

Sartopure® GA

Superior Non-Sterile Venting Filter Cartridges



Product Information

Sartopure® GA offers an outstanding flow rate at low differential pressure. Sartopure® GA and Sartofluor® GA are the ideal choice for air filtration in the biopharmaceutical industry.

Sartopure® 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.

Applications

Typical applications for Sartopure® GA air filters:

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

Introduction

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/cm² with *Bacillus subtilis var niger* spores and a LRV of 6.8/cm² 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 m² | 10", Sartopure® GA delivers a flow rate of nearly 40 m³/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 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 m² 7.5 ft²
Size 2	20"	1.4 m² 15.1 ft²
Size 3	30"	2.1 m² 22.6 ft²

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

Ordering Information

Retention

Aerosol tests: LRV/cm²
Bacillus subtilis var niger: 7/cm²
Brevundimonas diminuta: 6.8/cm²

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

Sterilization

In-Line Steam Sterilization

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

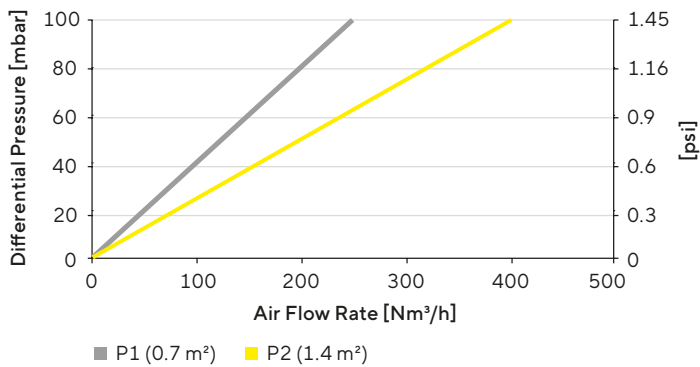
Autoclaving

134 °C, 2 bar | 29 psi, 30 min

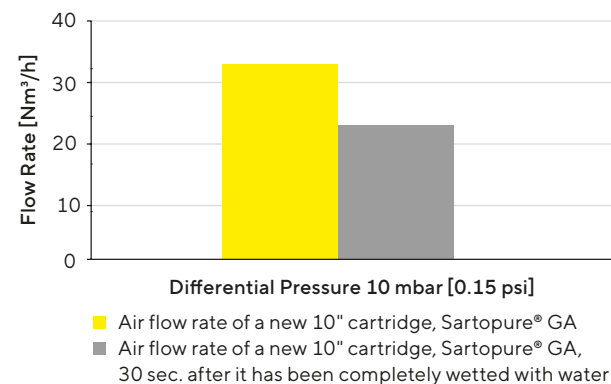
Sterilization Cycles

In-Line Sterilization: Min. 50

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 more information, visit
[sartorius.com](https://www.sartorius.com)