



Single Rounds

Mini & 1 Element Filter Housings

Product Information

There has been, and is, an increasing demand for quality filter cartridge systems for sterilizing and polishing filtration processes. A large emphasis has been placed on the integrity of construction of the filter cartridges. However, the filter cartridge housing is just as an important part of any filtration system. Without a proper housing the cartridge is useless.



Introduction

Even the best cartridge cannot do the job if enclosed in a housing that allows fluid to bypass the filter, has external leaks, are not chemically or mechanically compatible with the application. Quality gas or liquid filtration systems require both quality housings and quality filter cartridges. To meet this need, Sartorius Stedim Biotech has been producing a sanitary line of housings with quality as the primary objective.

Quality of Materials

Only 316L Stainless steel is used for all wetted surfaces to provide maximum durability. Supplied O-rings and gaskets are compounded only from FDA approved materials that meet the requirements for direct contact with food and pharmaceutical products.

Quality Surface Finishes

All Sartorius Stedim Biotech Sanitary housings come standard with internal finishes of at least 0.5 micron Ra and are electropolished. Electropolishing removes surface impurities in stainless steel left over from the machining and the finishing processes. Such impurities are sites for future initiation of corrosion and possible sources of contaminants leaching into the product.

Electropolishing also smoothes the microscratches left by mechanical polishing, thus reducing the total surface area the product will contact, and making it harder for bacteria or contaminants to lodge on the housing surface.

Finally, electropolishing leaves a highly corrosion resistant, passive film on the surface of the steel (passivation). Thus electropolishing is the recommended finish for all applications where cleanliness and corrosion resistance are critical.

Ease in Cleaning

Sartorius Stedim Biotech utilizes a unique filter cup design that is conducive for allowing a thorough cleaning.

Flexibility

Sartorius Stedim Biotech offers the widest range of housing sizes and design options to exactly match your flow rate and pressure differential requirements. Connections are available in many styles and sizes. Custom designs and unique configurations are available upon request.

Quality Control and Documentation

An important feature of pharmaceutical process validation is documentation. All our housings are given stringent inspections during and after manufacturing including dimensional checks, weld inspections, surface measurements and hydrostatic testing. Each housing is labeled by laser with a matching serial number on the bell and base. This serial number provides complete tractability for the Quality Control Certificate, Material Test Reports, and Weld Logs.

Ease of Installation

Sartorius Stedim Biotech housings are sold ready to install with all gaskets, o-rings and clamps. All that is required are the components needed to connect to your existing hardware.

PED 2014/68/EU Standard

Sartorius Stedim Biotech Stainless Steel Housings are designed and manufactured according to 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 specifications

Materials	
All Wetted Surfaces	316L
Clamps	304
Seals	Silicone (Fluoroelastomer or EPDM optional)

Available Heights	
Mini	5"
Single Round	5", 10", 20", 30", 40"

Surface Finishes	
Interior	Ra <0.5 µm EP
Exterior	Ra <1.6 µm EP

Housing Ratings	
Pressure	-1 + 10 bar
Temperature	-10 + 150°C

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--2009-e170806
Order No.: 85030-522-29
Ver. 08 | 2017