

Sartosteel

Removing Particles from
Liquids, Gas and Steam

Product Information

Sartosteel stainless steel depth filter cartridges contain sintered, homogeneous, 0.4 mm thick non woven stainless steel mesh layers, which are reinforced on both sides by mesh supports. These filters are used for removing particles from liquids and gases (steam). Sartosteel filter cartridges offer the user maximum security along with low filtration costs.

Applications

Sartosteel is applied in biopharmaceutical process such as:

- Steam filtration
- Condensate filtration
- Water filtration



Further Applications

Chemical Industrie

- Polymer filtration
- Catalyst retention
- Gas filtration ($\geq 80^{\circ}\text{C}$)
- Cleaning agents

Machine-building|Automotive Industries

- Fuel filtration
- Hydraulic oils

Product Benefits

- High dirt-handling capacity
- High mechanical stability
- Homogeneous material construction
- Absolutely leak-proof connections between end caps and filter unit

Quality Control

Sartosteel are designed, developed and manufactured in accordance with a DIN ISO 9001 certified Quality Management System.

Technical Data

Specifications

Materials

Sintered non woven stainless steel media, reinforced on both sides with sintered-on mesh

Filter Media AISI 316 L

Support Mesh AISI 304 | 316 L

Outer Support AISI 304 | 316 L

Core AISI 304 | 316 L

End Caps AISI 304 | 316 L

Gaskets silicone*

* standard: also available in Fluoroelastomer and EPDM

Retention Rate

3 µm

Available Sizes | Filtration Area

Size 9 (Mini Cartridge) 9" 200 cm² | 0.2 ft²

Size 1 10" 500 cm² | 0.5 ft²

Size 2 20" 1.000 cm² | 1.0 ft²

Size 3 30" 1.500 cm² | 1.5 ft²

Operation Parameters

Max. differential pressure:

≤ 20 bar, in the direction of filtration

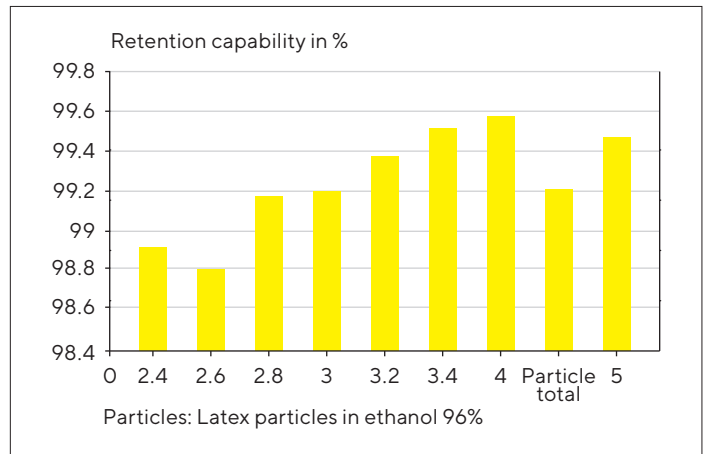
≤ 1 bar, opposite to the direction of filtration

Resistance | Compatibility

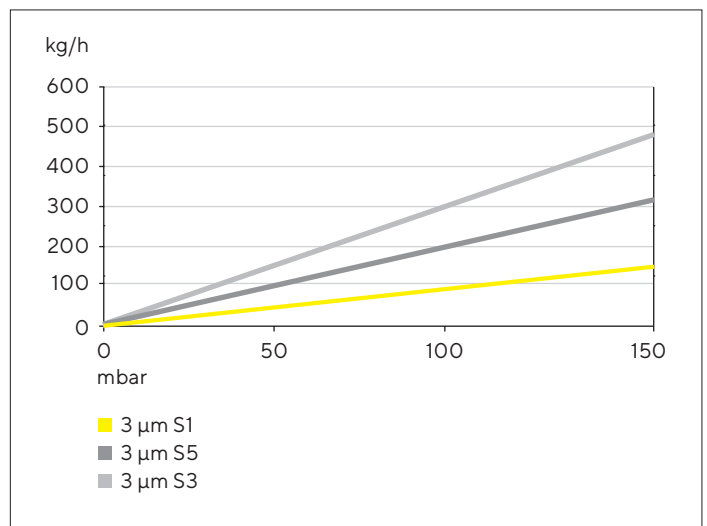
thermal up to 200°C (not with silicone sealing)

chemical inert to caustic solutions, solvents not compatible with aggressive and relatively high concentrations of acid (≥ 5%)

Particle Retention Rating 3 µm Sartosteel



Sartosteel – Sarturated Steam (T = 121°C, P = 1 Bar System Pressure)



Ordering Information

Order Code	Retention Rate [μm]	Adapter Type	Height	Pack Size [Pieces]
Mini Cartridge				
5701502S9	3	15*	9"	1
Cartridges				
5702502S1	3	25	10"	1
5702502S5	3	25	20"	1
5702502S3	3	25	30"	1
5702802S1	3	28	10"	1
5702802S5	3	28	20"	1
5702802S3	3	28	30"	1


* Bayonet adapter with internal O-ring

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