

F&B Filter Housings

For the Food and Beverage Industry



Introduction

Sartorius Stedim Biotech F&B Filter Housings are specifically designed for liquid filtration applications of the Food and Beverage Industry. Manufactured in an PED 97|23|EC certified facility, special attention has been paid to choice of materials, durability, cleanability, ease of use and quality control.

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 $32 \mu\text{inch}$ ($0.8 \mu\text{m}</math>) Ra mechanical polish.$

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 97|23|EC Standard

Sartorius Stainless Steel Housings are designed and manufactured according to the Pressure Equipment Directive PED 97|23|EC.
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 Specifications

Product Contact Surfaces	316L Surfaces
Gasket Materials	Silicone (Fluoroelastomer 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 <math><0.8 \mu\text{m}</math> internal MP Ra <math><1.6 \mu\text{m}</math> external MP

Standard Order Codes	Basic Description	In Out TC	Vent TC	Drain TC
810011G25ATT010L	1+10" F&B, T-type, Cd7, 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

Diaphragm Pressure Gauge, 1.5" TC, 0-10 bar, Glycerin	8ZML-0007
Ball Valve, 1.5" TC	8ZVL-0002
Sampling Valve, ½" (12 mm), 1.5" TC	8ZVL-0005
Butterfly Valve, 1" TC	8ZVL-0003
Butterfly Valve, 1.5" TC	8ZVL-0004
Sight Glass, 1.5" TC	8ZML-0009
Press. Gauge, Sampling Valve, Adapter KIT, 1.5" TC (10 bar)	8ZML-0021

Notes


1. Additional options may be available

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