



Sartocell Depth Filter Modules for Food & Beverages

Filter Sheets for Clarification
and Microbe-retentive filtration

Product Information

Sartocell filter Modules are used for clarification and microbe-retentive filtration in the food and beverage industry. The use of modules in closed filtration housings eliminates factors like dripping and contaminations from external sources. The sheets are manufactured from high grade, natural raw materials such as cellulose and diatomaceous earth and are available in various sizes.



Double layer

Typical applications often require a coarser filter grade immediately followed by a finer grade, both in separate housings. Sartocell depth filter modules combine the two most popular media combinations in a single depth filter module. This enables the combination of 2 process steps into a one. This reduces the work for exchanging the filters and minimizes the footprint.

Protection

The positively charged filters improve the clarifying effect by adsorption of negatively charged particles and colloids. This ensures that downstream filters are effectively protected and also lowers overall costs of consumables.

Housings

Sartocell filter housings are especially designed for the use of Sartocell modules in the food and beverages industries. The 12" and 16" diameter housings come with a self draining base plate design and a replaceable adapter for the use of O-ring and flat gasket adapters. A detailed datasheet for Sartocell filter housings is available (Publication No.: SR-2008-e140509).

Applications

The Sartocell F&B filter sheets are especially developed for Wine, Water, Beer, Fruit juices and Spirits applications.

Flexibility

Filter sheet systems can be optimally adapted to the flow rates required for each application by raising or lowering the number of sheets installed.

Quality

Sartocell F&B are manufactured according to a DIN EN ISO 9001:2000 certified quality management system. All Sartocell F&B products are according to the Quality standards for food production set by the German LFGB and US FDA.

Technical Data

Specifications

Materials	
Filter media	Cellulosic depth filter with inorganic filter aids
Gaskets	Silicone
Core	Stainless steel
Plastics	Polypropylene

Single layer	
Cells per module	16
Filtration area	12" : 1.9 m ² 16" : 3.7 m ²
Height	276 mm

Double layer	
Cells pro module	8
Filtration area	1.9 m ²
Height	185 mm

Operating Parameters	
Max. allowable differential pressure	2.0 Bar 36 psi
Max. allowable back pressure	0.03 Bar 0.4 psi

Sterilization

Steam sterilization
121°C at 1 bar for 60 min.

Hot water flush
80°C at 1 bar for 30 min.

Ordering Information

Grade	Nominal retention [µm]	Thickness [mm]	Ash content [%]	Water flow [L/m ² /min]	Diameter [mm]	Article number
12" Module with 16 Cells, 1 piece per package						
S7	0.2	3.9	50	40	297	291S7-1216--F
S5	0.3	3.8	50	52	297	291S5-1216--F
S3	0.5	3.8	49	84	297	291S3-1216--F
F7	0.8	3.8	50	155	297	291F7-1216--F
F4	1.5	3.8	50	240	297	291F4-1216--F
C8	2.5	3.8	50	425	297	291C8-1216--F
C5	3.5	3.8	46	1550	297	291C5-1216--F

16" Module with 16 Cells, 1 piece per package						
S7	0.2	3.9	50	40	402	291S7-1616--F
S5	0.3	3.8	50	52	402	291S5-1616--F
S3	0.5	3.8	49	84	402	291S3-1616--F
F7	0.8	3.8	50	155	402	291F7-1616--F
F4	1.5	3.8	50	240	402	291F4-1616--F
C8	2.5	3.8	50	425	402	291C8-1616--F
C5	3.5	3.8	46	1550	402	291C5-1616--F

Double layer grades	Grade combination	Diameter	Article number
CF5	C5 + F4	402	291CF51616--F
FS2	F4 + F7	402	291FS21616--F

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
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
Modcenterstrasse 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.p.A.
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

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 +34.902.110.935
Fax +34.91.358.96.23

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

LLC "Biohit"
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.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

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