

# Palletank<sup>®</sup> for Shipping

## Benefits

- ASTM qualification of Flexsafe<sup>®</sup> 3D bags with the Palletank<sup>®</sup> for Shipping ensures safe and robust liquid shipping applications
- Immobilisation of Flexsafe<sup>®</sup> I Flexel<sup>®</sup> 3D bag by compression ensures system integrity
- Dismountable construction improves space utilization on site before and after use

## Product Information

The Palletank<sup>®</sup> for Shipping is a qualified system that is specifically developed for safe and robust shipment of sterile bulk pharmaceutical fluids. It is available in volumes of 100 L, 200 L and 500 L. The system is composed of two basic elements, the Palletank<sup>®</sup> container and the associated shipping kit. The shipping kit acts as a compression unit to firmly secure the filled Flexsafe<sup>®</sup> I Flexel<sup>®</sup> 3D bag whilst in transit. Top and bottom tubing is protected to prevent damage and a rigid lid is locked securely before shipping. The Flexsafe<sup>®</sup> I Flexel<sup>®</sup> 3D bags are manufactured according to a qualified design that precisely fits the Palletank<sup>®</sup> to ensure highest safety during shipping.



## Safe Transportation of Critical Drug Substances

The superior strength and flexibility of Flexsafe® bags make them safe for critical liquid shipping applications by truck and air. The robustness of Flexsafe® bags for shipping is proven by applying extensive qualification regimes that combine the most stringent standard ASTM D4169 testing and real shipping conditions.

## Space-Saving

The Palletank® for Shipping can be stacked, when not in transit, up to three high to maximize the utilization of the available clean-room area. When not in use or on return, the Palletank® is foldable and stackable up to five high.

## Easy Cleanability

In order to simplify the cleaning | disinfection cycles, all movable parts of the Palletank® for Shipping can be dismantled (including door hinges).

## Palletank Family

The Palletank® for Shipping is part of the family of Palletank® containers which includes the following lines specifically developed for the various application requirements on fluid management in the biopharmaceutical industry:

- Palletank® for In-Process Fluid Handling
- Palletank® for Storage
- Plastic Palletank® for Storage
- Palletank® for Lev Mixer | Magnetic Mixer

## 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 Shipping are designed, developed and manufactured in accordance with a ISO 9001 certified Quality Management System. They undergo extensive testing before shipping.

# Technical Data

## Specifications

Description	Palletank® 100 L for Shipping	Palletank® 200 L for Shipping	Palletank® 500 L for Shipping
Bag volume(s)	100 L	200 L	500 L
Construction material	304 L		

## Dimensions (W × D × H)

Unfolded	800 × 600 × 688 mm (31.5 × 23.6 × 27.1 in)	800 × 600 × 935 mm (31.5 × 23.6 × 36.8 in)	1206 × 806 × 1050 mm (47.5 × 31.7 × 41.33 in)
Folded	800 × 600 × 510 mm (31.5 × 23.6 × 20.1 in)	800 × 600 × 510 mm (31.5 × 23.6 × 20.1 in)	1206 × 806 × 540 mm (47.5 × 31.7 × 21.26 in)
Weight (incl. lid)	63.5 kg (140 lb)	84 kg (185.18 lb)	129 kg (284.4 lb)
Bottom gates	1	1	1

## Stackability

Unfolded	3	3	2
Folded	5	5	3

## Accessories

Description	Shipping Kit 100 L	Shipping Kit 200 L	Shipping Kit 500 L
Weight	16 kg (35.27 lb)	16 kg (35.27 lb)	28 kg (61.8 lb)

## Ordering Information


Order Code	Description
FXC106362	Palletank® 100 L for Shipping
FXC103133	Palletank® 200 L for Shipping
FXC106609	Palletank® 500 L for Shipping
FXC102475	Palletank® 100 L shipping kit
FXC102477	Palletank® 200 L shipping kit
FXC102478	Palletank® 500 L shipping kit
FXS102254	Dolly for Palletank® 100 L   200 L for Shipping
FXS102256	Dolly for Palletank® 500 L for Shipping

**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

Publication No.: SPT2009-e | Order No.: 85034-534-88 | Status: 07 | 02 | 2021