

Jumbo Star Sartopure®

High Capacity Particle &
Bioburden Reduction
Filter Cartridges for Beverage
Applications

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.



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

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 Kieselgur 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
------------------------------	--------------------------

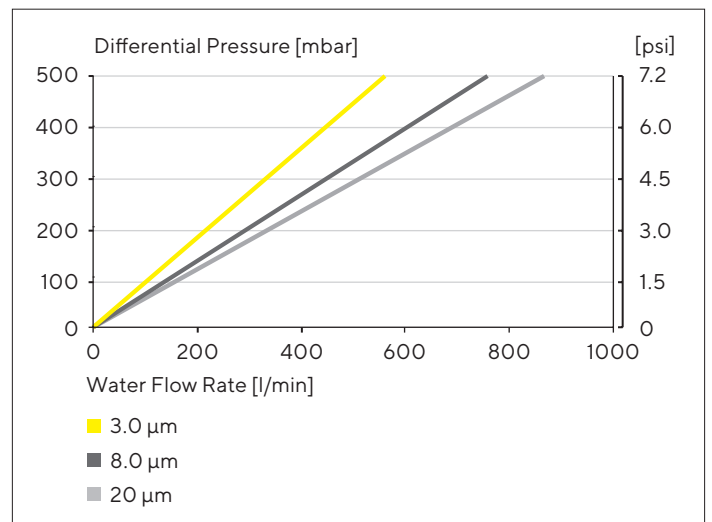
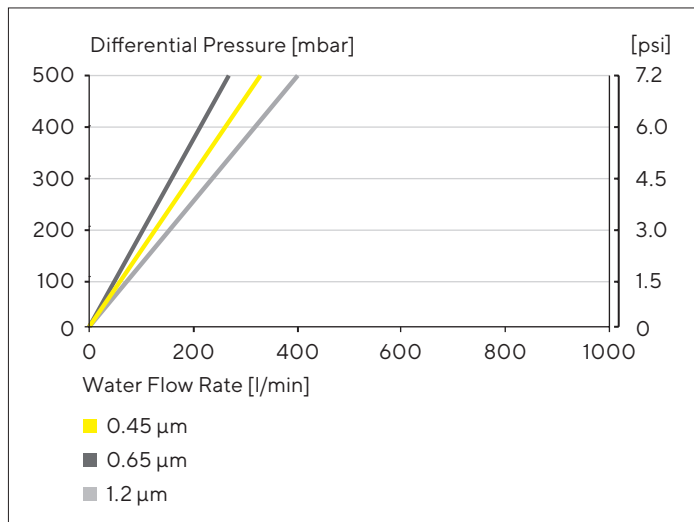
Sterilisation

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"

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