# SARTURIUS

# 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  $\mu m$ ) 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 minimizes 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 m² 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**

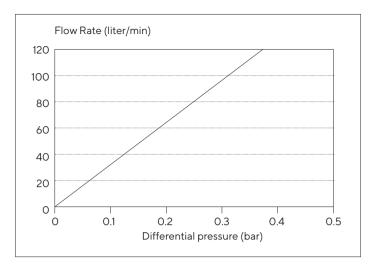
25, 27, 28

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
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	Saccharomcyes 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

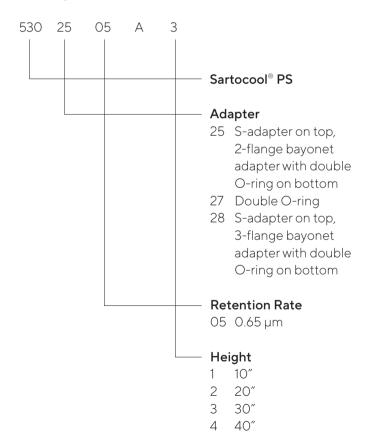
#### Flow rate for water

Sartocool® PS (0.65 µm) height 10"



# Ordering Information

## Cartridges



## 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