# Product Datasheet

# SARTURIUS

# 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 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.

## Order Information

Grade	Nominal Retention [μm]	Thickness [mm]	Ash Content [%]	Water Flow [L/m²/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.

#### 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

