

Sartopure[®] IND

Particle Reduction Filter Cartridges

Product Information

- Prefiltration of beverages
- Water filtration
- Industrial applications

Applications

Sartopure[®] IND cartridges are developed for a wide range of industrial prefiltration. The all polypropylene construct contains multiple layers of pleated polypropylene fleece material making the Sartopure[®] IND cartridges an effective prefilter for particle removal and the protection of process equipment. The brought range of retention rates allows the usage throughout all process steps.

Chemical compatibility and process economy

The all polypropylene construct is resistant to pH values from 1 to 14 and a variety of chemical cleaning agents. Sartopure[®] IND cartridges ensure a selective particle retention guaranteeing an optimal filtration result and protection of process equipment. This range of cartridges packaged by 25 units reduces the overall cost for filtration and process profitability.



Performance

Sartopure® IND filter cartridges combine a high dirt holding capacity, protective abilities and extremely high flow rates. Sartopure® IND cartridges are easy to clean for an efficient regeneration if necessary.

Documentation

Sartopure® IND cartridges are designed, developed and manufactured in accordance with a ISO 9001 certified Quality Management System. All the raw materials used to produce the Sartopure® IND cartridge comply with the European Union Regulation (EC) No. 1935/2004 as well as the Regulation (EU) No. 10/2011 concerning plastic materials and articles intended to come into contact with foodstuffs.

These guidelines for plastics allow the use in food and beverage applications.

All materials used meet the requirements of the FDA.

Technical Data

Specifications

Materials	
Filter Material	Multiple Polypropylene Layers
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone

Retention Rates	
0.45 µm, 0.65 µm, 1.2 µm, 3 µm, 5 µm, 8 µm, 20 µm, 50 µm	

Available Sizes | Filtration Area

Cartridges	
10"	0.6 m ² 6.5 ft ²
20"	1.2 m ² 12.9 ft ²
30"	1.8 m ² 19.4 ft ²
40"	2.4 m ² 25.9 ft ²

Available Adapters Cartridges

21, 25, 27, 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
------------------------------	------------------------

Sartopure® IND Cartridges are free of pork DNA and drinkable alcohol.

Sterilization

In-line Steam Sterilization

110°C, 30 minutes, ΔP max 0.5 bar | 7 psi

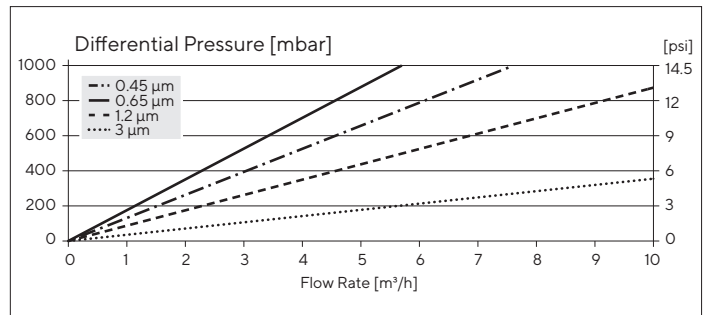
In-line hot water sterilization

85 – 95°C, 30 minutes, ΔP max 0.5 bar

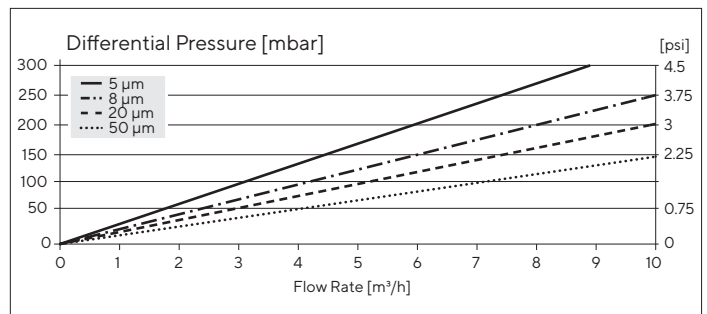
Sterilization Cycles

In-line Sterilization: minimum 100 cycles

Water Flow Rates for 10" Cartridges



Water Flow Rates for 10" Cartridges



Ordering Information

Order Code	Pore Size [μm]	Packaging [Units per box]
550**06Px-----G	0.45	25
550**05Px-----G	0.65	25
550**03Px-----G	1.2	25
550**02Px-----G	3	25
550**42Px-----G	5	25
550**01Px-----G	8	25
550**20Px-----G	20	25
550**50Px-----G	50	25
550**06Px-----C	0.45	6
550**05Px-----C	0.65	6
550**03Px-----C	1.2	6
550**02Px-----C	3	6
550**42Px-----C	5	6
550**01Px-----C	8	6
550**20Px-----C	20	6
550**50Px-----C	50	6

** : Adapter Style


x : Size

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