

Jumbo Star Sartopure® PP2

A "Giant", Ready-to-Use
Particulate Filtration &
Bioburden Reduction Filter



Description

Jumbo Star Sartopure® PP2 modular filter elements can be used for a wide range of prefiltration applications. Retention of hard, non deformable particles and reduction of bioburden from liquids is achieved through fractionated defined depth filtration. Jumbo Star Sartopure® PP2 combine multiple layers of progressively finer polypropylene depth filter fleeces in a pleated format.

Application

Jumbo Star Sartopure® PP2 are the ideal choice for prefiltration and clarification of:

- Plasma Fractions
- LVP Solution
- Ophthalmics
- WFI
- Process water

Security

Jumbo Star Sartopure® PP2 filter elements ensure selective and defined particle retention. They are a valuable protection for the final membrane filter. The completely polypropylene construction offers a broad chemical compatibility.

Economic Prefiltration

Highest dirt loading capacities in combination with high flow rates make Jumbo Star Sartopure® PP2 filter elements an ideal choice for a variety of large-scale filtration applications in the Pharmaceutical industry. Filtration costs are reduced to a minimum as these filters provide a long service life in many applications.

Reliable Operation

The new modular construction together with effective combination of nonwoven polypropylene achieves highest process reliability and reproducible results from batch to batch even under varying process conditions.

Cost Saving

The modular filter construction in combination with the expanded filter area results in a smaller filter housing, minimizing the required footprint. Jumbo Star filters are available in 4 different cartridge sizes to match a wide range of batch sizes.

Flexibility

Jumbo Star Sartopure® PP2 filter elements are available in a modular design from 7 m² filter area up to 28 m² filter area. This flexibility facilitates an easy adoption to your filtration process, depending on the batch size.

Documentation

Jumbo Star Sartopure® PP2 filter are designed, developed and manufactured in accordance with a ISO 9001 certified Quality Management System. A Validation Guide is available for compliance with regulatory requirements.

Specifications

Materials

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

Retention Rates

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 14.5 psi at 80 °C 0.5 bar 7.2 psi at 120 °C
Max. Allowable Back Pressure	1.5 bar 22 psi at 20 °C

Extractables

Jumbo Star Sartopure® PP2 filter meet, or exceed the requirements for WFI quality standards set by current USP without the need for a WFI flush, prior to use.

Regulatory Compliance

Non pyrogenic according to USP Bacterial Endotoxins Testing

Pass USP Plastic Class VI Test

Sterilization

Inline-Steam Sterilization

121 °C, 30 min. at max differential pressure of 0.5 bar | 7 psi

Autoclaving

134 °C, 30 min, 2 bar | 29 psi

Technical References:

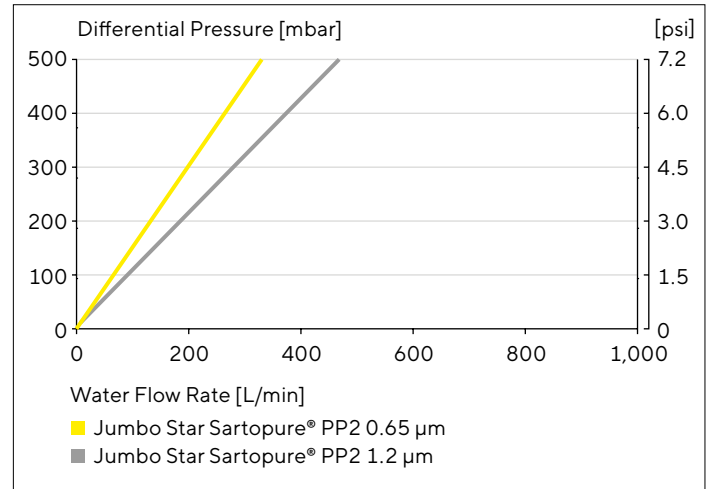
Validation Guide: DIR No.: 2930063

Ordering Information

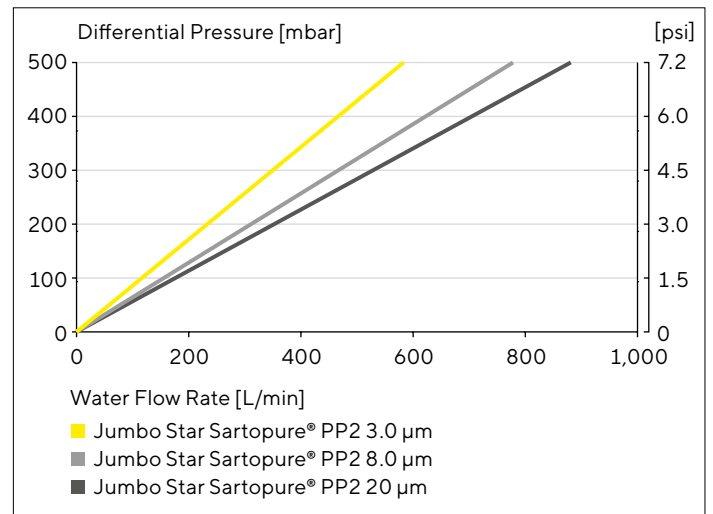
Order Code	Retention Rate
5594005JX	0.65 µm
5594003JX	1.20 µm
5594002JX	3.00 µm
5594042JX	5.00 µm
5594001JX	8.00 µm
5594020JX	20.0 µm

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

Water Flow 10" Cartridge



Water Flow 10" Cartridge




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)