

Sartocell Depth Filter Modules for Food and 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 two process steps into 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 beverage industry. 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 filter sheets are manufactured according to a DIN EN ISO 9001:2000 certified quality management system. All Sartocell F&B products are manufactured according to the quality standards for food production set by the German LFGB and the 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

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
www.sartorius.com