



# Aqua Star CA

Filter Cartridge for Particle & Bioburden Reduction



## Description

Aqua Star CA filter cartridges feature a high-performance, single-layer pleated cellulose acetate membrane. Specially developed for filtration of carbonated mineral water and non-carbonated drinks. Aqua Star CA filter cartridges guarantee the highest biological product safety in combination with an enhanced total throughput performance. Aqua Star CA filter cartridges are the ideal choice for filtration in the food and beverage industry.

## Application

Aqua Star CA membrane filter cartridges are used for final filtration prior to bottling for products, such as the following:

- Carbonated mineral water
- Non-carbonated water
- Near-water beverages

## Product Performance

Aqua Star CA filter cartridges incorporate a specially designed cellulose acetate membrane. This new cellulose acetate membrane has been developed to meet the demanding process requirements in terms of pressure pulsation and sterilization cycles in the food and beverage industry. In combination with an optimally designed prefilter, Aqua Star CA filter cartridges achieve maximum total throughput, and minimize filtration costs.

## Integrity-testable

The integrity of Aqua Star CA membrane filters cartridges can be easily checked by performing an integrity test before and after filtration to guarantee the highest biological safety.

## Quality

Aqua Star CA membrane filters are designed and manufactured according to an ISO 9001 certified quality management system. All materials used in the construction of Aqua Star CA meet FDA requirements for the use in food and beverage applications.

## Technical Specifications

Filter medium	Single-layer cellulose acetate membrane
Protective non-woven drainage layers	Polypropylene
Core   outer support   end caps   adapter	Polypropylene
O-rings	Silicone
Filtration area	0.7 m <sup>2</sup> per 10" element
Pore size	0.2 µm
Heights	10" (250 mm), 20" (500 mm), 30" (750 mm), 40" (1,000 mm)
Maximum differential pressure	5 bar   73 psi at 20°C 2 bar   29 psi at 80°C
Maximum back pressure	2 bar   29 psi at 20°C
Recommended flow rate	Approx. 3.000 liter/per hour/ per 30"

## Technical Specifications

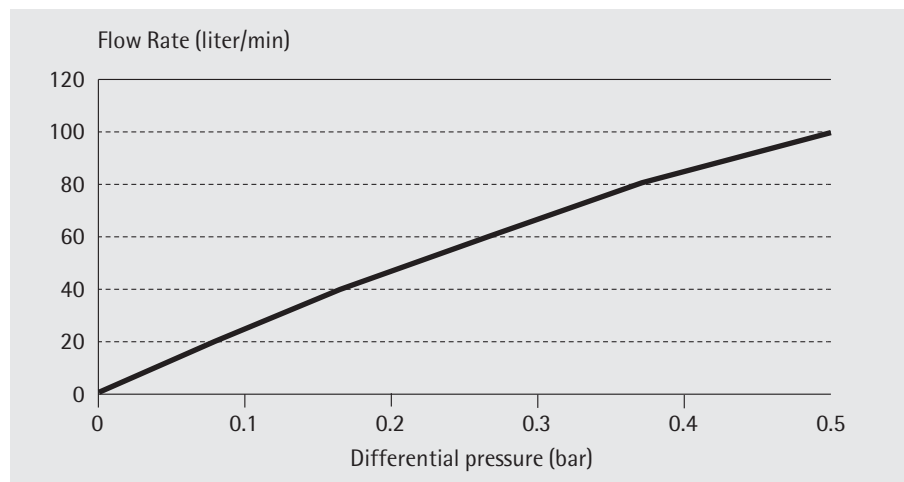
### Sterilization

Hot water	85 - 95°C, 30 minutes ΔP max. 0.5 bar
Steam	110°C, 30 minutes ΔP max. 0.5 bar for 100 cycles

### Integrity Test Data

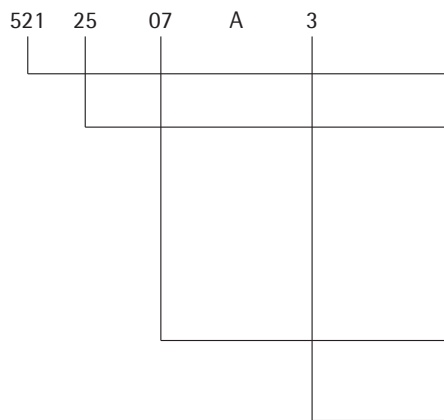
Pore size	0.2 μm
Test pressure	2.3 bar
Max. diffusion (10" element)	13 ml
Test time	10 minutes

## Flow rate for water Aqua Star CA (height 30")



## Ordering Information

### Cartridges



### Aqua Star CA

#### Adapter

- 21 Double open end
- 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

- 07 0.20 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 +34.90.2110935  
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.: SPK2077-e10073  
Order No.: 85030-523-59  
Ver. 07 | 2010