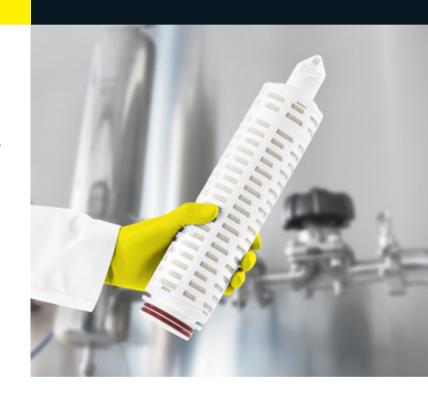
SARTURIUS

Sartofluor® HR

Heat Resistant Air | Gas Filter



Product Information

Sartofluor® has been synonomous with reliable air | gas filtration for decades. Employing the proven single-layer ePTFE membrane technology, this latest addition to the Sartofluor® broad offering of filter cartridges is specifically designed for long-term, high-temperature applications, such as Water For Injection (WFI) loop venting. Bacterial, phage and particle retention validation make Sartofluor® HR a valuable addition to the Sartofluor® product family.

Designed for High Temperature Applications

Sartofluor® HR filter cartridges can be used for an extended time period of at least six months at a temperature of 90°C. Incorporating the proven hydrophobic 0.2 μ m Sartofluor® ePTFE membrane (sterilizing grade according to current ASTM F-838 guideline), Sartofluor® HR also provides a validated retention of particles (sizes 0.005 μ m | 0.3 μ m), which is of the utmost importance for these types of applications. Integrity testing by Water Intrusion Test (WIT) allows in-situ integrity testing and circumvents the use of alcohol for wetting. Multiple in-line steaming | autoclaving cycles can be conducted without any loss of performance.

Applications

- Water For Injection (WFI) tanks | loops (80 90°C)
- Autoclaves
- Permanently heated filter housings
- All applications where hot air (approx. 90°C) is used over long periods

Inherent Quality Provides Full Confidence

The 0.2 µm hydrophobic ePTFE membrane is self dewetting and, therefore, provides high air flow rates even at very low differential pressures over an extended process time. Furthermore, the polyphenylene sulfide (PPS) drainage fleece is resistant to ozonated air.

Validated by various retention tests, Sartofluor® HR filter cartridges ensure sterile filtration of air | gas. The following indispensable requirements are fulfilled:

- Retention of ≥ 10⁷ Brev. diminuta/cm²
- Retention of ≥ 10⁷ MS-2 coliphage/cm²
- Retention of particles: sizes 0.005 µm | 0.3 µm
- Clear labeling with filter type, lot number, individual unit serial number, 2D data matrix code
- Developed, produced and distributed under a quality management system certified to DIN/ISO 9001.
- USP Plastic Class VI Test

Technical Data

Available Sizes	Filtration Area [m² ft²]	Max. WIT @ 2.5 bar 36 psi [mL/10 min]	Max. WFT @ 2.5 bar 36 psi [mL/min]	Max. Diffusion @ 0.7 bar 10 psi [mL/min]*	Min. Bubble Point [bar psi]*
Cartridges					
Size 1 (10")	0.56 6.03	12	0.34	10	1 14.5
Size 2 (20")	1.12 12.06	24	0.68	20	1 14.5
Size 3 (30")	1.68 18.09	36	1.03	30	1 14.5

^{*} Wetting agent IPA | Water (60 | 40).

Max. Differential Pressure for Air

In the direction of filtration:

Temperature [°C]	20	90
Pressure [bar]	5	2
Pressure [psi]	72.5	29

In the reverse direction of filtration:

Temperature [°C]	20	90
Pressure [bar]	2	1.5
Pressure [psi]	29	22

Sterilization

In-Line Steam Sterilization

134°C, 30 min. at max. differential pressure of 0.5 bar | 7.25 psi in forward and reverse direction Max. 100 Sterilization Cycles

or

Autoclaving

134°C, 2 bar, 30 min Max. 100 Sterilization Cycles

Technical Reference

Validation Guide SPK5814-e | 85037-554-85

Materials

Pore Size

Membrane Filter

0.2 µm

Single layer hydrophobic expanded polytetrafluoroethylene (ePTFE)

Drainage Layer

Polyphenylene Sulfide (PPS)

Core

Polypropylene (PP)

Outer Support

Polypropylene (PP)

End Caps

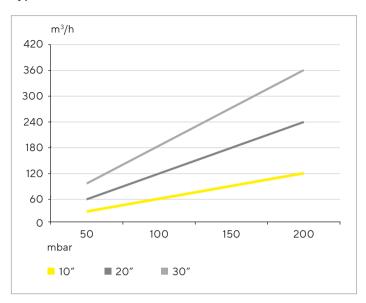
Polypropylene (PP)

O-Rings

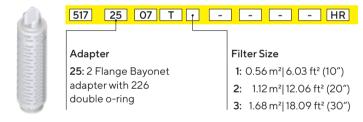
Silicone

Performance

Typical Air Flow



Ordering Information



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