

# Sartopure<sup>®</sup> GA

## Superior Non-Sterile Venting Filter Cartridges



### Applications

Typical applications for Sartopure<sup>®</sup> GA air filters:

- Prefiltration in front of Sartofluor<sup>®</sup> GA membrane filters or any other membrane air filter
- Venting of non-pressure resistant vessels
- Particle removal from air streams, e.g. pressure supplies

### Product Information

Sartopure<sup>®</sup> GA offers an outstanding flow rate at low differential pressure. Sartopure<sup>®</sup> GA and Sartofluor<sup>®</sup> GA are the ideal choice for air filtration in the biopharmaceutical industry. Sartopure<sup>®</sup> GA filters expand the service life time of sterilizing grade air filter systems by removal of particles from the air stream. In addition they can be used for all venting purposes that do not necessarily require an integrity testable membrane filter.

## Retention Efficiency

The excellent retention and therefore superior protection has been proven by particle retention filtration and bacteria challenge tests (aerosol tests) performed under worst case conditions. Sartopure® GA has got a LRV of 7/cm<sup>2</sup> with Bacillus subtilis var niger spores and a LRV of 6,8/cm<sup>2</sup> with Brevundimonas diminuta. Featuring a retention of 0.2 µm for gas, Sartopure® GA efficiently protects stored products, e.g. water, liquid sugar, oral solutions etc., in the pharmaceutical industry as well as the food and beverage industry.

## Flow Rate

Due to the large filter area of 0.7 m<sup>2</sup>|10", Sartopure® GA delivers a flow rate of nearly 40 m<sup>3</sup>/h at a differential pressure of 10 mbar. This means Sartopure® GA is the perfect product for high performance filling or draining of tanks and vessels.

## Optimized Filter Material

Sartopure® GA's hydrophobic material guarantees an air flow recovery of 60 – 80% within 30 seconds after the filter has been wetted with water. The water prevents high differential pressures, ensuring fast recovery of air flow rate e. g. after cleaning the tank with hot water|agents.

## Documentation

Sartopure® GA cartridges are designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System.

## Technical Data

### Specifications

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

### Pore Size

0.2 µm (nominal in Gases)

### Available Sizes | Filtration Area

Size 1	10"	0.7 m <sup>2</sup>   7.5 ft <sup>2</sup>
Size 2	20"	1.4 m <sup>2</sup>   15.1 ft <sup>2</sup>
Size 3	30"	2.1 m <sup>2</sup>   22.6 ft <sup>2</sup>

### Available Adapters Cartridges

25, 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
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone (EPDM or Fluoroelastomer optional)

### Regulatory Compliance

Filter material bacteria challenge tested with Bacillus subtilis var niger spores and Brevundimonas diminuta

Non pyrogenic according to USP Bacterial Endotoxins

Pass USP Plastic Class VI Test

Non fiber releasing according to 21 CFR

## Retention

Aerosol tests:	LRV/cm <sup>2</sup>
Bacillus subtilis var niger:	7/cm <sup>2</sup>
Brevundimonas diminuta:	6.8/cm <sup>2</sup>

## Sterilization

### In-Line Steam Sterilization

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

### Autoclaving

134°C, 2 bar | 29 psi, 30 min

### Sterilization Cycles

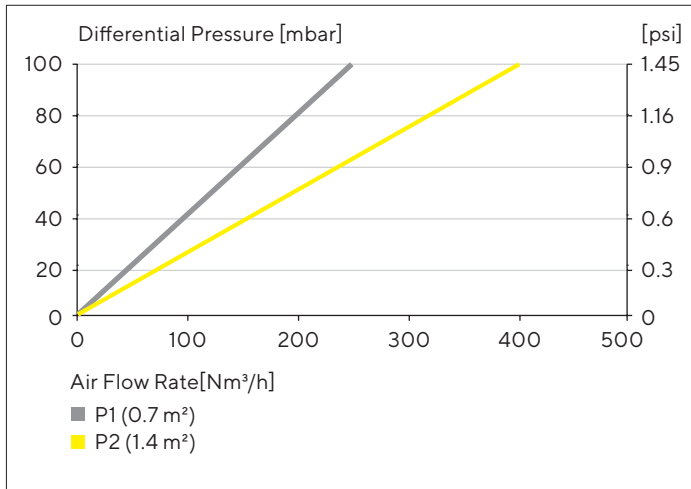
In-Line Sterilization: Min. 50

## Ordering Information

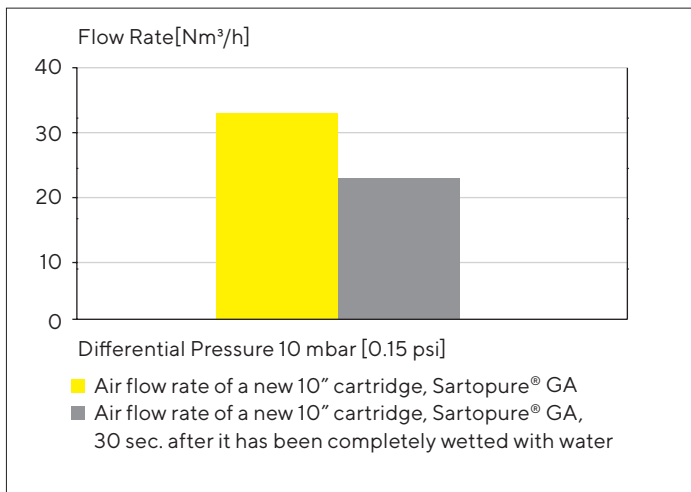
Order Code	Size	Pore Size [µm]	Pack Size [Pieces]
559**07P1--GA	1	0.2	1
559**07P2--GA	2	0.2	1
559**07P3--GA	3	0.2	1

\*\* : Adapter Type

## Air Flow Rates for 10" and 20" Cartridges



## Air Flow Recovery




**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)