



# 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.

## Specifications

### 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 <sup>2</sup>	4.3 ft <sup>2</sup>
Size 2	20"	0.8 m <sup>2</sup>	8.6 ft <sup>2</sup>
Size 3	30"	1.2 m <sup>2</sup>	12.9 ft <sup>2</sup>
Size 4	40"	1.6 m <sup>2</sup>	17.2 ft <sup>2</sup>

### 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

## Specifications

### 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

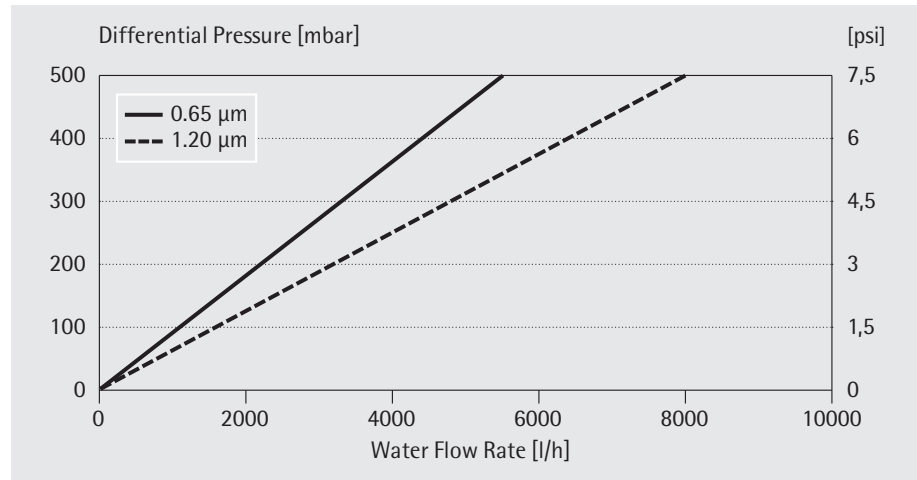
### Order Information

Order Code	Retentation Rate [µm]	Pack Size [pieces]
<b>Cartridges</b>		
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

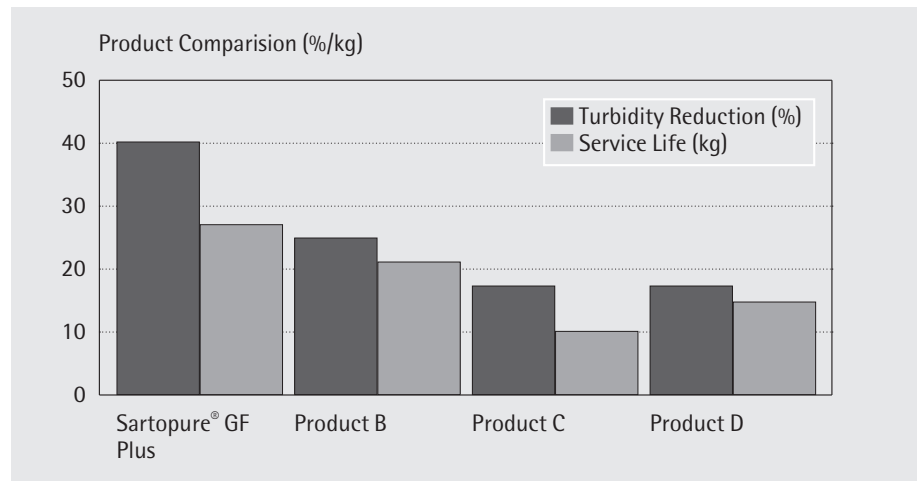
\*\* : Adapter Style

## Water Flow Rates for Sartopure® GF Plus

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



## Product Comparison Data



Sartorius Stedim Biotech GmbH

August-Spindler-Strasse 11  
37079 Goettingen, Germany

Phone +49.551.308.0

Fax +49.551.308.3289

www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178

UK +44.1372.737159

France +33.442.845600

Italy +39.055.63.40.41

Spain +34.90.2110935

Russian Federation +7.812.327.5.327

Japan +81.3.4331.4300

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH. | W  
Publication No.: SPK2064-e140508  
Order No.: 85030-521-97

Ver. 05 | 2014