

# Sartopure<sup>®</sup> GA

## Superior Autoclave Venting Filter

### Product Profile

Sartopure<sup>®</sup> GA filter cartridges are specially designed for autoclave venting. The filters have a protection layer to remove dust and particles from the air stream in order to avoid blockage of the filter. Sartopure<sup>®</sup> GA meets the requirements of the BS 3970/C14. Sartopure<sup>®</sup> GA offers the highest process security, even with high gas volume stream and extreme humidity.

### Performance

Based on the experiences, Sartopure<sup>®</sup> GA, offers an optimised ratio of air flow rate and retention.

### Retention Rate

The excellent retention capabilities and therefore superior protection for autoclave chambers in the laboratories has been proven by particle retention tests and bacteria challenge tests performed under worst case scenarios. Sartopure<sup>®</sup> GA retained 10 million *Bacillus subtilis* var niger

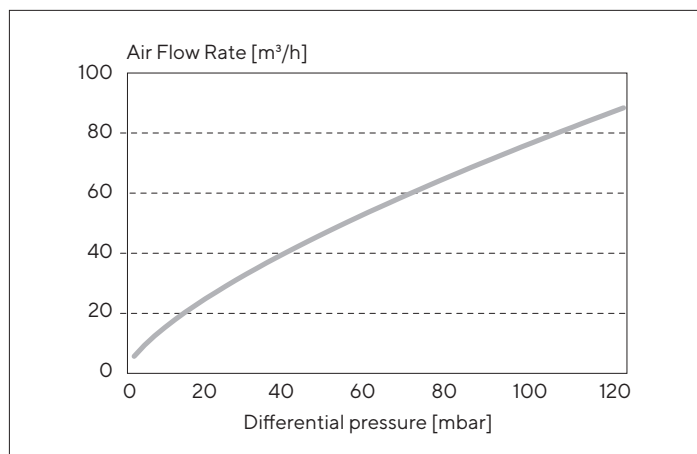


spores per cm<sup>2</sup> of filter area. Featuring a retention of 0.2 µm for gas, Sartopure<sup>®</sup> GA efficiently protects the autoclave chamber contents.

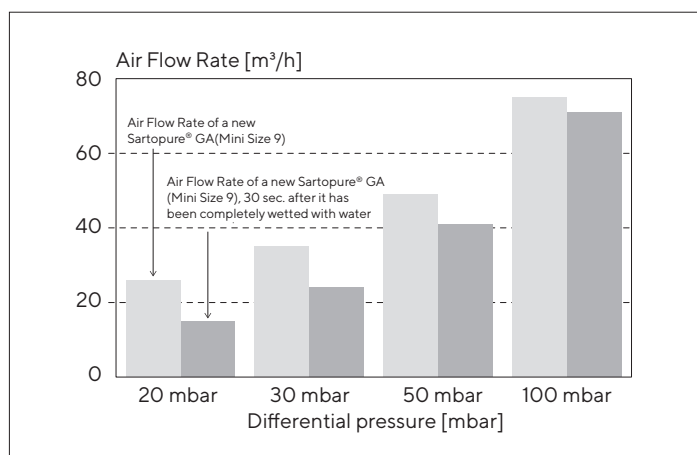
### Stability

Sartopure<sup>®</sup> GA can withstand high differential pressure either in the forward or reverse direction of flow. Its mechanical stability and the fleece structure are not affected by pulsation or high flow rates.

## Flow Rate for Mini Cartridge size 9



## Air Row Recovery



## Technical Data

Construction	hydrophobic glass fibre fleeces	
Retention rate (nominal)	0.2 µm in gases	
Filtration area	2000 cm²	
Adapter type	12 (½" Thread connector)	
Max. allowable differential pressure	max. 5 bar   73 psi at 20°C forward direction of flow max. 2 bar   29 psi at 20°C reverse direction of flow	
Sterilization	in-line steam sterilization 20 min at 134°C Autoclaving max. 134°C, 20 min	
Sterilization cycles	min. 50	
Material	Support fleece	Polypropylene
	Filtration fleece	hydrophobic glass fibre
	Core and end caps	Polypropylene
	Gaskets	Silicone

## Order Information

Order Code	Packaging
5591207P9----GAK	50 per box

### 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)