



Hydrosart® Ultrafiltration Cassettes

Protein purification,
concentration and diafiltration

Product Information

Hydrosart® is characterized by ultra-low protein adsorption, easy cleaning, higher yields, and long product life. Even with repeated use, Hydrosart® ultrafiltration membrane maintains its performance without fouling or loss of retention.



Description

Hydrosart® Membrane is a stabilized cellulose based membrane that has been optimized for the biopharmaceutical industry. Hydrosart® is stable across a broad pH range, is extremely hydrophilic, making it non-protein binding and virtually non-fouling. Membrane regeneration, storage and depyrogenation can be accomplished by using NaOH even at elevated temperatures. These features make Hydrosart® an ideal membrane for biological applications. Hydrosart® ultrafiltration cassettes are available in the following nominal molecular weight cutoffs: 2 kD | 5 kD | 10 kD | 30 kD | 100 kD | 300 kD

Applications

Hydrosart® ultrafiltration membranes are designed for use in the biotechnological and pharmaceutical industries. They can be used for the following applications:

- Oligonucleotide
- Proteins
 - Albumin, even with 40% EtOH
 - Hemoglobin

- Coagulation factors
 - Factor VIII
 - Factor III
- Vaccines
 - Tetanus
 - Diphtheria
- Monoclonal antibodies

Feature	Benefits
Non-adsorptive	No loss of proteins, easy to clean, sustained flux
Non-protein binding	High product yield
Wide pH and temperature range	More choices in sanitizing agents
High flow rates	Economical filtration runs
Self sealing cassette	No gaskets needed
Silicone sealing compound	Low Extractables
Enlarged inlet and outlet holes	Lower system pressure drops

Better solvent resistance than Polyethersulfone and Cellulose Triacetate

Technical Data

Specifications

Materials of construction	
Membrane	Hydrosart® (stabilized cellulose based membrane)
Gaskets	PVDF
Spacer	Polypropylene
Sealing compound	Silicone white

Pore size | Retention rate

Hydrosart® ultrafiltration cassettes are available in a choice of the following nominal molecular weight cut offs:
2 kD | 5 kD | 10 kD | 30 kD | 100 kD | 300 kD

Available sizes

Sartorius Stedim Biotech Crossflow Cassettes are available in Standard Cassette size for pilot- | production scale and in Sartococon® Slice format for reduced volume handling.

Available filter holder

Sartorius Stedim Biotech Crossflow Cassettes are designed for Sartorius Stedim Biotech filter holders like Sartococon® Slice (0.1 m² Cassettes only), Sartococon® 2 Plus, and different SARTOFLOW® holder.

Filtration area

Filter area Sartococon® Cassette	0.6 m ²
Filter area Sartococon® Slice Cassette	0.1 m ²

Operating parameters	
Feed pressure, P _{in}	58 psi 4 bar maximum
Operating temperature	50°C maximum, at 20°C
Max. air diffusion rates at P _{in} = 15 psi 1 bar	15 mL air/min for 0.6 m ² filter area 5 mL air/min for 0.1 m ² filter area
Cleaning	NaOH, 1 M, 40°C
Disinfection	NaOH, 1 M, 40°C, 30 min
Storage	NaOH, 0.1 M

Sterilization

NaOH, 1 M, 40°C, 30 min

Regulatory compliance

All materials have passed the USP Biological Test. The filtrate meets or exceeds the currently valid USP and EP for sterile Water for Injection, with respect to particulate matter, extractable substances, oxidizable substances, pH dependent conductivity, Ammonia, Chloride, Sulfate, Calcium and Bacteria Endotoxins.

Quality control

Each filter cassette is individually assigned a serial number, integrity tested and certified.

It complies with cGMP requirements for non-fiber-releasing filters and is filed under the Drug Master File Number DMF 5967 by the Food and Drug Administration, Washington, DC. Validation information is available upon request.

If you use holding devices from other suppliers, please contact our Applications Department. A different torque might be needed due to specific variations in design.

For further assistance, please contact your local Sartorius Stedim Biotech field engineer or our Goettingen-based Applications Department in Germany.

Technical references

Validation Guide
Publication No.: SPC5705-e

Directions for Use
(Sartococon® Cassettes and Sartococon® Slice Cassettes)
Publication No.: 1000035949

Average Dynamic Water Flux*

Nominal molecular weight cutoff (kD)	2 kD	5 kD	10 kD	30 kD	100 kD	300 kD
Permeate Flow Hydrosart® (L/h/m ²)	10	18	50	136	370	860

* (Feed pressure, P_{feed} = 29 psi | 2.0 bar; Retentate pressure, P_{ret} = 7.3 psi | 0.5 bar; P_{rem} = open valve)

Retention rates Hydrosart®

Substance	Approx. Mol. Wt.	2 kD	5 kD	10 kD	30 kD	100 kD	300 kD
Vitamin B12	1,200	≥ 88%	-	-	-	-	-
Inulin	5,000	-	> 96%	-	-	-	-
Cytochrome C	12,400	-	-	> 97.5%	-	-	-
Albumin	67,000	-	-	-	> 97.5%	≤ 60%	-
γ Globulin	169,000	-	-	-	-	≥ 96%	-
Blue Dextran	500,000	-	-	-	-	-	< 90%

Ordering Information

Available types and order numbers

Cut Off	Sartocon® Cassettes 0.6 m ² Filter Area	Sartocon® Slice Cassettes 0.1 m ² Filter Area
2 kD	302 144 19 06 E--SW*	305 144 19 01 E--SW*
5 kD	302 144 29 06 E--SW*	305 144 29 01 E--SW*
10 kD	302 144 39 06 E--SW*	305 144 39 01 E--SW*
30 kD	302 144 59 06 E--SW*	305 144 59 01 E--SW*
100 kD	302 144 68 06 E--SW*	305 144 68 01 E--SW*
300 kD	302 144 79 06 E--SW	305 144 79 01 E--SW

* For grey silicone cassettes material number ends with "...--SG"

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49.551.308.0

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen
Phone +49.5665.407.0

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex
Phone +33.442.845600

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex
Phone +33.442.845600

Austria

Sartorius Stedim Austria GmbH
Modectcenterstrasse 22
1030 Vienna
Phone +43.1.7965763.18

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.80

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi
Phone +36.23.457.227

Italy

Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41

Netherlands

Sartorius Stedim Netherlands B.V.
Phone +31.30.60.25.080
filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.647.38.40

Russian Federation

LLC "Sartorius Stedim RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone +34.913.586.098

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen
Phone +41.52.354.36.36

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ
Phone +44.1372.737159

Ukraine

LLC "Sartorius Stedim RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Africa

Republic of South Africa

Sartorius South Africa (Pty) Ltd
Unit 4, Alphen Square South
853 16th Road
Midrand 1685, RSA
Phone +27.11.315.5444

Americas

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1.800.368.7178

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Beijing Branch Office
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Guangzhou Branch Office
Room 1105
Xing Guang Ying Jing Building
No. 119, Shui Yin Road
Yue Xiu District, Guangzhou 510075
Phone +86.20.3836.4193

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.4331.4300

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Rd
The Alpha #02-13/14
Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, BunDang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700



▶ www.sartorius-stedim.com