

# Plastic Palletank<sup>®</sup> for Storage



## Benefits

### PP and ABS containers

Low cost and robust storage solution

### Stackable when full and collapsible when empty

Minimize storage volume

### Smooth scratch-resistant plastic surface

Clean-room compatible

### Standard product

Available from stock

## Product Information

The plastic Palletank<sup>®</sup> for Storage represents a cost-effective containment solution for the storage and processing of biopharmaceutical fluids contained in Flexsafe<sup>®</sup> | Flexel<sup>®</sup> 3D bags. Their flat, smooth surfaces and rounded corners facilitate cleaning and sanitization for clean-room operations. The plastic Palletank<sup>®</sup> for Storage are foldable for minimal storage space when empty. They are also stackable for optimal floor space utilization when filled with biopharmaceutical solutions. They are available as standard products with volumes of 200 L and 500 L that perfectly fit with 100 L | 200 L and 500 L Flexsafe<sup>®</sup> | Flexel<sup>®</sup> 3D bags.

## Applications

The plastic Palletank® for Storage with Flexsafe® | Flexel® 3D bags have been designed for the safe storage and processing of a wide range of biopharmaceutical fluids in a variety of applications such as:

- Buffers and media storage
- Bulk harvest
- Product pooling
- Fraction collection

## Optimal Space Utilization

The plastic Palletank® for Storage allows for secure stacking, up to three high, to meet the complex in-process and storage requirements, while maximizing the utilization of the available clean-room area. It saves up to 50% of the space required for cylindrical drums. When not in use, the collapsible tank can be folded to approximately half of its height for intermediate storage.

| Description |  |   |              |
|-------------|--|---|--------------|
| Volume      | Floor Space Required   |   | Space Saving |
|             | Plastic Palletank® for Storage                               | Drum                                      |              |
| 200 L       | 9 × 200 L stacked three high 1.5 m <sup>2</sup> (16.1 sq ft) | 9 × 200 L 3.2 m <sup>2</sup> (34.4 sq ft) | 50 %         |
| 500 L       | 6 × 500 L stacked two high 3 m <sup>2</sup> (32.3 sq ft)     | 6 × 500 L 3.7 m <sup>2</sup> (40 sq ft)   | 20 %         |

## Easy Setting and Use

One person can easily prepare a plastic Palletank® for Storage; the walls are raised and locked in place by hand. Their low tare weight ensures easier operator manipulation and facilitates intra-facility transfer.

## Hygienic Design

The plastic Palletank® for Storage is not subject to rust and corrosion due to all plastic construction. The containers are made from a robust, scratch-resistant plastic material. The flat, smooth walls and the rounded corners facilitate cleaning operations. The plastic Palletank® for Storage is compatible with cleaning and decontaminating agents commonly used in the biopharmaceutical industry.

# Technical Data

## Specifications

### 1. Plastic Palletank® for Storage

| Description                    |                   | STD Plastic Palletank® for Storage     |  |
|--------------------------------|-------------------|--|--|
|                                |                   | 200 L                                  | 500 L                                    |
| Part number                    |                   | FXC116215                              | FXC116216                                |
| Bag volume(s)                  |                   | 100 L   200 L                          | 500 L <sup>1</sup>                       |
| Main construction material     |                   | PP                                     | ABS                                      |
| Dimensions (W × D × H) Approx. | Unfolded          | 805 × 607 × 969 mm<br>32 × 24 × 38 in. | 1215 × 805 × 1086 mm<br>48 × 31 × 43 in. |
|                                | Folded            | 805 × 607 × 510 mm<br>32 × 24 × 20 in. | 1215 × 805 × 572 mm<br>48 × 31 × 22 in.  |
| Bottom gate                    |                   | 1                                      | 1  |
| Stackability -static           | Folded            | 5                                      | 3  |
|                                | Unfolded and full | 3                                      | 2  |

### 2. Ancillary Products

#### 2.1. Dolly

| Description            | STD Palletank® Accessory Dolly        |  |
|------------------------|---------------------------------------|--|
|                        | 200 L                                 | 500 L                                  |
| Part number            | FXA116630                             | FXA116631                              |
| Construction material  | 304 L and polyamide (wheels)          |  |
| Surface finishing      | Glass bead blasted                    |  |
| Dimensions (W × D × H) | 807 × 607 × 214 mm<br>32 × 24 × 8 in. | 1207 × 807 × 214 mm<br>47 × 32 × 8 in. |
| Weight (approx.)       | 10 kg (22 lb)                         | 20 kg (44 lb)                          |



<sup>1</sup>The maximum filling volume of the Flexel® 3D bag in the 500 L plastic Palletank® for Storage is 490 L.

## 2.2. Weighing Platform

| Description                              |  |  |
|--|--|--|
|  | IFS4-300LL-I                           | IFS4-600RN-I                           |
| Weighing capacity                        | 300 kg   1322 lb                       | 600 kg   1322 lb                       |
| Platform size (D + W)                    | 1000 × 1000 mm<br>39 × 39 in.          | 1500 × 1250 mm<br>59 × 49 in.          |
| Height                                   | Standard:<br>35 mm   1.4 in.           | Standard:<br>45 mm   1.8 in.           |
| Load plate                               | AISI304   1.4301 (V2A)<br>bead blasted | AISI304   1.4301 (V2A)<br>bead blasted |
| Resolution                               | 30.000 d                               | 30.000 d                               |
| Readability                              | 10 g                                   | 20 g                                   |
| Suitable with plastic Palletank® storage | 200 L                                  | 500 L                                  |

Please contact your sales representative for detailed ordering information.



## Ordering Information


| Order Code | Description   |
|------------|---|
| FXC116215  | STD Plastic Palletank® Storage 200 L                      |
| FXC116216  | STD Plastic Palletank® Storage 500 L                      |
| FXA116630  | STD Palletank® Accessory Dolly 200 L<br>(Plastic Storage) |
| FXA116631  | STD Palletank® Accessory Dolly 500 L<br>(Plastic Storage) |

**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.: SPT2022-e | Order No.: 85032-538-70 | Status: 07 | 02 | 2021