

Sartovent P

Protection for Sanitary Tanks

Sartovent P Air Filter: Maximum Filtration Reliability

Sartovent is completely made of polypropylene. And naturally it won't affect the taste or odor of your products because it's approved by food safety laws.

It Ensures That Your Products Will Remain Microbe-Free

- Basic ingredients
- Fruit juices
- Liquid sugar
- Unsweetened must for wine



“Microbe-Proof” Design

Thanks to its special construction, Sartovent retains particles and microbes reliably.

Heat Stable

Sartovent can be steam-sterilized while mounted in place on top of a container or autoclaved separately.

Leaktight Fit

A few turns with a common hook wrench are all it takes to securely install Sartovent on a tank, thus ensuring a leaktight, secure fit.

Wide Selection of Beverage Threads

Sartovent is available in various beverage thread sizes to fit most tanks on the market: DN 25, 32, 40.

Specifications

Materials	Polypropylene (PP) housing Nonwoven PP filter material
Retention rating	0,2 µm, nominal
Connectors	Beverage thread acc. to DIN 11.851, DN 25, DN 32, DN 40
Height	80 mm
Outer diameter	110 mm Ø

No Investment Costs

Sartovent eliminates the need for conventional stainless steel housings with filter elements. Why? Sartovent is an all-in-one unit made of plastic – unbeatably affordable.

Tamper-Proof

Sartovent is lead-sealed to the tank to prevent unauthorized removal that may go unnoticed.

Practical Design

The air inlet and outlet openings are configured to ensure optimum flow geometry in the filter medium. The integrated splash guard prevents premature blockage of the membrane by stopping liquid from splashing the pleated filter material, especially during transportation of containers.

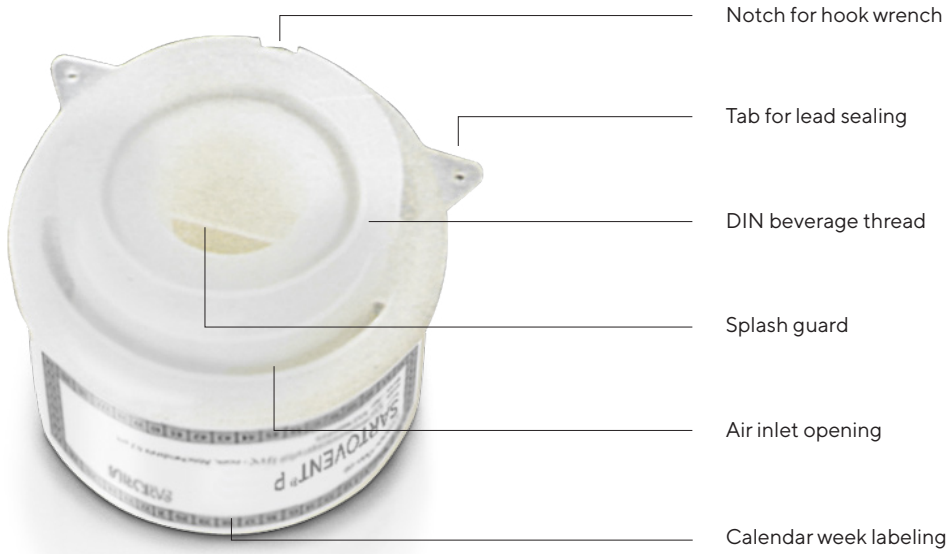
Convenient Label for 100% Traceability

Sartovent's convenient label has numbered squares representing the calendar weeks so you can keep track of its initial use by marking the appropriate square. Moreover, just one glance at the label and you'll know immediately when Sartovent has to be reesterilized or exchanged.



Order Information

5521607P25-----B	DN 25 Thread
5521607P32-----B	DN 32 Thread
5521607P40-----B	DN 40 Thread




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