



# Filtration Solutions

For Food and Beverage  
Applications Only

Simplifying Progress

**SARTORIUS**

# Shift your F & B Process to a New Level of Quality

Specifically designed for your requirements, Sartorius offers a broad range of filtration solutions to make your processes faster, easier and more economical. With flow rates from 100 L/hr to 100,000 L/hr or even higher, our fully automated systems ensure:

- Significantly reduced filtration cost
- Enhanced flow rates
- Increased total throughput
- Optimized bioburden
- Stabilized process flow
- Minimized non-productivity and downtime

## Areas of Application

The broad range of products fulfills all requirements for the filtration of:

- Wine | Sparkling wine
- Beer
- Water
- Alcohol
- Soft drinks
- Gases | Steam

## Worldwide Services

### Support

- On-site assistance
- Filtration optimization trials
- Training seminars
  - Filtration technology
  - Process designing
  - Validation aspects

### Quality Assurance

- Compliant with relevant FDA and EU Regulations for food contact (refer to Validation Guide)
- Every membrane filter is integrity tested to ensure its 100% integrity

# Our Portfolio for Your Applications

## Prefilter Cartridges

The broad variety of different prefilter materials combine unmatched total throughput performance with a level of clarification as never seen before. Increase the total throughput of your final filtration run and protect your processes from premature blockage with the right choice of prefilter.



### Jumbo Star Cartridges

Unique Jumbo filter cartridges for high-volume flow rates and maximum throughput and featuring the smallest footprint.

## Membrane Filter Cartridges

Wherever and whenever 100% retention ratings for beverage-spoilage microorganisms are essential: These membrane filters are generally used in the final filtration stage and installed directly upstream of the filling unit. Sartorius Stedim Biotech supplies the following membrane types for filtration of liquids and gases:

- Polyethersulfone (PES)
- Polytetrafluoroethylene (PTFE)

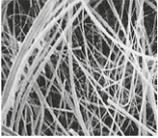
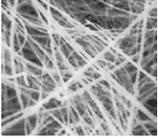
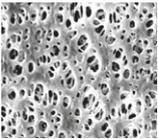
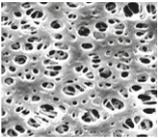
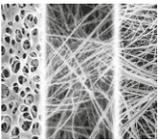


## Integrity Testing

Ensure 100% reliable filling by testing all final membrane filters with our testing system Sartocheck® mini.

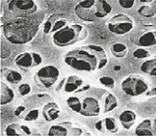
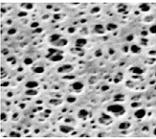
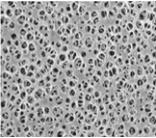
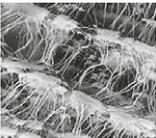
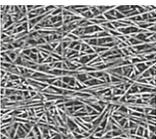


# Technical Specifications

	Applications	Cartridge Construction	Available Heights	Adapter Types (**)
<p><b>Sartopure® IND</b> The prefilter cartridge with particle-removing with polypropylen filter material</p> 	Retention of particles; reduction of microorganisms by fractionized filter fleeces	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ Nonwoven polypropylene filter layers</li> <li>▪ Nonwoven polypropylene drainage layer (heat-sealed, non-fiber-releasing)</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> <li>▪ Delivered in packages of 25 cartridges</li> </ul>	10", 20", 30", 40"	21, 25, 27, 28
<p><b>Sartopure® GF Plus</b> The fleece based filter cartridge with adsorptive nonwoven glass fiber materials for maximum protection</p> 	Retention of particles and colloids, reduction of microorganisms by fractionized filter fleeces	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ Nonwoven glass fiber layers</li> <li>▪ Nonwoven polypropylene drainage layer (heat-sealed, non-fiber-releasing)</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> </ul>	10", 20", 30", 40"	21, 25, 27, 28
<p><b>Sartocool® PS</b> The high-performance membrane cartridge for cold sterilization of beer</p> 	Retention of yeasts and beer-spoilage bacteria	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ Single-layer PES membrane</li> <li>▪ Nonwoven polypropylene drainage layer</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> </ul>	10", 20", 30", 40"	25, 27, 28
<p><b>Vinosart® PS</b> The high-performance membrane filter cartridge for wine and sparkling wine filtration</p> 	Specially developed for the retention of microorganisms, particles and colloids in all types of wine and sparkling wine	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ Single-layer PES membrane</li> <li>▪ Nonwoven polypropylene drainage layer</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> </ul>	10", 20", 30", 40"	25, 27, 28
<p><b>Jumbo Star</b> The large-area filter cartridge for easy handling and high flow rates</p> 	Retention of particles; reduction of colloids and bioburden in water, wine and beer and in alcohol production; also for venting large tanks	Pleated construction with filter areas of up to 30 m <sup>2</sup> ; pleats made of polypropylene or glass fiber material*	10", 20", 30", 40"	40

\* Polypropylene 559xxx | Glass fiber 555xxx

The pictures shown are for reference only, the actual product may differ.

	Applications	Cartridge Construction	Available Heights	Adapter Types (**)
<b>Sartobev® PS</b> The classic membrane filter cartridge for filtration of wine and sparkling wine 	For the retention of microorganisms, particles and colloids in wine, sparkling wine and water	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ PES membrane</li> <li>▪ Nonwoven polypropylene drainage layer</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> </ul>	20", 30", 40"	25, 28
<b>Aquasart® PS</b> The high-performance membrane filter cartridge for water filtration 	For the retention of microorganisms, particles and colloids in mineral water	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ Single-layer PES membrane</li> <li>▪ Nonwoven polypropylene drainage layer</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> </ul>	10", 20", 30", 40"	25, 27, 28
<b>Aquasart® Plus</b> The innovative membrane filter cartridge for liquids to guarantee the lowest filtration costs 	Best retention of microorganisms, particles and colloids in mineral water as well as near-water beverages	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ Double layer PES</li> <li>▪ Nonwoven polypropylene drainage layer</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone</li> </ul>	20", 30", 40"	25, 28
<b>Aerosart®</b> The hydrophobic PTFE membrane filter cartridge 	Sterile venting of tanks fermenters and bioreactors. Sterile filtration of inlet and outlet air, gases, solvents and aggressive media	<ul style="list-style-type: none"> <li>▪ Protective nonwoven polypropylene layer</li> <li>▪ PTFE membrane</li> <li>▪ Nonwoven polypropylene drainage layer</li> <li>▪ Outer support, core and end caps: polypropylene</li> <li>▪ O-rings: silicone or EPDM</li> <li>▪ Optional: fluoroelastomer</li> </ul>		25, 27, 28
<b>Sartosteel®</b> Stainless steel filter cartridge for filtration of steam 	For the removal of particles from steam and gases	<ul style="list-style-type: none"> <li>▪ Sintered nonwoven stainless steel mesh: AISI 316L</li> <li>▪ Reinforced on both sides with sintered-on mesh filter support: AISI 316L</li> <li>▪ Core and outer support: AISI 316L</li> <li>▪ O-rings: silicone</li> </ul>	10", 20", 30"	21, 25, 27, 28

The pictures shown are for reference only, the actual product may differ.

# Ordering Information

	Retention Rate	Height 1 (10")	Height 2 (20")	Height 3 (30")	Height 4 (40")
<b>Sartopure® IND</b>	0.45 µm	550**06P1	550**06P2	550**06P3	550**06P4
	0.65 µm	550**05P1	550**05P2	550**05P3	550**05P4
	1.2 µm	550**03P1	550**03P2	550**03P3	550**03P4
	3 µm	550**02P1	550**02P2	550**02P3	550**02P4
	5 µm	550**42P1	550**42P2	550**42P3	550**42P4
	8 µm	550**01P1	550**01P2	550**01P3	550**01P4
	20 µm	550**20P1	550**20P2	550**20P3	550**20P4
	50 µm	550**50P1	550**50P2	550**50P3	550**50P4
<b>Sartopure® GF Plus</b>	0.65 µm	555**05P1	555**05P2	555**05P3	555**05P4
	1.2 µm	555**03P1	555**03P2	555**03P3	555**03P4
<b>Sartocool® PS</b>	0.45 µm	530**06A1	530**06A2	530**06A3	530**06A4
	0.65 µm	530**05A1	530**05A2	530**05A3	530**05A4
<b>Vinosart® PS</b>	0.45 µm	542**06A1	542**06A2	542**06A3	542**06A4
	0.65 µm	542**05A1	542**05A2	542**05A3	542**05A4
<b>Jumbo Star</b>	0.45 µm	-	5504006J2	5504006J3	5504006J4
	0.65 µm	5594005J1	5594005J2	5594005J3	5594005J4
		5554005J1	5554005J2	5554005J3	5554005J4
	1.20 µm	5594003J1	5594003J2	5594003J3	5594003J4
		5554003J1	5554003J2	5554003J3	5554003J4
	3.00 µm	5594002J1	5594002J2	5594002J3	5594002J4
		5634002J1	5634002J2	5634002J3	5634002J4
	5.00 µm	5594042J1	5594042J2	-	5594042J4
8.00 µm	5594001J1	5594001J2	5594001J3	5594001J4	
20.0 µm	5594020J1	5594020J2	5594020J3	5594020J4	

	Retention Rate	Height 1 (10")	Height 2 (20")	Height 3 (30")	Height 4 (40")
Sartobev® PS	1.00 µm	-	531**24A2	531**24A3	531**24A4
Aquasart® PS	0.2 µm	541**07A1	541**07A2	541**07A3	541**07A4
Aquasart® Plus	0.2 µm	-	538**07H2	538**07H3	538**07H4
	0.1 µm	-	-	538**58K3	-
Aerosart®	0.2 µm	515**07T1---E--C	515**07T2---E--C	515**07T3---E--C	-
Sartosteel®	3 µm	570**02S1	-	570**02S3	-

Replace the \*\* in the order number with the code for the adapter of your choice.

**Pack Size**  
blank box of 1  
C box of 6  
G box of 25

**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](http://www.sartorius.com)