

# Sartopure® GF Plus

Adsorptive High Capacity  
Filter Cartridges  
(Food & Beverage)



## Applications

Sartopure® GF Plus are ideal for the prefiltration of:

- Water
- Wine
- other beverages

## Description

Sartopure® GF Plus cartridges are optimized for the protection of downstream membrane filter systems and clarification in food and beverage applications. Due to their high adsorptive power by glass fiber fleeces, they are the ideal solution for removal of colloids from beverages.

## High Adsorption

Based on electrostatic forces of attraction, the negatively charged glass fiber material retains positively charged, small particles that would otherwise pass through the filter. These particles can even be smaller than the nominal filter retention rate.

## Defined Retention Rates

The key feature of Sartopure® GF Plus cartridges is their defined and accurate retention rating >99.9%!

## High Profitability

The cartridges' high flow rates at a low differential pressure reduce overall costs as less energy is used. Moreover, a high total throughput guarantees the maximum in process profitability.

## Flexibility

Sartopure® GF Plus filters are available as standard filter cartridges.

## Documentation

Sartopure® GF Plus cartridges are designed, developed and manufactured in accordance with a ISO 9001 certified Quality Management System.

## 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

### Cartridges

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

## Sterilization

### In-Line Steam Sterilization

134°C, 20 min. at max differential pressure of 0.5 bar|7 psi

### Sterilization Cycles

In-Line Sterilization: Min. 25

## Technical References

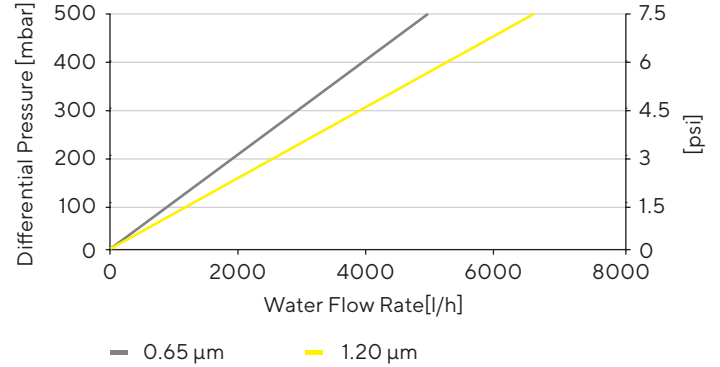
Validation Guide SPK5762-e

# Order Information

Order Code	Retentation Rate [ $\mu\text{m}$ ]	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 10" Standard Cartridges 0.65 $\mu\text{m}$ , 1.2 $\mu\text{m}$




### 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](http://www.sartorius.com)

Specifications subject to change without notice.  
© 2021 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

DIR No.: 2672589-000-00 | Order No.: 85030-519-91 | Status: 01 | 13 | 2021