



## Sartocon<sup>®</sup> Slice 50

Small scale process development disposable ultrafilters for biopharmaceutical applications

### Product Information

Sartocon<sup>®</sup> Slice 50 is available with Hydrosart<sup>®</sup> and Polyethersulfone membrane. It can easily be interfaced with the SARTOFLOW<sup>®</sup> Smart system.



### Sartocon<sup>®</sup> Slice 50

The Sartocon<sup>®</sup> Slice 50 is a ready to use 50 cm<sup>2</sup> fully disposable crossflow module that is ideally suited for membrane screening and process development, available in a wide variety of MWCO's and polymers. The Sartocon<sup>®</sup> Slice 50 can easily be interfaced with the SARTOFLOW<sup>®</sup> Smart system as well as many other laboratory scale crossflow systems.

### Flow Channel Geometry

Sartocon<sup>®</sup> Slice 50 is available in two flow channel geometry versions. E-Screen filters are dominantly used for protein concentrations up to above 20% as well as viscosity conditions greater than 3 cp. ECO-Screen devices mainly used for water based protein solutions of below 20% protein concentration as well as lower than 3 cp viscosity.

### The Hydrosart<sup>®</sup> Membrane

Hydrosart<sup>®</sup> is a stabilized cellulose based membrane that has been optimized for the biotechnological and pharmaceutical industry. The Hydrosart<sup>®</sup> membrane is a stable polymer that can operate in a broad pH range. Hydrosart<sup>®</sup> is also extremely hydrophilic, making it non-protein binding, virtually non-fouling, and has extremely high flux. Membrane depyrogenation can be accomplished by using NaOH even at elevated temperatures. These features make Hydrosart<sup>®</sup> an ideal membrane for biological applications.

Feature	Benefits
Available in all membrane materials and cut offs	Flexibility in membrane screening
Identical flow dynamics with larger modules	True linear scale up
Self contained unit	Ready to use, no compression holder needed
Low adsorption	Minimal product loss
Low protein-binding	High flux and product yield
DataMatrix code	Fast access to order code and shelf life data.
Luerlock Female ports	Quick connection to existing equipment ready to use on SARTOFLOW <sup>®</sup> Smart system
Low hold up volume	High concentration factor

### The Polyethersulfone Membrane

The polyethersulfone membrane (PESU) is a membrane polymer that is well established in the biotechnological and pharmaceutical industries. The PESU membrane is a stable polymer that can operate in a broad pH and temperature range. Polyethersulfone membranes are designed for use in the biotechnological and pharmaceutical industries.

## Technical Data

### Specifications Materials of Construction

Membrane	Hydrosart®   Polyethersulfone
Sealing and Gaskets	Silicone, Epoxide, Polypropylene
Spacer	Polypropylene, PVDF
Encapsulation	Reinforced ABS
Connector	ABS

### Available Filterholder

No special filter holder is required to operate a Sartocan® Slice 50. For ease of handling in vertical orientation a matching holder (order number 3ZG---0108) can be used.

### Rejection Coefficient (Characterization of Membrane)

#### Retention rates Hydrosart®

Substance	Approx. Mol. Wt.	2 kD [%]	5 kD [%]	10 kD [%]	30 kD [%]
Vitamin B12	1,200	≥ 88	-	-	-
Inulin	5,000	-	≥ 96	-	-
Cytochrome C	12,400	-	-	≥ 97,5	-
Albumin	67,000	-	-	-	≥ 97,5
γ-Globulin	169,000	-	-	-	-

#### Retention rates Polyethersulfone (PESU)

Substance	Approx. Mol. Wt.	1 kD [%]	5 kD [%]	8 kD [%]	10 kD [%]	30 kD [%]	50 kD [%]	100 kD [%]	300 kD [%]
Vitamin B12	1,200	> 70	50 – 80	-	-	-	-	-	-
Cytochrome C	12,400	-	-	≥ 99	> 95	60 – 90	-	-	-
Albumin	67,000	-	-	-	-	-	> 95	< 80	-
γ-Globulin	169,000	-	-	-	-	-	> 99	≥ 98	< 70
Dextran	2,000,000	-	-	-	-	-	-	-	> 95

### Operating Parameters

Feed pressure  $P_{in}$  3.5 bar maximum

Operating temperature range 2 – 40°C,

Air diffusion rates ≤ 3.0 mL/min at 2.5 bar

pH stability Hydrosart 2 – 14 | PESU 1 – 14

Hold up volume 1 – 2 ml

Recommended cross flow rate E Screen 50 – 120 ml/min

Recommended cross flow rate ECO Screen 20 – 60 ml/min

Cleaning NaOH, 1 M, 40°C

Disinfection NaOH, 1 M, 40°C, 30 min

Storage NaOH, 0.1 M

### Sterilization

No thermal sterilization possible.

### Regulatory Compliance

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

### Quality Control

Each Sartoco<sup>®</sup> Slice 50 is individually assigned a serial number, integrity tested and certified. It complies with cGMP requirements 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. For further assistance, please contact your local Sartorius Stedim Biotech Application Department.

### Technical References

Directions for Use	SPC6068
Validation Guide	SPC6067

## Ordering Information

Sartocon Slice 50 available as "E"-Screen and "ECO"-Screen Filter

Membrane Polymer <sup>™</sup>	Cut-Off	Material Number "E"-Screen Filter	Material Number "ECO"-Screen Filter
Hydrosart <sup>®</sup>	2 kD	3D91441950ELLPU	3D91441950MLLPU
Hydrosart <sup>®</sup>	5 kD	3D91442950ELLPU	3D91442950MLLPU
Hydrosart <sup>®</sup>	10 kD	3D91443950ELLPU	3D91443950MLLPU
Hydrosart <sup>®</sup>	30 kD	3D91445950ELLPU	3D91445950MLLPU
PESU	1 kD	3D91460950ELLPU	3D91460950MLLPU
PESU	5 kD	3D91462950ELLPU	3D91462950MLLPU
PESU	8 kD	3D91463450ELLPU	3D91463450MLLPU
PESU	10 kD	3D91463950ELLPU	3D91463950MLLPU
PESU	30 kD	3D91465950ELLPU	3D91465950MLLPU
PESU	50 kD	3D91465050ELLPU	3D91465050MLLPU
PESU	100 kD	3D91466850ELLPU	3D91466850MLLPU
PESU	300 kD	3D91467950ELLPU	3D91467950MLLPU

### Available Crossflow System

The SARTOFLOW<sup>®</sup> Smart is a modular small scale benchtop crossflow system optimized for concentration and diafiltration applications used in many downstream processes, such as purification of vaccines, monoclonal antibodies and recombinant proteins. The system is suitable for flexible use in laboratory environments for process development and clinical trials as well as for cGMP environments.



SARTOFLOW<sup>®</sup> Smart including options

# Sales and Service Contacts

For further contacts, visit [www.sartorius-stedim.com](http://www.sartorius-stedim.com)

## Europe

### Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289

Sartorius Stedim Systems GmbH  
Robert-Bosch-Strasse 5 – 7  
34302 Guxhagen

Phone +49.5665.407.0  
Fax +49.5665.407.2200

### France

Sartorius Stedim FMT S.A.S.  
ZI des Paluds  
Avenue de Jouques – CS 91051  
13781 Aubagne Cedex

Phone +33.442.845600  
Fax +33.442.845619

Sartorius Stedim France SAS  
ZI des Paluds  
Avenue de Jouques – CS 71058  
13781 Aubagne Cedex

Phone +33.442.845600  
Fax +33.442.846545

### Austria

Sartorius Stedim Austria GmbH  
Modecenterstrasse 22  
1030 Vienna

Phone +43.1.7965763.18  
Fax +43.1.796576344

### Belgium

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

Phone +32.2.756.06.80  
Fax +32.2.756.06.81

### Hungary

Sartorius Stedim Hungária Kft.  
Kagyló u. 5  
2092 Budakeszi

Phone +36.23.457.227  
Fax +36.23.457.147

### Italy

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

Phone +39.055.63.40.41  
Fax +39.055.63.40.526

### Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080  
Fax +31.30.60.25.099

[filtratie.nederland@sartorius-stedim.com](mailto: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  
Fax +48.61.879.25.04

### Russian Federation

LLC "Sartorius Stedim RUS"  
Uralskaya str. 4, Lit. B  
199155 St. Petersburg

Phone +7.812.327.53.27  
Fax +7.812.327.53.23

### Spain

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

Phone +34.913.586.098  
Fax +34.913.589.623

### Switzerland

Sartorius Stedim Switzerland AG  
Ringstrasse 24 a  
8317 Tagelswangen

Phone +41.52.354.36.36  
Fax +41.52.354.36.46

### U.K.

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

Phone +44.1372.737159  
Fax +44.1372.726171

### Ukraine

LLS "Sartorius RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine

Phone +380.44.411.4918  
Fax +380.50.623.3162

## Americas

### USA

Sartorius Stedim North America Inc.  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Toll-Free +1.800.368.7178  
Fax +1.631.254.4253

### Argentina

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

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962

São Bernardo do Campo  
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900  
Fax +55.11.4362.8901

### 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  
Fax +52.55.5562.2942

[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

### Peru

Sartorius Peru S.A.C.  
Av. Emilio Cavenecia 264 San Isidro  
15073 Lima, Perú

Phone +51.1.441 0158  
Fax +51.1.422 6100

## Asia | Pacific

### Australia

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

Phone +61.3.8762.1800  
Fax +61.3.8762.1828

### China

Sartorius Stedim Biotech (Beijing) Co. Ltd.  
No. 33 Yu'an Road  
Airport Industrial Park Zone B  
Shunyi District, Beijing 101300

Phone +86.10.80426516  
Fax +86.10.80426580

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  
Fax +86.21.6878.2882

Sartorius Stedim Biotech (Beijing) Co. Ltd.  
Guangzhou Representative Office

Unit K, Building 23  
Huihua Commerce & Trade Building  
No. 80 Xianlie Middle Road  
Guangzhou 510070

Phone +86.20.37618687 | 37618651  
Fax +86.20.37619051

### India

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

Phone +91.80.4350.5250  
Fax +91.80.4350.5253

### 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  
Fax +81.3.4331.4301

### 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  
Fax +60.3.8996.0755

### Singapore

Sartorius Stedim Singapore Pte. Ltd.  
1 Science Park Road,  
The Capricorn, #05-08A,  
Singapore Science Park II  
Singapore 117528

Phone +65.6872.3966  
Fax +65.6778.2494

### 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  
Fax +82.31.622.5799



▶ [www.sartorius-stedim.com](http://www.sartorius-stedim.com)