



## Palletank for shipping

### Benefits

- ASTM qualification of Flexsafe® 3D Bags with the Palletank for shipping ensures safe and robust liquid shipping applications.
- Immobilisation of Flexsafe® | Flexel® 3D Bag by compression assures system integrity
- Dismountable construction improves space utilization on site before and after usage



### Product Information

The Palletank 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 container and the associated shipping kit. The shipping kit acts as a compression unit to firmly secure the filled Flexsafe® | Flexel® 3D Bag whilst in transit. Top and bottom tubing is protected to prevent damage and a rigid lid is locked securely prior to shipping. The Flexsafe® | Flexel® 3D Bags are manufactured according to a qualified design that precisely fits the Palletank to assure 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 3 high to maximize the utilization of the available clean room area. When not in use or upon return, the Palletank is foldable and stackable up-to 5 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 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 LevMixer® | 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 assures 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

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen, Germany  
Phone +49.551.308.0  
[www.sartorius-stedim.com](http://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. Printed and copyrighted by Sartorius Stedim Biotech GmbH.  
Publication No.: SPT2009-e180205  
Order No.: 85034-534-88  
Ver. 02 | 2018