

# Aquasart® PS 1 µm

Membrane Filter Cartridge  
for the Food & Beverage  
Industry



## Product Information

Aquasart® PS filter cartridges are integrity testable membrane filters especially developed for filtration of water in the Food & Beverage Industry. They are characterised by a unique hydrophilic Polyethersulfone membrane offering a broad chemical compatibility, exceptional high flow rates and total throughputs.

## Applications

Aquasart® PS filter cartridges can be used especially in applications which require a chemical disinfection or cleaning with caustic or acid, due to the broad chemical compatibility from pH 1 to pH 14.

## Total Throughput

The asymmetric pore structure of the Polyethersulfone membrane of the Aquasart® PS filter cartridges provides outstanding total throughputs and high flow rates. Economic system design for filtration processes is assured. The total throughput of the final filters can be significantly increased by an optimised prefilter system reducing the filtration costs overall.

## Flow Rates

The high effective filtration area of Aquasart® PS filter cartridges allows highest flow rates and economic filtration system design.

## Stability

Aquasart® PS filter cartridges withstand more than 5.000 pulsation cycles with differential pressure up to 5 bar | 75 psi and multiple steam sterilisation cycles up to 134°C without loss of filter integrity.

## Safety

Aquasart® PS filter cartridges membrane can be checked before and after every use with an integrity test.

## Quality Control

Each individual element is tested for integrity during production assuring highest product quality.

## Documentation

Aquasart® PS filter cartridges are designed, developed and produced under a ISO 9001 certified Quality Management System.

## Technical Data

### Specifications

---

Materials	
Membrane	Highly asymmetric, single layer, hydrophilic Polyethersulfone membrane
Fleeces	Polypropylene
Outer Cage	Polypropylene
Inner Core	Polypropylene
Endkaps	Polypropylene
O-Rings	Silicone

---

### Pore Size

1 µm absolute

---

### Available Sizes | Filtration Area

---

Size 3	30"	2.1 m <sup>2</sup>   22.6 ft <sup>2</sup>
--------	-----	---

---

### Available Adapters

25, 27, 28

### Operating Parameters

---

Max. Allowable Differential Pressure	5 bar (72.5 psi) at 20°C 2 bar (29 psi) at 80°C
Max. Allowable Back Pressure	2 bar (29 psi) at 20°C
Sterilisation	Hot Water 85–95°C, 30 min, Δp max. 0.5 bar   7 psi In-Line-Steam at 110°C, 30 min, Δp max. 0.5 bar   7 psi
Sterilisation cycles	min. 100

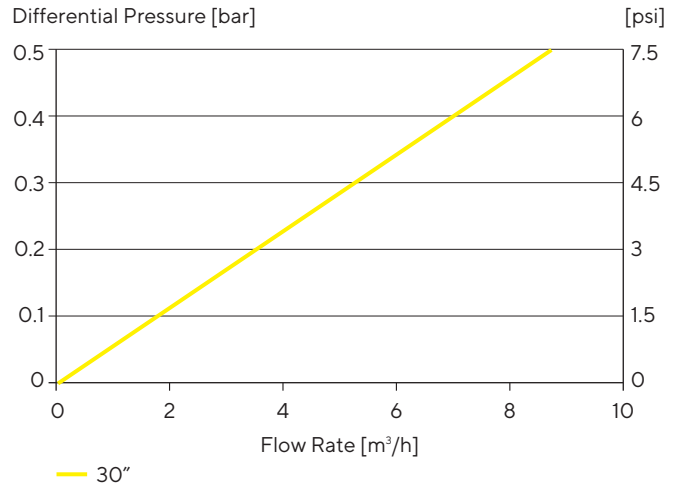
---

## Compatibility

Aquasart filter cartridges meet the requirements laid down in the Framework Regulation (EC) N° 1935/2004 and Commission Regulation (EU) N° 10/2011 of 14th January 2011 on plastic materials and articles intended to come into contact with food. These guidelines for plastics allow the use in food and beverage applications. All materials used meet the requirements of the FDA. pH-Compatibility: pH 1 to pH 14.

Aquasart® PS filter cartridges are free of pork DNA and drinkable alcohol.

## Water Flow Rates for 30" Cartridges



## Ordering Information


Order Code	Pore Size [µm]	Filtration area [m²]
541**24A3	1	2.1

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Straße 11  
37079 Göttingen  
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)