



Jumbo Star Sartopure®

High Capacity Particle & Bioburden Reduction
Filter Cartridges for Beverage Applications

Benefits

- Extend the life time of the final membrane filter before filling if used as prefilter before bottling
- Reduce filter consumption
- Highest filtration area on smallest footprint: reduce hardware investment costs
- Robustness & configuration adapted to backwash
- Quick installation



Product Information

With the new Jumbo Star Sartopure® 0.45 µm Sartorius introduces the finest available fleece material on the market to the food and beverage industry. The new fleece material shows an unmatched total throughput, particle retention, protective abilities of the final membrane filtration before filling at an extreme flow rate. The all polypropylene construction makes it resistant to all common cleaning chemicals and numerous sterilization cycles.

Jumbo Star Sartopure® cartridges combine the highest filtration area on the smallest footprint on the market. Particle-rich beverages can be filtered with Jumbo Star cartridges that are easily and efficiently backwashed with water. This allows for a significant reduction of filtration area, filter consumption and overall costs for filtration. Sartopure® Jumbo Star filter cartridges are available in the widest range of retention rates and their modular design allows the adaptation to all production requirements.

Applications

- Prefiltration upstream of the final membrane filtration before filling (water, wine and beer applications)
- Trapfiltration downstream of the diatomaceous earth | PVPP filter for beer industry
- Clarification step in winery and distillery
- Retention of tartar crystals after cold treatment of wine
- Retention of particles, reduction of colloids and bioburden in water, wine and beer and in alcohol production, also for venting large tanks
- Possibility of use on raw wine as an alternative to Kieselghur filter or Cross-flow filter or as cellar filter after tartaric stabilization

Documentation

Jumbo Star Sartopure® filters are designed and manufactured according to an ISO 9001 certified quality management system. All raw materials which are used to produce Jumbo Star Sartopure® are in accordance with EU regulation 1935/2004/EC or are listed in the Regulations (EU) No. 10/2011 relating to plastic materials and article intended to come into contact with food. All used materials comply with FDA requirements.

Flexibility

The modular concept allows the ideal adaptation to production requirements from 7 m² to 28 m² of filtration area.

Technical Data

Specifications

Materials	
Filter Material	Multiple Polypropylene layers
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone

Retention Rates
0.45 µm, 0.65 µm, 1.2 µm, 3 µm, 5 µm, 8 µm, 20 µm

Available Sizes | Filtration Area

Size 1	10"	7 m ²
Size 2	20"	14 m ²
Size 3	30"	21 m ²
Size 4	40"	28 m ²

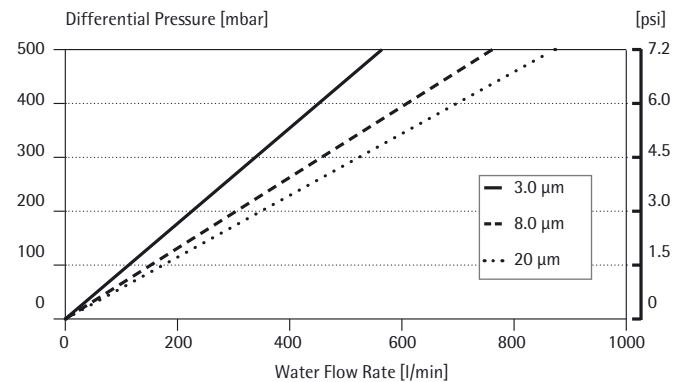
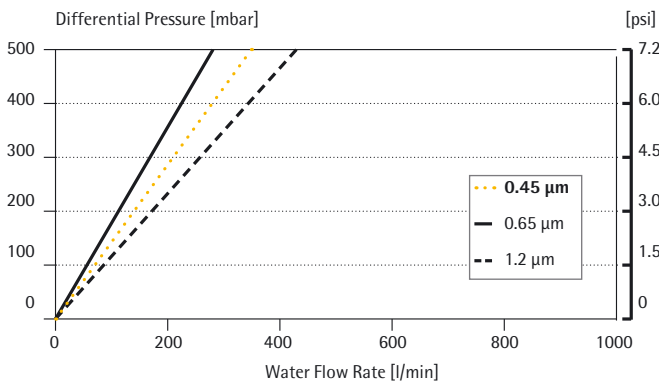
Available Adapter

40

Operating Parameters	
Max. Allowable Differential Pressure	4 bar 58 psi at 20°C 1 bar 15 psi at 80°C 0.5 bar 7.5 psi at 120°C
Max. Allowable Back Pressure	1.5 bar 22 psi at 20°C

Sterilisation	
Hot water at	85–95°C, 30 minutes, ΔP max 0.5 bar
Steam at	110°C, 30 minutes, ΔP max 0.5 bar
Sterilization cycles	min. 100

Water Flow Rates for 10" Cartridges



Ordering Information

Order Code	Retention Rate [μm]
5504006JX	0.45 μm
5594005JX	0.65 μm
5594003JX	1.2 μm
5594002JX	3.0 μm
5594042JX	5.0 μm
5594001JX	8.0 μm
5594020JX	20.0 μm

X = 1 Size 10"
X = 2 Size 20"
X = 3 Size 30"
X = 4 Size 40"

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

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen
Phone +49.5665.407.0

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex
Phone +33.442.845600

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex
Phone +33.442.845600

Austria

Sartorius Stedim Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965763.18

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.80

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi
Phone +36.23.457.227

Italy

Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41

Netherlands

Sartorius Stedim Netherlands B.V.
Phone +31.30.60.25.080
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

Russian Federation

LLC "Sartorius Stedim RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone +34.913.586.098

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen
Phone +41.52.354.36.36

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ
Phone +44.1372.737159

Ukraine

LLC "Sartorius Stedim RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Toll-Free +1.800.368.7178

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

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

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Beijing Branch Office
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Guangzhou Branch Office
Room 1105
Xing Guang Ying Jing Building
No. 119, Shui Yin Road
Yue Xiu District, Guangzhou 510075
Phone +86.20.3836.4193

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

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

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

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

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



▶ www.sartorius-stedim.com