

Vacusart

Reliable Retention of Bacteria, Particles and Water

Product Description

Vacusart is a ready-to-connect filtration unit for reliable venting of vacuum systems. The permanently hydrophobic PTFE membrane incorporated into this unit reliably retains bacteria, particles and water from gas streams. Vacusart offers the highest possible reliability even at high flow rates and for high moisture content.

Applications

Vacusart filter units have been specially designed for reliable venting of vacuum lines and tubing. Vacusart is especially reliable as a water or liquid barrier upstream of valves or vacuum pumps.

Performance

PTFE is the most hydrophobic material available that is used in gas filtration. Even after undergoing several autoclaving cycles, the PTFE membrane still retains its outstanding hydrophobic properties.



Technical Specification

Materials

Membrane:	Reinforced PTFE
Housing:	Polypropylene
Membrane Area:	20 cm ²
Housing Diameter:	64 mm
Pore Size:	0.45 µm
Connectors:	Inlet Outlet Stepped hose barb 6-12 mm (interior thread fits luer male connector)

Operating Pressure

3 bar max. (~ 45 psi max.)

Autoclaving Temperature

134°C max.

Ordering Information

Vacusart	17804-----M
Quantity per package	3 units

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

Specifications subject to change without notice.
© 2021 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

DIR No.: 2729660-000-00 | Order No.: 85030-509-18 | Status: 03 | 22 | 2021