

# Sartoclear® Depth Filter Sheets for Food and 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 are 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 filter sheets are manufactured according to a DIN EN ISO 9001 certified quality management system. All materials of construction used to produce Sartoclear® F&B products are selected in compliance with the EU and US requirements for products intended to come in contact with foodstuff:

- Meet the requirements for European Food Contact Materials Regulation as detailed in Regulation (EC) No 1935/2004
- Meet the United States Food and Drug Administration (US FDA) requirements as detailed in Code of Federal Regulations (CFR), Title 21, paragraphs 170–189. The FDA Compliance for indirect food additives is based on the information provided by our raw material suppliers.

## Order Information

Grade	Nominal Retention [µm]	Thickness [mm]	Ash Content [%]	Water Flow [L/m <sup>2</sup> /min]	Size [mm]	Article Number	Quantity [pieces]
S7	0.2	3.9	50	40	400 × 400	290S7-040040FN	100
S5	0.3	3.8	50	52	400 × 400	290S5-040040FN	100
S3	0.5	3.8	49	84	400 × 400	290S3-040040FN	100
F7	0.8	3.8	50	155	400 × 400	290F7-040040FN	100
F4	1.5	3.8	50	240	400 × 400	290F4-040040FN	100
C8	2.5	3.8	50	425	400 × 400	290C8-040040FN	100
C5	3.5	3.8	46	1,550	400 × 400	290C5-040040FN	100
C3	4.0	3.8	42	1,900	400 × 400	290C3-040040FN	100

S7	0.2	3.9	50	40	600 × 615	290S7-060060FK	50
S5	0.3	3.8	50	52	600 × 615	290S5-060060FK	50
S3	0.5	3.8	49	84	600 × 615	290S3-060060FK	50
F7	0.8	3.8	50	155	600 × 615	290F7-060060FK	50
F4	1.5	3.8	50	240	600 × 615	290F4-060060FK	50
C8	2.5	3.8	50	425	600 × 615	290C8-060060FK	50
C5	3.5	3.8	46	1,550	600 × 615	290C5-060060FK	50
C3	4.0	3.8	42	1,900	600 × 615	290C3-060060FK	50

Special formats are available on request.

**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  
[sartorius.com](https://www.sartorius.com)