



Sartoclear® Depth Filter Sheets for Food & Beverages

Filter Sheets for Clarification and
Microbe-retentive filtration

Product Information

Sartoclear® filter sheets are used for clarification and microbe-retentive filtration in the food and beverage industry.

The sheets are manufactured from high grade, natural raw materials such as cellulose and diatomaceous earth and can be available in various sizes.



Total Throughput

The structure of the cellulose forms a three-dimensional matrix. The Sartoclear® F&B filter sheets have a large void volume of up to 85%. These features ensure outstanding total throughput and thereby cost-effective filtration.

Protection

The positive Zeta potential of the Sartoclear® F&B filter sheets strengthens the mechanical retention capacity and clarifying effect, allowing additional adsorption of minute negatively charged particles and colloids. This ensures that downstream filters are effectively protected and also lowers overall costs of consumables.

Applications

The Sartoclear® F&B filter sheets are especially developed for Wine, 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

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

Ordering Information

Grade	Nominal retention (µm)	Thickness (mm)	Ash content (%)	Water flow (L/m ² /min)	Size (mm)	Article number
S7	0.2	3.9	50	40	400 × 400	290S7-040040FN
S5	0.3	3.8	50	52	400 × 400	290S5-040040FN
S3	0.5	3.8	49	84	400 × 400	290S3-040040FN
F7	0.8	3.8	50	155	400 × 400	290F7-040040FN
F4	1.5	3.8	50	240	400 × 400	290F4-040040FN
C8	2.5	3.8	50	425	400 × 400	290C8-040040FN
C5	3.5	3.8	46	1550	400 × 400	290C5-040040FN
S7	0.2	3.9	50	40	600 × 615	290S7-060060FK
S5	0.3	3.8	50	52	600 × 615	290S5-060060FK
S3	0.5	3.8	49	84	600 × 615	290S3-060060FK
F7	0.8	3.8	50	155	600 × 615	290F7-060060FK
F4	1.5	3.8	50	240	600 × 615	290F4-060060FK
C8	2.5	3.8	50	425	600 × 615	290C8-060060FK
C5	3.5	3.8	46	1550	600 × 615	290C5-060060FK

Special formats are available on request.

Sartorius Stedim Biotech GmbH
 August-Spindler-Strasse 11
 37079 Goettingen, Germany
 Phone +49.551.308.0
 Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
 Argentina +54.11.4721.0505
 Brazil +55.11.4451.6226
 Mexico +52.5555.62.1102
 UK +44.1372.737159
 France +33.442.845600
 Italy +39.055.63.40.41
 Spain +34.913.586.098
 Russian Federation +7.812.327.5.327
 Japan +81.3.4331.4300
 China +86.21.68782300

Specifications subject to change without notice.
 Copyright Sartorius Stedim Biotech GmbH. Printed
 in the EU on paper bleached without chlorine.
 Publication No.: SR-2010-e160605
 Order No.: 85030-532-02
 Ver. 06 | 2016