

Palletank[®] for Storage



Benefits

Qualified System

Perfect fit and protection of the Flexsafe[®] | Flexel[®] 3D bag in its Palletank[®]

Stackable Version

Space saving containment system

Technology Integration Support

For a successful single-use manufacturing implementation and validation.

Product Information

The Palletank[®] for Storage are stainless steel containers designed for the safe and robust storage of biopharmaceutical fluids contained in Flexsafe[®] | Flexel[®] 3D bags. They are available in volumes of 50 L, 200 L and 500 L. The Flexsafe[®] | Flexel[®] 3D bags are manufactured according to a qualified design that precisely fits the Palletank[®].

Applications

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

- Buffers and media storage
- Bulk harvest
- Product pooling
- Fraction collection
- Sample collection
- Bulk intermediate hold

Space-Saving

The stackable version of the Palletank® for Storage enables users to meet the complex in-process demands as well as the high requirements for storage, while maximizing the utilization of the available clean-room area. It saves up to 50% of the space required for cylindrical drums.

Safety

Flexsafe® | Flexel® 3D bags coupled with the rigid structure of Palletank® provide a stable and secure solution for processing, storage and in-house transportation of buffers, media, intermediates and final bulk products.

Palletank® Family

Besides the Palletank® for Storage, the product range of Palletank® container includes the following lines specifically developed for the various application requirements on fluid management in the biopharmaceutical industry:

- Palletank® for Shipping
- Palletank® for In-Process Fluid Handling
- Plastic Palletank® for Storage
- Palletank® for Mixing

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

The Palletank® for Storage are designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System. They undergo extensive testing before shipping.

Technical Data

Specifications

1. Palletank® for Storage

Description	Palletank® 200 L for Storage	Palletank® 500 L for Storage
Bag volume(s)	100 L or 200 L	500 L
Construction material	304 L stainless steel	304 L stainless steel
Surface finishing	Bead blasted	Bead blasted
Dimensions (W × D × H)	792 × 592 × 892 mm (31 × 23 × 35 in)	1192 × 792 × 1008 mm (46 × 31 × 39 in)
Weight	48 kg (106 lb)	100 kg (220 lb)
Bottom gate	1	1
Stackability	No	No

2. Palletank® for Storage Stackable

Description	Palletank® 50 L for Storage Stackable ¹	Palletank® 200 L for Storage Stackable	Palletank® 500 L for Storage Stackable
Bag volume(s)	50 L	100 L or 200 L	500 L
Construction material	304 L stainless steel	304 L stainless steel	304 L stainless steel
Surface finishing	Bead blasted	Bead blasted	Bead blasted
Dimensions (W × D × H)	490 × 490 × 750 mm (19 × 19 × 29 in)	792 × 592 × 929 mm (31 × 23 × 36 in)	1135 × 792 × 1044 mm (44 × 31 × 41 in)
Weight	25 kg (55 lb)	70 kg (154.3 lb)	110 kg (242 lb)
Bottom gate sliding gate	1	1	1
Stackability (Static)	3 high	3 high	2 high

¹ The 50L volume only available with Flexel® 3D bag

3. Ancillary Products

3.1. Dolly

Description	Dolly for Palletank® 50 L for Storage and Shipping	Dolly for Palletank® 200 L for Storage and Shipping	Dolly for Palletank® 500 L for Storage and Shipping
Construction material	304L and polyamide (wheels)	304L and polyamide (wheels)	304L and polyamide (wheels)
Surface finishing	Glass bead blasted	Glass bead blasted	Glass bead blasted
Dimensions (W × D × H)	490 × 490 × 185 mm (19.3 × 19.3 × 7.3 in)	815 × 615 × 188 mm (32.1 × 24.2 × 6.4 in)	1215 × 815 × 161 mm (47.8 × 32.1 × 6.4 in)
Weight (approx.)	7 kg (15.4 lb)	10 kg (22 lb)	18 kg (39.7 lb)



3.2. Weighing Platforms

The IFS4 flat-bed scales are entirely constructed of stainless steel and have an extremely low height, making it ideally suited for floor installation without a pit or anchoring.

The ramp is securely attached to the scale using special retainers for prevention of force shunt. This high-quality platform can be connected to any of a wide range of indicators, for use as a Class III legal measuring instrument or without legal verification. The CIS1 Combics 1 indicator allows strain gauge weighing with flat bed scales as well as with load cells to be connected.

	IFS4-150GG-I	IFS4-300LI-I	IFS4-1000RN-I
Weighing capacity	150 kg (330.7 lb)	300 kg (661.4 lb)	1000 kg (2204.6 lb)
Platform size	600 × 600 mm (23.6 × 23.6 in)	1000 × 800 mm (39.3 × 31.5 in)	1500 × 1250 mm (59 × 49.2 in)
Height	Standard: 35 mm	Standard: 35 mm	Standard: 45 mm
Load plate	AISI304 1.4301 (V2A) bead-blasted	AISI304 1.4301 (V2A) bead-blasted	AISI304 1.4301 (V2A) bead-blasted
Resolution	30.000 d	30.000 d	30.000 d
Readability	5 g	10 g	50 g
Suitable with Palletank® Storage and Storage Stackable	50 L	200 L	500 L

Refer to specific Sartorius datasheet for Combics indicators ranges, printers and other accessories specifications and ordering information.



Integrated Features

Features	Benefits	Palletank® for Storage	Palletank® for Storage Stackable
Level marks	Allow rapid visual monitoring of the fluid level in the bag	■	■ *
Integrated pallet	Allows easy carriage by pallet-jack or forklift	■	■
Tubing and fitting tray	Simplifies fluid handling operations and provides a convenient and secure place for inlet and outlet tubing assemblies during transport	■	■
Lid	Protects the bag against dust and light		■ *
Bottom gate sliding gate	Allows passage of large bore tubing, 1,5" tri clamps, QC bags and filters; facilitates bag positioning and maintain in position	■	■
Stacking corner	Enables the stacking of the Palletank® in order to maximize the utilization of available clean-room area		■
Dolly (accessory)	Facilitates the movement of material throughout a facility	■	■

* Except 50 L volume

Ordering Information


Order Code	Description
FXC110888	Palletank® 200 L for Storage
FXC110889	Palletank® 500 L for Storage
FXC113946	Palletank® 50 L for Storage Stackable
FXC110733	Palletank® 200 L for Storage Stackable
FXC110734	Palletank® 500 L for Storage Stackable
FXA113988	Dolly for Palletank® 50 L for Storage
FXS102254	Dolly for Palletank® 100 L 200 L for Storage and Shipping
FXS102256	Dolly for Palletank® 500 L for Storage and 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

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.: SPL2002-e | Order No.: 85030-533-48 | Status: 07 | 02 | 2021