

Sartopure[®] GF Plus

The New Generation of Adsorptive Depth Filters



Description

Sartopure[®] GF Plus adsorptive depth filters are designed for removal of contaminants like colloids, lipids, protein aggregates (Host Cell Protein) and particles from biopharmaceutical fluids. They are used for protection of membrane filters, chromatography columns and ultrafiltration systems in pharmaceutical and biotechnological production processes.

Applications

Sartopure[®] GF Plus adsorptive depth filters are the ideal choice for prefiltration and clarification of:

- Cell Culture fluids after cell harvest
- Fermentation broths
- Serum free or serum containing cell culture media
- Serum
- Highly viscous ophthalmic and LVP solutions
- All media containing lipids and colloids as contaminants

Effective Clarification

Sartopure® GF Plus adsorptive depth filters feature highly charged glass fiber layers for effective clarification of fluid streams based on the combination of adsorptive and mechanical retention.

Economic Prefiltration

The 3-dimensional filter matrix of Sartopure® GF Plus adsorptive depth filters assures outstanding total throughput performance of the filters thus ensuring most economic design of your prefiltration scheme.

Reliable Operation

The high and defined particle retention capability of Sartopure® GF Plus allows reliable operation and reproducible results from batch to batch even under varying process conditions.

Cost Saving

The efficient protection of downstream membrane filters and purification equipment saves filter costs and helps to increase the yield of biotech production processes.

Flexibility

Sartopure® GF Plus filters are available as standard cartridges and Maxicaps®. Cartridges are strong and robust and designed for maximum pressure differentials and multiple steaming cycles. Disposable Maxicaps are designed for single use and are integral component of disposable manufacturing lines.

Documentation

Sartopure® GF Plus adsorptive depth filters are designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System. A Validation Guide is available for compliance with regulatory requirements.

Technical Data

Materials	
Filter Material:	Glass Fiber Fleeces
Support Fleece:	Polypropylene
Core:	Polypropylene
End Caps:	Polypropylene
O-Rings:	Silicone (optional EPDM or Fluoroelastomer)

Retention Rates

0.65 µm, 1.2 µm

Available Sizes|Filtration Area (nominal)

Size 1	10"	0.4 m ² 4.3 ft ²
Size 2	20"	0.8 m ² 8.6 ft ²
Size 3	30"	1.2 m ² 12.9 ft ²
Size 4	40"	1.6 m ² 17.2 ft ²

Available Adapters Cartridges

21, 25, 27, 28

Operating Parameters

Max. Allowable Differential Pressure	5 bar 72.5 psi at 20°C 2 bar 29 psi at 80°C
Max. Allowable Back Pressure	2 bar 29 psi at 20°C

Extractables

Sartopure® GF Plus cartridges meet, or exceed the requirements for WFI quality standards set by the current USP.

Regulatory Compliance

Non pyrogenic according to USP Bacterial Endotoxins

Pass USP Plastic Class VI Test

Non fiber releasing according to 21 CFR

Sterilization

In-line Steam Sterilization:
134°C, 20 min. at max differential pressure of 0.5 bar

Note

Maxicaps® cannot be in-line steam sterilized!

Autoclaving:
134°C, 2 bar, 30 min

Sterilization Cycles

In-Line sterilization Min. 25
Autoclaving: Min. 25

Technical References

Validation Guide SPK5762-e

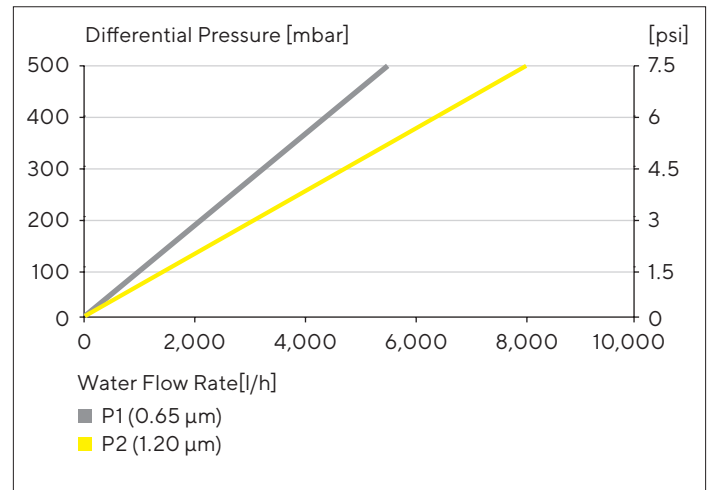
Order Information

Order Code	Retention Rate [µm]	Pack Size [Pieces]
Cartridges		
555**05P1	0.65	1
555**05P2	0.65	1
555**05P3	0.65	1
555**05P4	1.2	1
555**03P1	1.2	1
555**03P2	1.2	1
555**03P3	1.2	1
555**03P4	1.2	1

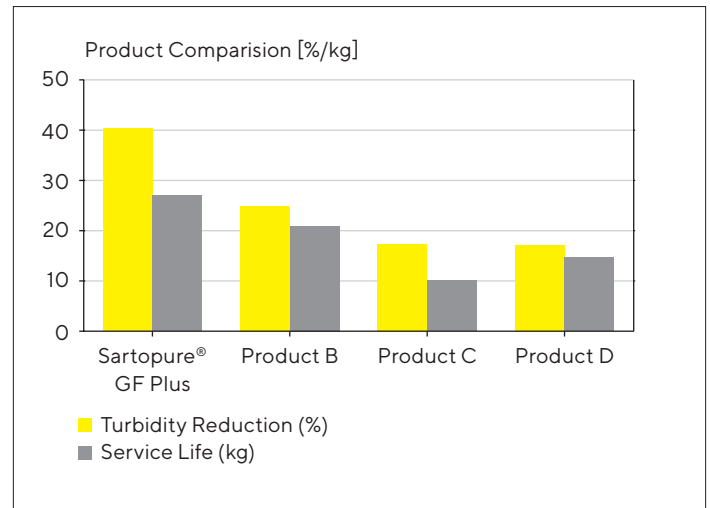
** Adapter Style

Water Flow Rates for Sartopure® GF Plus

Sartopure® GF Plus 10" Standard Cartridges 0.65 µm, 1.2 µm



Product Comparison Data




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