

Jumbo Star Sartopure[®] GA

The Giant-Sized
Venting Cartridge



Benefits

- Air filter cartridge with extremely high throughput
- Highest filtration area on smallest footprint

Product Information

The Jumbo Star Sartopure[®] GA cartridge is designed for all venting applications that do not necessarily require an integrity testable membrane filter cartridge.

Application

Featuring a retention of 0.2 µm for gas, Jumbo Star Sartopure® GA efficiently protects stored products in the food and beverage industry. Ideally designed for special reservoirs and storage tanks with a high capacity.

Security

The Jumbo Star Sartopure® GA filter elements ensure selective and defined particle and bioburden retention.

Performance

Jumbo Star Sartopure® GA offers an optimized ratio of air flow rate and retention as well as outstanding hydrophobicity.

Flexibility

The modular concept allows the ideal adaptation to customer requirements. Jumbo Star Sartopure® GA filter cartridges are available in two different heights.

Documentation

Jumbo Star Sartopure® GA filter cartridges are designed, developed and manufactured in accordance with a ISO 9001 Quality Management System.

Technical Data

Specifications

Materials	
Filter Material	Glass fibre and polypropylene
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone

Retention Rates

0.2 µm

Available Sizes | Filtration Area

Size 1	10"	15 m ²
Size 2	20"	30 m ²

Available Adapter

40

Operating Parameters

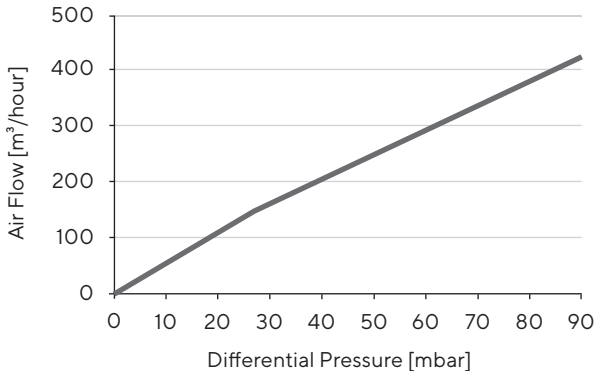
Max. Allowable Differential Pressure	4 bar 58 psi at 20°C
	1 bar 15 psi at 80°C
	0.5 bar 7.5 psi at 120°C

Max. Allowable Back Pressure	1.5 bar 22 psi at 20°C
------------------------------	--------------------------

Sterilisation

Hot Water	85 – 95°C	30 min
In-line Steaming	110°C	30 min

Air Flow Rates Jumbo Star Sartopure® GA (height 10")



Ordering Information


Order Code	Size	Retention Rate
5594007J1----GA	10"	0.2 µm
5594007J2----GA	20"	0.2 µm

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