



# Sartocool PS (0.65 µm)

Filter Cartridge for Particle & Yeast Cell Retention



## Description

Sartocool PS filter cartridges feature a high-performance, single-layer pleated PES membrane. The ideal choice for the filtration in the food and beverage industry, especially developed for the filtration of beer and sparkling wine. Sartocool PS (0.65 µm) guarantee the retention of particles and yeast cells. Enhanced total throughput performance in combination with highest product safety is achieved.

## Application

Sartocool PS membrane filter cartridges are used for the final filtration prior to bottling or as an integrity testable prefilter for such beverages as the following:

- Beer
- Wine
- Sparkling Wine

## Product Performance

Sartocool filter cartridges incorporate a PES membrane. This new polyethersulfon membrane has been developed to meet the demanding process requirements in terms of pressure pulsation and sterilization cycles in the food and beverage industry. Sartocool PS achieves maximum total throughput to minimize the filtration costs. The membrane is chemically resistant within the pH-range of 1–14. Caustic soda and acids can be used for a chemical regeneration.

## Integrity-testable and Quality

The integrity of Sartocool PS membrane filters cartridges can be easily checked by performing an integrity test before and after filtration to guarantee the highest biological safety. This membrane filter is designed and manufactured according to an ISO 9001 certified quality management system. All materials used in the construction of Sartocool PS meets FDA requirements for use in food and beverage applications.

## Technical Specifications

### Material

Filter medium	Single-layer PES membrane
Protective non-woven drainage layers	Polypropylene
Core   outer support   end caps   adapter	Polypropylene
O-rings	Silicone

### Pore Size

0.65 µm

### Filtration Area

0.78 m<sup>2</sup> per 10" element

### Available Sizes | Filtration Area

Size 1	10"	0.78 m <sup>2</sup>
Size 2	20"	1.56 m <sup>2</sup>
Size 3	30"	2.34 m <sup>2</sup>
Size 4	40"	3.12 m <sup>2</sup>

### Available Adapter

25, 27, 28

### Operating Parameters

Maximum differential pressure	5 bar   73 psi at 20°C
	2 bar   29 psi at 80°C
	0.5 bar   7 psi at 121°C
Maximum back pressure	2 bar   29 psi at 20°C

### Technical Specifications

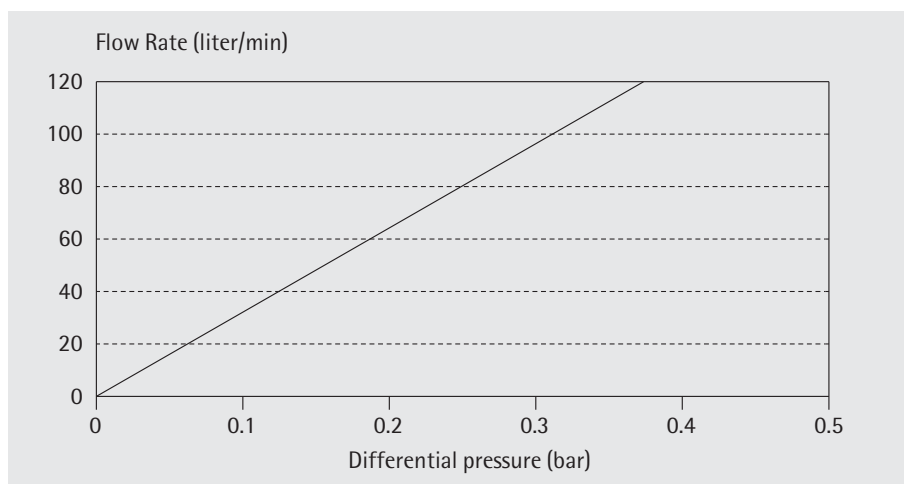
Recommended flow rate	10 to 12 hl beer/h/30" 3.000 liter water/h/30"
Sterilization	Hot water (85-95°C) or Steam (110°C), at least 100 cycles
Microbiological recovery rate	Saccharomyces cerevisiae > 10 <sup>7</sup> /cm <sup>2</sup>
ph stability	ph 1-14

### Integrity Test Data

Test pressure	0.8 bar
Max. diffusion (10" element)	8 ml
Test time	10 minutes
Bubble point	> 0.8 bar

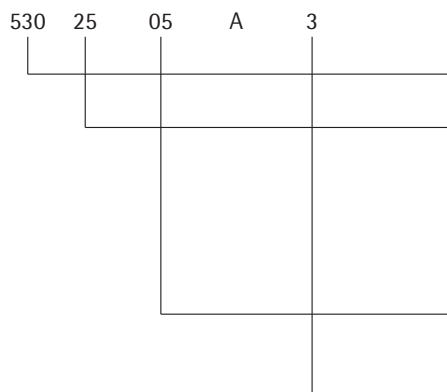
### Flow rate for water

Sartocool PS (0.65 µm) height 10"



### Ordering Information

#### Cartridges



#### Sartocool PS

#### Adapter

- 25 S-adapter on top, 2-flange bayonet adapter with double O-ring on bottom
- 27 Double O-ring
- 28 S-adapter on top, 3-flange bayonet adapter with double O-ring on bottom

#### Retention Rate

- 05 0.65 µm

#### Height

- 1 10"
- 2 20"
- 3 30"
- 4 40"

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen, Germany  
Phone +49.551.308.0  
Fax +49.551.308.3289  
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178  
UK +44.1372.737159  
France +33.442.845600  
Italy +39.055.63.40.41  
Spain +34.91.3586102  
Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH  
W/sart-000 - G  
Publication No.: SPK2086-e08051  
Order No.: 85034-535-34