



F&B Filter Housings

For the Food and Beverage Industry

Product Information

Sartorius Stedim Biotech F&B Filter Housings are specifically designed for liquid filtration applications of the Food and Beverage Industry. Manufactured in an PED 2014/68/EU certified facility, special attention has been paid to choice of materials, durability, cleanability, ease of use and quality control.



Introduction

Sartorius Stedim Biotech F&B Filter Housings are made to support the high expectations and standards of our customers.

Applications

Sartorius Stedim Biotech F&B Filter Housings are ideally suited for liquid filtration, including:

- Beer
- Wine & spirits
- Mineral | Bottled water
- Carbonated Soft drinks
- Non-carbonated Soft drinks

Quality of Material

Only 316L grade stainless steel is used for all wetted surfaces to provide maximum durability and resistance to corrosion. All supplied gaskets and O-rings meet FDA regulatory requirements.

Quality of Surface Finishes

Sartorius F&B Filter Housings come standard with internal finishes of at least <math><32 \mu\text{inch}</math> (0.8 μm) Ra mechanical polish.

Ease in Cleaning

Thanks to their optimized design, our F&B housings are exceptionally easy to clean for thorough results and can be in-line steam sterilized as a complete unit, incl. gaskets.

Quality Control and Documentation

Standard documentation includes: User manual, Declaration of Conformity, Certificate for Manufacture and Pressure Testing.

Ease of Installation

Sartorius Stedim Biotech Filter Housings are sold ready-to-install with the gasket(s), O-Ring(s) and clamp included.

CE PED 2014/68/EU Standard

Sartorius Stainless Steel Housings are designed and manufactured according the Pressure Equipment Directive PED 2014/68/EU. Our manufacturing process follows the highest quality standards and is monitored by an internal quality control system as well as by independent notified bodies on a regular basis.

Technical Data

Technical Specifications

Product Contact Surfaces	316L
Gasket Materials	Silicone (Viton or EPDM available)
CFR Compliance	All gasket materials comply with the FDA regulations 21 CFR 17.2600
Closure System	Clamp closure
Height	10", 20", 30", 40"
Adapter	25 Code 7 (-226)
Surface Finishes	Ra <0.8 µm internal MP, Ra <1.6 µm external MP

Ordering Information

Standard Order Codes	Basic Description	In Out TC	Vent TC	Drain TC
810011G25ATT010L	1 × 10" F&B, T-type, Cd 7, Silicone	1"	1.5"	1"
810012G25ATT010L	1 × 20" F&B, T-type, Cd 7, Silicone	1"	1.5"	1"
810013G25ATT010L	1 × 30" F&B, T-type, Cd 7, Silicone	1"	1.5"	1"
810014G25ATT010L	1 × 40" F&B, T-type, Cd 7, Silicone	1"	1.5"	1"
810032G25ATT015L	3 × 20" F&B, T-type, Cd 7, Silicone	1.5"	1.5"	1.5"
810033G25ATT015L	3 × 30" F&B, T-type, Cd 7, Silicone	1.5"	1.5"	1.5"
810034G25ATT015L	3 × 40" F&B, T-type, Cd 7, Silicone	1.5"	1.5"	1.5"
810052G25ATT020L	5 × 20" F&B, T-type, Cd 7, Silicone	2"	1.5"	1.5"
810053G25ATT020L	5 × 30" F&B, T-type, Cd 7, Silicone	2"	1.5"	1.5"
810054G25ATT020L	5 × 40" F&B, T-type, Cd 7, Silicone	2"	1.5"	1.5"
810082G25ATT025L	8 × 20" F&B, T-type, Cd 7, Silicone	2.5"	1.5"	1.5"
810083G25ATT025L	8 × 30" F&B, T-type, Cd 7, Silicone	2.5"	1.5"	1.5"
810084G25ATT025L	8 × 40" F&B, T-type, Cd 7, Silicone	2.5"	1.5"	1.5"
810123G25ATT025L	12 × 30" F&B, T-type, Cd 7, Silicone	2.5"	1.5"	1.5"
810124G25ATT025L	12 × 40" F&B, T-type, Cd 7, Silicone	2.5"	1.5"	1.5"
810183G25ATT030L	18 × 30" F&B, T-type, Cd 7, Silicone	3"	1.5"	1.5"
810243G25ATT040L	24 × 30" F&B, T-type, Cd 7, Silicone	4"	1.5"	1.5"
810303G25ATT040L	30 × 30" F&B, T-type, Cd 7, Silicone	4"	1.5"	1.5"

Notes

1. Additional versions are available on request

Accessories

Order Code	Description
8ZML--0007	Diaphragm Pressure Gauge, 1.5" TC, 0-10 bar, Glycerin
8ZVL--0002	Ball Valve, 1.5" TC
8ZVL--0005	Sampling Valve, 1/2" (12 mm), 1.5" TC
8ZVL--0003	Butterfly Valve, 1" TC
8ZVL--0004	Butterfly Valve, 1.5" TC
8ZML--0009	Sight Glass, 1.5" TC
8ZML--0021	Press. Gauge, Sampling Valve, Adapter KIT, 1.5" TC (10 bar)

Notes

1. Additional options may be available

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
Argentina +54.11.4721.0505
Brazil +55.11.4362.8900
Mexico +52.55.5562.1102
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.913.586.098
Russian Federation +7.812.327.53.27
Japan +81.3.4331.4300
China +86.21.6878.2300

Specifications subject to change without notice.
Copyright Sartorius Stedim Biotech GmbH.
Printed in the EU on paper bleached without chlorine.
Publication No.: S--2094-e160903
Order No.: 85037-542-07
Ver. 09 | 2016