

Biopharma Pressure Vessel Type 385

The Simple Solution

Product Information

Sartorius pharmaceutical-grade vessels are pressure-resistant tanks for transporting, storing, and distributing liquids.

Features and Benefits

- Scalable volumes (0.1 L to 200 L)
- Simple design – simple usage
- Suitable for dangerous media
- Optional trolleys



Introduction

Our vessels are manufactured in facilities that are certified for compliance with the European Pressure Equipment Directive 2014/68/EU. The key features of our pressure vessels are their ease-of-use and thorough cleanability. Sartorius pressure vessels are designed to fulfill our customers' quality standards and meet their needs.

Use

Once filled with the liquid to be filtered, the vessel can be drained by feeding in compressed air or gas. To do this, connect a pressure hose to a pressure source and another pressure hose the outlet side of the tank and a filtration system (e.g. 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.

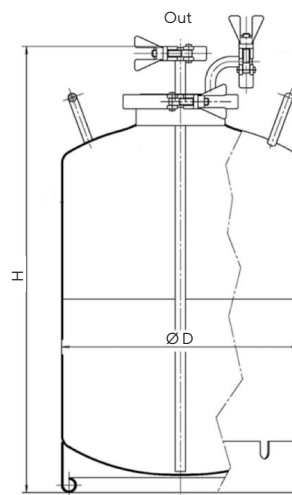
Technical Specifications

Vessel

Material	Vessels: 316 L (1.4404) Trolleys: 1.4301
Surface:	- Inside and outside electro-polished - Ra-inside ≤ 0.8 µm
Gasket material	EPDM
Documentation	Manual Drawing Declaration of conformity Manufacturer certificate Certificate of marking control and traceability material and gasket certificates
Autoclaving Conditions	+121 °C at 3.6 bar
Fluid Group acc. to PED 2014-68-EU:	1 (dangerous)
Inlet Outlet tube:	½" - 12.7 mm × 1.65 mm
Inlet Outlet TC (0.1 L and 0.25 L)	TC 25
Inlet Outlet TC (1 L - 200 L)	TC 50.5

Ordering Information

Order Number	Net Volume [L]	Operation Pressure [bar]	Operating Temperature [°C]	Weight [kg]	Height (H) [mm]	Outer Diameter (D) [mm]	Main TC Diameter (Lid) [mm]
385AA0001LM001IC	0.1	-1 +5	+50	1.5	120	67	64
385AA0002LM001IC	0.25	-1 +5	+50	2	200	67	64
385AA0005LM001IC	0.5	-1 +5	+50	2.5	230	89	91
385AA0010LM001IC	1	-1 +5	+50	3	250	108	106
385AA0020LM001IC	2	-1 +5	+50	3	325	140	119
385AA0050LM001IC	5	-1 +5	+50	3,5	380	181	119
385AA0100LM001IC	10	-1 +5	+50	6	430	229	119
385AA0200LM001IC	20	-1 +5	+50	8	500	278	119
385AA0300LM001IC	30	-1 +5	+50	13	620	324	119
385AA0500LM001IC	50	-1 +5	+50	16	700	363	119
385AA1000LM001IC	100	-1 +5	+50	32	900	450	119
385AA2000LM001IC	200	-1 +4	+50	45	1,000	600	119



Spare Parts and Accessories

Clamps

Vessel Size	Vessel Main Clamp	Order Code
Volume	Clamp Size	
0.1 - 0.25	2" - TC 64	7ZSB--0202
0.5	3" - TC 91	7ZSB--0203
1	3.5" - TC 106	7ZSB--0204
2 - 200	4" - TC 119	7ZSB--0205



Vessel Size	Inlet Outlet Clamp	Order Code
Volume	Clamp Size	
0.1 - 0.25	TC 25	7ZSB--0200
0.5	TC 25	7ZSB--0200
1	TC 50.5	7ZSB--0201
2 - 200	TC 50.5	7ZSB--0201



Gaskets

Gasket Material	Gasket Size [in]	TC-Diameter [mm]	Order Code
EPDM	4" - main gasket	119	7EDECV0007
Silicone	4" - main gasket	119	7EDSCV0007
EPDM	3.5" - main gasket	106	7EDECV0183
Silicone	3.5" - main gasket	106	7EDSCV0183
EPDM	3" - main gasket	91	7EDECV0006
Silicone	3" - main gasket	91	7EDSCV0006
EPDM	2" - main gasket	64	7EDECV0004
Silicone	2" - main gasket	64	7EDSCV0004
EPDM	1.5" - in out	50.5	7EDECV0003
Silicone	1.5" - in out	50.5	7EDSCV0003
EPDM	½" - in out	25	7EDECV0001
Silicone	½" - in out	25	7EDSCV0001



Trolleys With a Swivel Castor and Fixed Castor (PA-Conductive)

Vessel Net Volume [L]	Outer Diameter Vessel [mm]	Size (L × W × H) [mm]	Weight [kg]	Order Code (Trolley)
10	229	240 × 240 × 150	5	1ZNI--0001
20	278	280 × 290 × 150	5.3	1ZNI--0002
30	324	330 × 330 × 150	5.7	1ZNI--0003
50	363	380 × 380 × 150	6.1	1ZNI--0004
100	450	460 × 460 × 150	7.1	1ZNI--0005
200	600	610 × 610 × 150	8.8	1ZNI--0006



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 more information, visit

www.sartorius.com