



Pressure Vessels

Pressure vessels with Capacities
from 5–100 Liters

Product Information

Sartorius Stedim Biotech pressure vessels are pressure-resistant vessels for the transport, storage and distribution of liquids. Our pressure vessels are manufactured in facilities certified for compliance with the European Pressure Equipment Directive 2014/68/EU. Flexibility, thorough cleanability and ease of use are the major features of our pressure vessels.

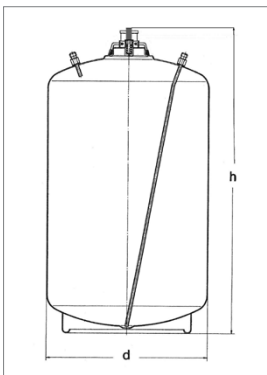


Use

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

Quality of Materials

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



Ease of Cleaning

The vessels are designed to ensure thorough cleaning. Vessels can be sterilized, along with the pressure gauge in place, by autoclaving at 121°C (30 min.). The maximum operating temperature is 95°C.

Easy Installation

Sartorius Stedim Biotech pressure vessels are supplied as ready-to-use units along with a preset safety valve, a pressure gauge and gaskets. A wide variety of accessories, such as tubing adapters, is available to choose from.

European Pressure Equipment Directive

The design and manufacture of Sartorius Stedim Biotech 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 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. The vessel and the safety valve are certified acc. to PED 2014/68/EU for media of fluid group 2.

Technical specifications

Product-contact surfaces	316 L
Gasket materials	Silicone (Standard)
CFR-Compliance	Gasket materials comply with the FDA regulations code 21 CFR 17.2600
Surface finishes	Ra < 0.8 µm internal Ra < 1.6 µm external
Surface quality	E-polished

Please note that the pressure vessels may not be operated under vakuum.



Safety valve TÜV.SV.2055 CE 0036 type examined
Ø DN 8 Brass nickel-plated

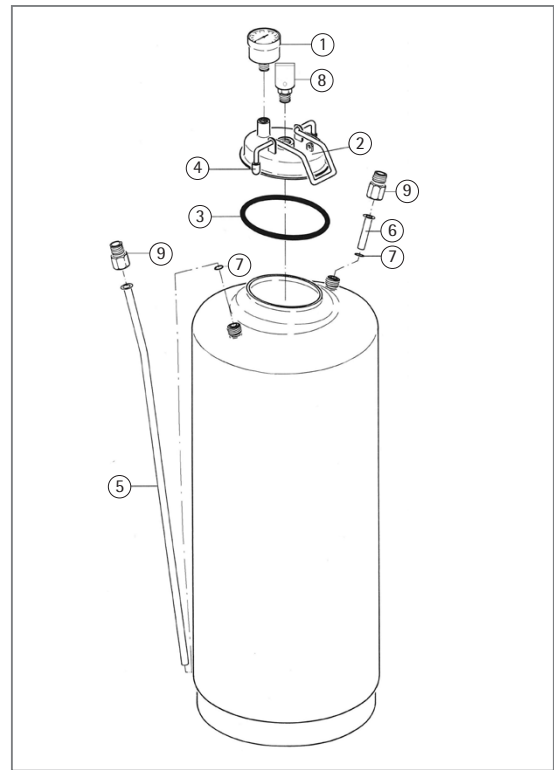
Ordering Information

Replacement parts and Accessories

Type Ordercode	17530	17531	17532	17533	17534	17535	17536
Volume	5 ltr.	10 ltr.	20 ltr.	40 ltr.	60 ltr.	80 ltr.	100 ltr.
Max. op. pressure	7 bar	7 bar	7 bar	5 bar	3 bar	2 bar	2 bar
Max. op. temperature	95°C	95°C	95°C	95°C	95°C	95°C	95°C
Weight	3.9 kg	5.35 kg	8.15 kg	11.8 kg	15.2 kg	18.4 kg	21.7 kg
Diameter Ø	234 mm	234 mm	234 mm	300 mm	400 mm	400 mm	400 mm
Height h ± 8 mm	265 mm	390 mm	630 mm	735 mm	673 mm	832 mm	992 mm

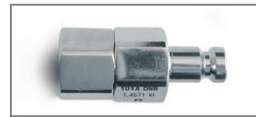
Replacement parts and Accessories

Pos.	Replacement part	Order no.																																	
1A	Pressure gauge 7 bar	6980390																																	
1B	Pressure gauge 5 bar	6980415																																	
1C	Pressure gauge 2 bar	6986119																																	
1D	Pressure gauge 3 bar	6986137																																	
2A	Lid with valve and pressure gauge 7 bar	6986130																																	
2B	Lid with valve and pressure gauge 5 bar	6986129																																	
2C	Lid with valve and pressure gauge 2 bar	6986131																																	
2D	Lid with valve and pressure gauge 3 bar	6986138																																	
3A	Silicone o-ring for lid acc. FDA	6986110																																	
3B	EPDM o-ring for lid acc. FDA	6986111																																	
3C	Fluoroelastomer o-ring for lid acc. FDA	6980389																																	
4	PTFE cap for lid clamp (pack of 2)	6985131																																	
5A	Outlet tube 5 ltr.	6986112																																	
5B	Outlet tube 10 ltr.	6986113																																	
5C	Outlet tube 20 ltr.	6986114																																	
5D	Outlet tube 40 ltr.	6986115																																	
5E	Outlet tube 60 ltr.	6986116 </tr <tr> <td>5F</td> <td>Outlet tube 80 ltr.</td> <td>6986117</td> </tr> <tr> <td>5G</td> <td>Outlet tube 100 ltr.</td> <td>6986118</td> </tr> <tr> <td>6</td> <td>Inlet tube</td> <td>6980395</td> </tr> <tr> <td>7A</td> <td>O-ring for in- outlet tubes Silicone</td> <td>6986132</td> </tr> <tr> <td>7B</td> <td>O-ring for in- outlet tubes EPDM</td> <td>6986133</td> </tr> <tr> <td>7C</td> <td>O-ring for in- outlet tubes Fluoroelastomer</td> <td>6980396</td> </tr> <tr> <td>8A</td> <td>Safety valve 1/4" 2 bar acc. PED brass nickel-plated</td> <td>6986701</td> </tr> <tr> <td>8B</td> <td>Safety valve 1/4" 5 bar acc. PED brass nickel-plated</td> <td>6986702</td> </tr> <tr> <td>8C</td> <td>Safety valve 1/4" 7 bar acc. PED brass nickel-plated</td> <td>6986703</td> </tr> <tr> <td>8D</td> <td>Safety valve 1/4" 3 bar acct. to 97/23/EC nickel-plated</td> <td>6986704</td> </tr> <tr> <td>9</td> <td>Connector 19/32 18-NS-2 female. thread G 3/8" male thread</td> <td>6980420</td> </tr>	5F	Outlet tube 80 ltr.	6986117	5G	Outlet tube 100 ltr.	6986118	6	Inlet tube	6980395	7A	O-ring for in- outlet tubes Silicone	6986132	7B	O-ring for in- outlet tubes EPDM	6986133	7C	O-ring for in- outlet tubes Fluoroelastomer	6980396	8A	Safety valve 1/4" 2 bar acc. PED brass nickel-plated	6986701	8B	Safety valve 1/4" 5 bar acc. PED brass nickel-plated	6986702	8C	Safety valve 1/4" 7 bar acc. PED brass nickel-plated	6986703	8D	Safety valve 1/4" 3 bar acct. to 97/23/EC nickel-plated	6986704	9	Connector 19/32 18-NS-2 female. thread G 3/8" male thread	6980420
5F	Outlet tube 80 ltr.	6986117																																	
5G	Outlet tube 100 ltr.	6986118																																	
6	Inlet tube	6980395																																	
7A	O-ring for in- outlet tubes Silicone	6986132																																	
7B	O-ring for in- outlet tubes EPDM	6986133																																	
7C	O-ring for in- outlet tubes Fluoroelastomer	6980396																																	
8A	Safety valve 1/4" 2 bar acc. PED brass nickel-plated	6986701																																	
8B	Safety valve 1/4" 5 bar acc. PED brass nickel-plated	6986702																																	
8C	Safety valve 1/4" 7 bar acc. PED brass nickel-plated	6986703																																	
8D	Safety valve 1/4" 3 bar acct. to 97/23/EC nickel-plated	6986704																																	
9	Connector 19/32 18-NS-2 female. thread G 3/8" male thread	6980420																																	

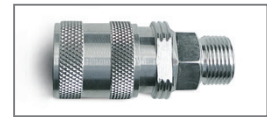


Accessories for pressure filtration (optional)

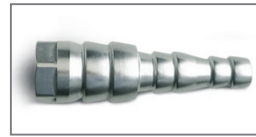
Pos.	Article	Order no.
1	PTFE pressure hose 1.5 m stainless steel armoured G 3/8"	16999
2	PTFE pressure hose 0.8 m stainless steel armoured G 3/8"	16823
3	Sanitary flange 1"-1½" TC to 18-NS2 female thread	17070
4	Hose nipple 18-NS2 female thread for tubes DN10, DN 13 und DN19	16863
5	Sanitary flange 18-NS2 female thread to quick connect nipple	17170
6	Quick connect nipple to G 3/8" male thread	6985128
7	Sanitary flange 1"-1½" TC to quick connect nipple	17154
8	Sanitary flange 1"-1½" TC to quick connector nipple Stäubli RBE03	7ZML-0012
9	Sanitary flange 1"-1½" TC to G 3/8" male	17019
10	Pressure pump 220 Volt Pmax. 7 bar, R 3/8"	16617
11	Compressed air pistol with quick connector nipple	6980407
12	Trolley for 40, 60, 80, und 100 ltr. container	17640
13	Clamp closure 1"-1½" (50.5 mm) for sanitary flange	7ZSB-0009
14	Gasket silicone for sanitary flange 1"-1½"	7EDSCV0003



17170



6985128



16863



17070



17019



16999



17640

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--2067-e161005
 Order No.: 85032-537-76
 Ver. 10 | 2016