

Sartopure[®] GF Plus

Adsorptive Depth Filters



Product Information

Sartopure[®] GF Plus Cartridges, Maxicaps[®] and Midicaps[®] are depth filter elements utilizing two layers of pleated glass fleece. The highly adsorptive non fiber releasing material is ideally suited for the removal of contaminants like colloids, lipids, protein aggregates and particles. The unique construction results in an unmatched combination of total throughput and turbidity reduction.

Applications

Sartopure[®] GF Plus filter elements are the ideal choice for prefiltration and clarification of:

- Cell Cultures
- Fermentation broths
- Serum free or serum containing cell culture media
- All media containing lipids, colloids and protein aggregates.

Effective Clarification

Sartopure® GF Plus Cartridges, Maxicaps® and Midicaps® feature highly adsorptive glass fiber layers for effective clarification of fluid streams based on the combination of adsorptive and mechanical retention.

Robust Processes

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.

Economic Filtration

The efficient protection of downstream equipment saves filtration costs and increases the total throughput of biotech production processes.

Technical Data

Specifications

Materials	
Filter Matrial	Glass fiber fleeces double layer
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
Capsule Housing	Polypropylene
O-Ring	Silicone (optional EPDM or Fluoroelastomer)

Retention Rates	
0.65 µm	
1.2 µm	

Flexibility

Sartopure® GF Plus Cartridges, Maxicaps® and Midicaps® are available with various filtration areas from 400 cm² | 0.4 ft² up to 1.6 m² | 17.2 ft² for easy adoption to any filtration process independent from the batch size.

Documentation

Sartopure® GF Plus filter elements 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.

Available Sizes | Filtration Area (nominal)

Cartridges Maxicaps®	
10"	0.4 m ² 4.3 ft ²
20"	0.8 m ² 8.6 ft ²
30"	1.2 m ² 12.9 ft ²
40"	1.6 m ² 17.2 ft ² (only Cartridges)

Midicaps®	
Size 7	0.04 m ² 0.4 ft ²
Size 8	0.08 m ² 0.9 ft ²
Size 9	0.12 m ² 1.3 ft ²
Size 0	0.25 m ² 2.7 ft ²

Available Adapters Cartridges

- 21: Double open end with flat gasket
- 25: 2 Flange bayonet adapter with 226 double o-ring
- 27: Bayonet adapter with 222 double o-ring
- 28: 3 Flange bayonet adapter with 222 double o-ring

Available Connectors Midicaps®

- S: 1½" Tri-Clamp (Sanitary)
- O: ½" single stepped hose barb
- F: ¾" Tri-Clamp (Sanitary)
- H: ¼" Multiple stepped hose barb (only size 7)

Available Connectors Maxicaps®

- S: 1½" Tri-Clamp (Sanitary)
- O: ½" single stepped hose barb

Operating Parameters

Max. Allowable Differential Pressure	5 bar 72.5 psi at 20°C (Cartridges, Maxicaps®, Midicaps®) 3 bar 43.5 psi at 50°C (Maxicaps®, Midicaps®) 2 bar 29 psi at 80°C (Cartridges)
--------------------------------------	--

Max. Allowable Back Pressure	2 bar 29 psi at 20°C
------------------------------	------------------------

Extractables

Sartopure® GF Plus Cartridges, Maxicaps® & Midicaps® 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

for Cartridges

In-Line Steam Sterilization

134 °C, 20 min. at max differential pressure of 0.5 bar

or

Autoclaving

134°C, 2 bar, 30 min

for Maxicaps® & Midicaps®

Autoclaving

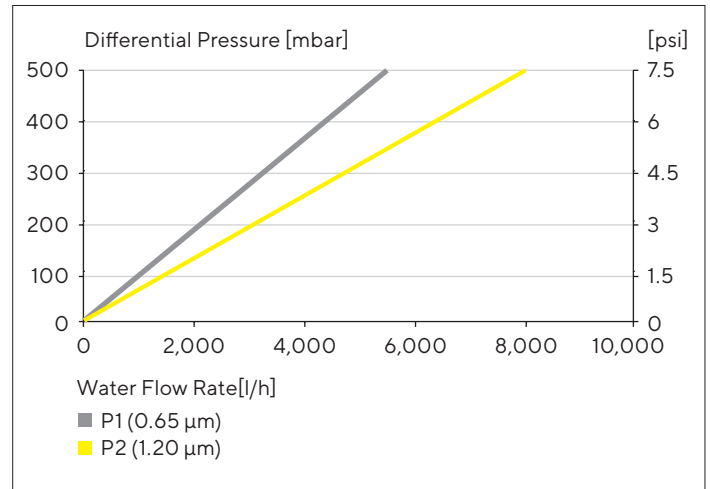
134°C, 2 bar, 30 min

Technical References

Validation Guide

SPK5762-e

Water Flow Rates for Sartopure® GF Plus 10" Standard Cartridges



Ordering Information

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

Maxicaps®

5557305P1-**	0.65	1
5557305P2-**	0.65	1
5557305P3-**	0.65	1
5557303P1-**	1.2	1
5557303P2-**	1.2	1
5557303P3-**	1.2	1

** Connector|Adapter Styles


Order Code	Retentation Rate [μm]	Pack Size [Pieces]
Midicaps®		
5555305P7-**-A	0.65	4
5555305P8-**-A	0.65	4
5555305P9-**-A	0.65	4
5555305P0-**-V	0.65	2
5555303P7-**-A	1.2	4
5555303P8-**-A	1.2	4
5555303P9-**-A	1.2	4
5555303P0-**-V	1.2	2

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