SVISCISVS

Product Datasheet

Jumbo Star Sartopure GF Plus

A "Giant" Step Forward in Pleated Depth Filters



Description

Jumbo Star Sartopure GF Plus modular filter elements are ideal for removal of contaminates like colloids, lipids, protein aggregates (Host Cell Protein) and bioburden from biopharmaceutical fluids. They can be used for large-scale cell harvest clarification in lieu of lenticular filters and for aggregates removal in large-scale Protein Pool filtrations. They offer excellent protection to the membrane filters & chromatography columns in downstream processing. These filters do not contain any DE embedded in a loose cellulosic matrix. As a result, these GF Plus filters typically have significantly less Extractables compared to the Lenticular filters.

Application

Jumbo Star Sartopure GF Plus are the ideal choice for large-scale prefiltration and clarification of:

- Harvested Cell Culture fluids
- Microbial Fermentation broths
- Serum free or serum containing cell culture media
- Process Intermediates containing lipids, colloids and protein aggregates as contaminants.

Effective Clarification

Jumbo Star Sartopure GF Plus feature glass fiber layers for an effective clarification of fluid streams based on the combination of adsorption and sieve retention.

Economic Prefiltration

Based on a cutting-edge pleating technology; extremely high filter area is incorporated in each 10" filter element. In addition, the 3-dimensional filter matrix of Sartopure GF Plus depth filters ensures outstanding total throughput performance thus ensuring most economic design of your prefiltation scheme.

Reliable Operation

The new modular construction, coupled with effective combinations of nonwoven polypropylene and glass fibre layers, achieves the 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 GF Plus filter elements are available in a modular design from 5 m² up to 20 m² of filter area. This flexibility facilitiates an easy adoption to your filtration process, depending on the batch size.

Documentation

Jumbo Star Sartopure GF Plus 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 glass fibre layers
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone

Retention Rates

0.65 µm 1.2 µm

Available Sizes | Filtration Area

Size 1	10″	5 m²
Size 2	20″	10 m ²
Size 3	30″	15 m²
Size 4	40″	20 m ²

Available Adapter

40

Operating Parameters

Max. Allowable	4 bar 58 psi at 20°C
Differential Pressure	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 GF Plus meet or exceed the requirements for WFI quality standards set by current USP after WFI flush.

Regulatory Compliance

Non pyrogenic according to USP Bacterial Endotoxins Testing

Passed USP Plastic Class VI Test

Non-fiber releasing according to 21 CFR

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:

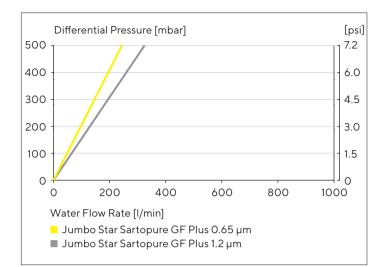
Validation Guide: SPK5774-e09021 85034-537-52

Ordering Information

Order Code	Retention Rate
5554005JX	0.65 µm
5554003JX	1.20 µm

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

Water Flow 10" Cartridge



Germany

USA

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen Phone +49 551 308 0 Sartorius Stedim North America Inc. 565 Johnson Avenue Bohemia, NY 11716 Toll-Free +1 800 368 7178

For further contacts, visit www.sartorius.com

Specifications subject to change without notice. Copyright Sartorius Stedim Biotech GmbH. DIR No.: 2648617-000-00 Order No.: 85034-537-62 Status: 11 | 2020