



Jumbo Star Sartopure[®] GA

The Giant Sized Venting Cartridge

Benefits

- Air filter cartridge with extrem high throughput
- Highest filtration area on smallest footprints



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 & beverage industry. Ideally designed for special reservoirs and storage tanks with a high capacity.

Security

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

Performance

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

Flexibility

The modular concept allows the ideal adaptation to the 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	

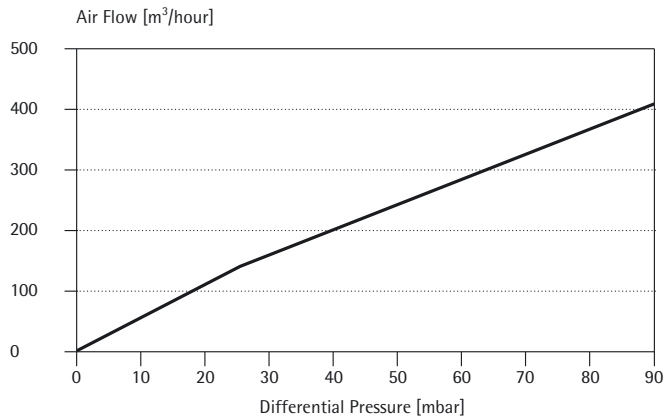
Available Sizes | Filtration Area

Size 1	10"	15 m ²
Size 2	20"	30 m ²

Available Adapter

40

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

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

Sartorius Stedim Biotech GmbH
 August-Spindler-Strasse 11
 37079 Goettingen, Germany
 Phone +49.551.308.0
 www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
 Argentina +54.11.4721.0505
 Brazil +55.11.4362.8900
 Mexico +52.55.5562.1102
 UK +44.1372.737159
 France +33.442.845600
 Italy +39.055.63.40.41
 Spain +34.913.586.098
 Russian Federation +7.812.327.53.27
 Japan +81.3.4331.4300
 China +86.21.6878.2300

Specifications subject to change without notice.
 Copyright Sartorius Stedim Biotech GmbH. Printed
 in the EU on paper bleached without chlorine.
 Publication No.: SGJ2004-e181204
 Order No.: 85030-521-72
 Ver. 12 | 2018