

# Biopharma Pressure Vessel Type 380



## Product Information

Sartorius pharmaceutical-grade vessels are pressure-resistant tanks for the transportation, storage and distribution of liquids.

### Introduction

Our pressure vessels are manufactured in facilities certified for compliance with the European Pressure Equipment Directive 2014/68/EU. Thorough cleanability and ease of use are the major features of our pressure vessels. Sartorius pressure vessels are designed to fulfill our customers' quality standards and thus completely meet market expectations.

### Use

The vessel filled with the liquid to be filtered is drained by feeding in compressed air or gas. For this purpose, a pressure hose is connected to a pressure source and another pressure hose is connected to the outlet side of the tank and a filtration system (such as a filter holder or capsule). When pressure is applied, liquid flows into the outlet tube of the tank and through the connected pressure hose into the filtration system.

## Ease of Cleaning

The vessels are designed to ensure thorough cleaning. These pharmaceutical-grade vessels can be sterilized by autoclaving at 134 °C.

## Accessories (optional)

### Trolley

- 3ZGLA-0106 – Trolley for 5 L and 10 L
- 3ZGLA-0101 – Trolley for 20 L to 100 L

### TC-Gasket DN40 (41 mm × 1.5 mm TC 50.5 mm for Inlet)

- 7EDECV0170 - EPDM
- 7EDSCV0170 - Silicone
- 7EDVCV0170 - FKM (Fluoroelastomer)
- 7EDWCV0170 - PTFE with FKM-core (Fluoroelastomer)

### TC-Gasket DN25 (29 mm × 1.5 mm TC 50.5 mm for Outlet)

- 7EDECV0132 - EPDM
- 7EDSCV0132 - Silicone
- 7EDVCV0132 - FKM (Fluoroelastomer)
- 7EDWCV0132 - PTFE with FKM-core (Fluoroelastomer)

### TC-Gasket DN 6.625" (168.3 mm × 2.6 mm TC 183 mm for lid)

- 7EDECV0008 - EPDM
- 7EDSCV0008 - Silicone
- 7EDVCV0008 - FKM (Fluoroelastomer)
- 7EDWCV0008 - PTFE with FKM-core (Fluoroelastomer)

### Safety valve

- 3ZVIG-0134 – Safety valve spring loaded 4 bar DN25 TC 50.5 mm; Medium group I, type examined for gases, steam and fluids, 316L
- 3ZVIG-0133 – Safety valve spring loaded 3 bar DN25 TC 50.5 mm; Medium group I, type examined for gases, steam and fluids, 316L
- 3ZVIG-0132 – Safety valve spring loaded 2 bar DN25 TC 50.5 mm; Medium group I, type examined for gases, steam and fluids, 316L
- 3ZVIG-0131 – Safety valve spring loaded 1 bar DN25 TC 50.5 mm; Medium group I, type examined for gases, steam and fluids, 316L

### Manometer

- 7ZMA-0024 - Manometer 0–10 bar, TC 50.5 mm, autoclavable at 134 °C

### Diaphragm valve

- 7ZVD-0111 – Diaphragm valve DN25, 2 × TC 50.5 mm, material 316L, EPDM

## Quality of Materials

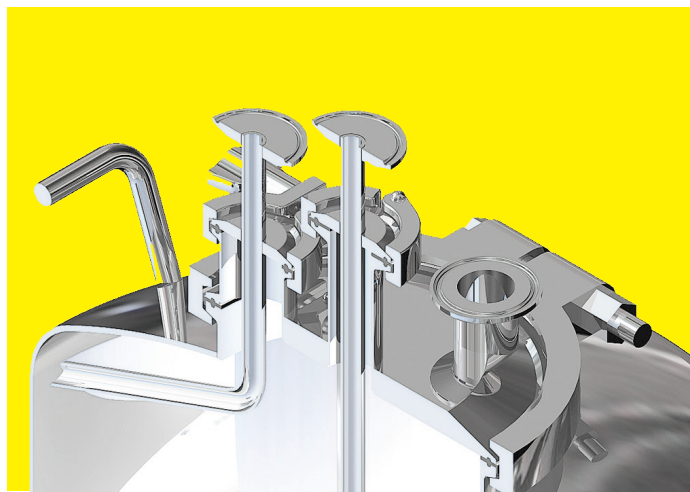
Only 316L stainless steel is used for all product-contact surfaces in order to provide maximum durability. The gaskets supplied with the vessel are FDA and USP Class VI approved materials that meet the requirements for direct contact with food and pharmaceutical products.

## Quality Control and Documentation

An important part of pharmaceutical process validation is documentation. All our vessels undergo stringent testing during and after manufacturing. Each vessel is labeled by laser with a matching serial number on the body. This serial number provides complete traceability to the Quality Control Certificate and Material Test Reports.

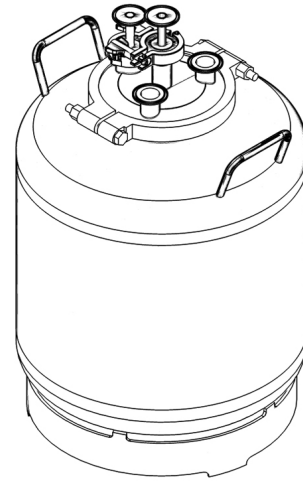
## European Pressure Equipment Directive

The design and manufacture of Sartorius Stedim Biotech pharmaceutical-grade pressure vessels meet the requirements of the European Pressure Equipment Directive 2014/68/EU. Our production facilities fulfill the highest quality standards and are audited and monitored at regular intervals by both our in-house quality control system and by independent, accredited auditors. The optional safety valve we use on our pressure vessels has been approved and certified according to AD2000 and is optimally matched for each specific vessel size.



# Technical Specifications

Product-contact surfaces	316L; others available on request
Gasket materials	Silicone, Fluoroelastomer, EPDM FEP, FFKM, etc.
CFR compliance	Gasket materials comply with the FDA Code 21 CFR 172.600 and USP Class VI
Surface finishes	Ra <0.5 µm interior Ra <1.6 µm exterior
Surface quality	Electropolished
Max. pressure	
5 L, 10 L, 20 L, 30 L, 40 L	4 bar
60 L, 80 L	2 bar
50 L	3 bar
100 L	1 bar



## Ordering Information

Order Number with Standard N-documentation	Volume (Net)	Operation pressure	Weight	Height vessel	Height lid	Diameter
380AA0005ID006LN	5 L	-1...+4 bar	9.0 kg	200 mm	170 mm	219.1 mm
380AA0010ID006LN	10 L	-1...+4 bar	11.0 kg	350 mm	170 mm	219.1 mm
380AA0020FD002LN	20 L	-1...+4 bar	19.0 kg	290 mm	100 mm	406.4 mm
380AA0030FD002LN	30 L	-1...+4 bar	21.4 kg	370 mm	100 mm	406.4 mm
380AA0040FD002LN	40 L	-1...+4 bar	23.8 kg	450 mm	100 mm	406.4 mm
380AA0050FD022LN	50 L	-1...+3 bar	26.3 kg	530 mm	100 mm	406.4 mm
380AA0060FD002LN	60 L	-1...+2 bar	28.7 kg	610 mm	100 mm	406.4 mm
380AA0080FD002LN	80 L	-1...+2 bar	33.5 kg	770 mm	100 mm	406.4 mm
380AA0100FD022LN	100 L	-1...+1 bar	41.4 kg	1030 mm	100 mm	406.4 mm

Order Number with Premium Z-documentation	Volume (Net)	Operation pressure	Weight	Height vessel	Height lid	Diameter
380AA0005ID006LZ	5 L	-1...+4 bar	9.0 kg	200 mm	170 mm	219.1 mm
380AA0010ID006LZ	10 L	-1...+4 bar	11.0 kg	350 mm	170 mm	219.1 mm
380AA0020FD002LZ	20 L	-1...+4 bar	19.0 kg	290 mm	100 mm	406.4 mm
380AA0030FD002LZ	30 L	-1...+4 bar	21.4 kg	370 mm	100 mm	406.4 mm
380AA0040FD002LZ	40 L	-1...+4 bar	23.8 kg	450 mm	100 mm	406.4 mm
380AA0050FD022LZ	50 L	-1...+3 bar	26.3 kg	530 mm	100 mm	406.4 mm
380AA0060FD002LZ	60 L	-1...+2 bar	28.7 kg	610 mm	100 mm	406.4 mm
380AA0080FD002LZ	80 L	-1...+2 bar	33.5 kg	770 mm	100 mm	406.4 mm
380AA0100FD022LZ	100 L	-1...+1 bar	41.4 kg	1030 mm	100 mm	406.4 mm


\* Standard design: Material: 316L; silicone gasket; electropolished internal and external surfaces; incl. fill and drain tube (13 x 1.5 mm d) with sanitary TC.

**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](http://www.sartorius.com)