

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.



Features and Benefits

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

Introduction

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 ($\geq 80\text{ }^{\circ}\text{C}$)
- Cleaning agents

Machine-Building | Automotive Industries

- Fuel filtration
- Hydraulic oils

Quality Control

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

Technical 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 cm^2 0.2 ft^2
Size 1	10"	500 cm^2 0.5 ft^2
Size 2	20"	1,000 cm^2 1.0 ft^2
Size 3	30"	1,500 cm^2 1.5 ft^2

Operation Parameters

Max. differential pressure:

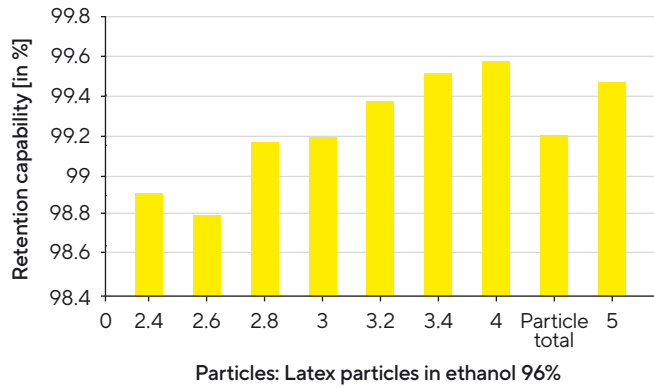
- ≤ 20 bar, in the direction of filtration
- ≤ 1 bar, opposite to the direction of filtration

Resistance | Compatibility

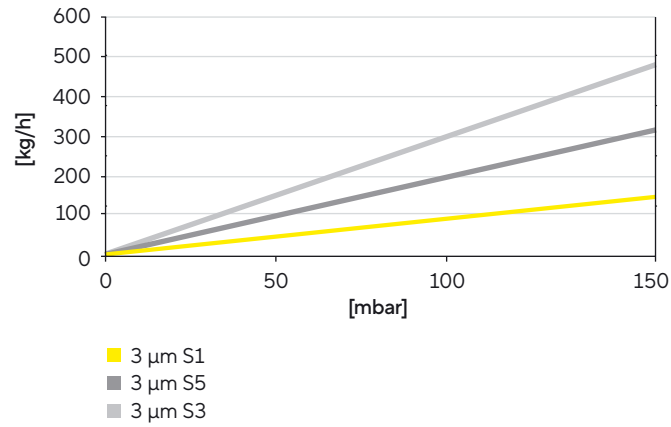
Thermal	Up to 200 $^{\circ}\text{C}$ (not with silicone sealing)
Chemical	Inert to caustic solutions, solvents not compatible with aggressive and relatively high concentrations of acid ($\geq 5\%$)

Ordering Information

Particle Retention Rating 3 μm Sartosteel[®]



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



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 more information, visit

[sartorius.com](https://www.sartorius.com)