocoon® SP and Sartoflow® SP are sold under a sublicense from Cytiva US LLC under U.S. Patent No. 7,384,549 B2 and/or its US and foreign equivalents. The aforementioned patent rights are owned by SPF Innovations, LLC.

By purchasing this product, customer is granted the right to use said product. However, the right to use does not include the right to make, to have made, to import, to sell, to resell or to have sold said product covered by one or more claims of the aforementioned patent rights.

Ordering Information

Available Sartocan® SP Types and Order Numbers

Cut-Off Pore Size	Connectors	Hydrosart® 0.2 m² Filter Area
30 kDa	TC 25 – ½"	3T514459M1465902
30 kDa	AseptiQuik® S	3Q514459M1465902

Related Products

Sartocan® SP cassettes are designed to be used with the Sartoflow® SP TFF system, Sartoflow® SP flow path, and Sartocan® Slice SC hydraulic filter holder.

Related Product	Order Code
Sartoflow® SP TFF system (including flow path)	SFSP-M-SYS
Sartoflow® SP flow path	SFSP-M-TUSET
Sartocan® Slice SC hydraulic holder	3D540SQ01-01

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
[sartorius.com](https://www.sartorius.com)