



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.

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.

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	

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®

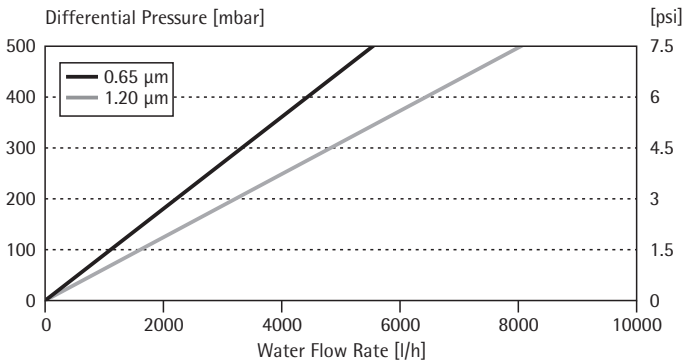
Autoclaving

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	Retention 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

Order Code	Retention Rate [µm]	Pack Size [Pieces]
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

Order Code	Retention 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

** Connector | Adapter Styles

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5 – 7
34302 Guxhagen

Phone +49.5665.407.0
Fax +49.5665.407.2200

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques – CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Modcenterstrasse 22
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi

Phone +36.23.457.227
Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.p.A.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080
Fax +31.30.60.25.099

filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.647.38.40
Fax +48.61.879.25.04

Russian Federation

LLC "Sartorius ICR" and LLC "Biohit"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone +34.902.110.935
Fax +34.91.358.96.23

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen

Phone +41.52.354.36.36
Fax +41.52.354.36.46

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax + 55.11.4362.8901

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Stedim Biotech (Beijing) Co. Ltd.
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim (Shanghai)
Trading Co. Ltd
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, China

Phone +86.21.68782300
Fax +86.21.68782332 | 68782882

Sartorius Stedim Biotech (Beijing) Co. Ltd.
Guangzhou Representative Office
Unit K, Building 23
Huihua Commerce & Trade Building
No. 80 Xianlie Middle Road
Guangzhou 510070

Phone +86.20.37618687 | 37618651
Fax +86.20.37619051

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300
Fax +81.3.4331.4301

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, BunDang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799



▶ www.sartorius-stedim.com