

# Palletank<sup>®</sup> for In-Process Fluid Handling

## Benefits

- Easy operation due to continuous processing without movement of the containers
- Easy access through double hinged front door with safety latches
- Optimal bag unfolding and filling through a lifting system for large size containers

## Product Information

The Palletank<sup>®</sup> for In-Process Fluid Handling are specifically developed for users who wish to leave the containers in close proximity to process equipment. Palletank<sup>®</sup> for in-process fluid handling are available in 200 L, 500 L, 1000 L, 1500 L, 2000 L, 2500 L and 3000 L volumes.<sup>1</sup> Flexsafe<sup>®</sup> I Flexel<sup>®</sup> 3D bags are manufactured according to a qualified design that precisely fits the Palletank<sup>®</sup>.

<sup>1</sup>1500 L volume and above are only available with Flexel<sup>®</sup> 3D bags.



## Easy Operation

The double hinged front doors allow bag replacement, installation and manipulation whilst Palletank® are stacked up in the process area (up to 1000 L), thus enabling continuous processing without movement of the containers. As a result, a forklift is not routinely required in the manufacturing area. A specially designed dolly, up to 1000 L (accessory), facilitates the convenient movement of material throughout the facility.

## Ready for Large Volumes

The unique lifting system for large volume tanks provides optimal bag unfolding and filling. A specially designed bottom gate allows safe passage of large bore tubing, 1.5" tri clamps, QC bags and filters. The front and side bin access also eases the bag deployment in the container when filling.

## Palletank® Family

Besides the Palletank® for In-Process Fluid Handling, the product range of Palletank® containers includes the following lines, specifically developed for the various application requirements of fluid management in the biopharmaceutical industry:

- Palletank® for Shipping
- Plastic and stainless steel Palletank® for Storage
- Palletank® for Lev Mixer | Magnetic Mixer

## Technology Integration Support

Sartorius Stedim Biotech supports users from the design and implementation phase of a new production facility with the most comprehensive support program, that ensures successful design implementation and validation of single-use manufacturing.

## Security of Supply

Sartorius Stedim Biotech has established multiple manufacturing sites with consistent industrial processes. The expertise of designing single-use solutions based on collaborative supplier management and customer demand planning ensures a state-of-the-art and robust supply chain that can cope with strong market growth.

## Quality Assurance

Palletank® for In-Process Fluid Handling are designed, developed and manufactured in accordance with a ISO 9001 certified Quality Management System. They undergo extensive testing before shipping.

# Technical Data

## Specifications

### Palletank® for In-Process Fluid Handling

Description	Palletank® 200 L for In-Process Fluid Handling	Palletank® 500 L for In-Process Fluid Handling	Palletank® 1000 L for In-Process Fluid Handling
Bag volume(s)	100 L or 200 L	500 L	1000 L
Construction material	304 L stainless steel		
Surface finishing	Glass bead blasted		
Dimensions (W×D×H)	860 × 660 × 929 mm (33.9 × 26 × 36.6 in)	1260 × 860 × 989 mm (49.6 × 33.9 × 38.9 in)	1260 × 1060 × 1443.5 mm (49.6 × 41.7 × 59.3 in)
Weight (approx.)	50 kg (110.2 lb)	89 kg (196.2 lb)	120 kg (264.6 lb)
Bottom gate	1	1	2
Stackability	3 high	3 high	2 high
Total weight (approx.)	68 kg (150.2 lb)	111 kg (244.7 lb)	144.5 kg (317.9 lb)

Description	Palletank® 1500 L for In-Process Fluid Handling	Palletank® 2000 L for In-Process Fluid Handling	Palletank® 2500 L for In-Process Fluid Handling	Palletank® 3000 L for In-Process Fluid Handling
Bag volume(s)	1500 L	2000 L	2500 L	3000 L
Construction material	304 L stainless steel			
Surface finishing	Glass bead blasted			
Dimensions (W×D×H)	1200 × 1000 × 2043 mm (47.2 × 39.4 × 80.4 in)	1200 × 1000 × 2543 mm (47.2 × 39.4 × 100.1 in)	1200 × 1000 × 3043 mm (47.2 × 39.4 × 119.8 in)	1200 × 1000 × 3543 mm (47.2 × 39.4 × 139.5 in)
Weight (approx.)	300 kg (661.4 lb)	350 kg (771.6 lb)	400 kg (881.8 lb)	450 kg (992 lb)
Bottom gate	3			
Tubing access in front of Palletank®	3			
Stackability	No			

# Specifications

Palletank® for In-Process Fluid Handling

Description	Dolly for Palletank® 200 L for In-Process Fluid Handling	Dolly for Palletank® 500 L for In-Process Fluid Handling	Palletank® 1000 L for In-Process Fluid Handling
Weight	18 kg (40 lb)	22 kg (48.5 lb)	24.5 kg (54 lb)

## Ordering Information


Order Code	Description
FXC106230	Palletank® 200 L for In-Process Fluid Handling
FXC106235	Palletank® 500 L for In-Process Fluid Handling
FXC106223	Palletank® 1000 L for In-Process Fluid Handling
FXC109797	Palletank® 1500 L for In-Process Fluid Handling
FXC109798	Palletank® 2000 L for In-Process Fluid Handling
FXC109799	Palletank® 2500 L for In-Process Fluid Handling
FXC109764	Palletank® 3000 L for In-Process Fluid Handling
FXS102255	Dolly for Palletank® 200 L for In-Process Fluid Handling
FXS102257	Dolly for Palletank® 500 L for In-Process Fluid Handling
FXS102259	Dolly for Palletank® 1000 L for In-Process Fluid Handling

### 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)

Specifications subject to change without notice.

© 2021 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France