

# Flexboy® Tray and Rack System



## Description

The Flexboy® Tray and Rack systems are designed to facilitate handling of both individual and manifold Flexboy® single-use bioprocessing bags (5 L-20 L) within biopharmaceutical manufacturing processes.

## Applications

The Flexboy® Tray and Rack systems are designed to support Flexboy® bags for the safe processing and transfer of a wide range of biopharmaceutical fluids in applications such as:

- Buffer and media storage
- Bulk harvest
- Product pooling
- Fraction collection
- Sample collection
- Bulk intermediate hold
- Final product transport

## Flexibility

The modular Tray and Rack system offers the possibility to expand the system according to process needs. Swivel wheels integrated in the rack allow individual or manifold Flexboy® bags to be easily moved around the facility.

Additional rack systems focused on specific application demands are available on request.

## Safety

Smooth finish and integrated handles make the system both user friendly and safe. This allows reliable stacking of additional modules by maintaining highest safety levels.

## Specifications Tray and Lid

Material	Stainless steel 304 L
Surface finish	Bead blasted
Volumes	5 L, 10 L and 20 L

## Dimensions

5 L	475 × 330 × 117 mm (18.7" × 12.9" × 4.6")
10 L	760 × 330 × 117 mm (29.9" × 12.9" × 4.6")
20 L	770 × 404 × 170 mm (30" × 15.9" × 6.7")

## Weight of Tray | Bag and Lid (Bag Filled to Nominal Volume)

5 L	8.6 kg
10 L	15.7 kg
20 L	28.6 kg

## Technology Integration Support

Sartorius Stedim Biotech supports users already at 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

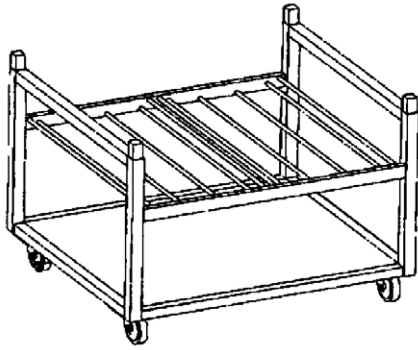
Sartorius Stedim Biotech has established multiple manufacturing sites with consistent industrial processes. The expertise of designing single-use solutions based on collaborative supplier management and customer demand planning ensures a state of the art and robust supply chain that can cope with strong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

## Quality Assurance

The Flexboy® Tray and Rack systems are designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System. They undergo extensive testing before shipping.

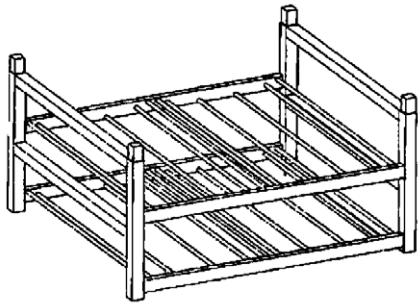
## Available From Stock

Flexboy® Tray and Rack systems are available from stock.



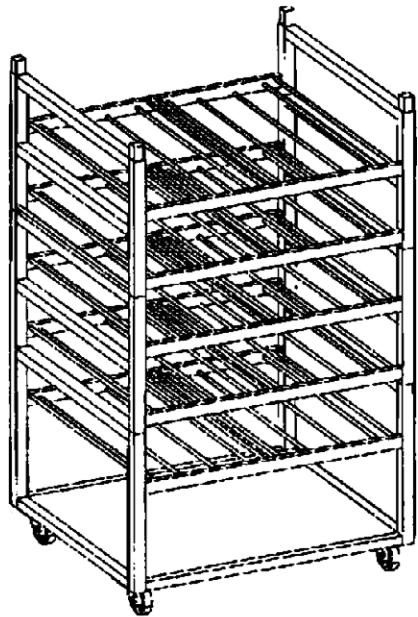
**Base module 1**

H: 772 mm × L: 887 mm × W: 775 mm  
(30.4" × 34.9" × 30.5")



**Module 2**

H: 640 mm × L: 887 mm × W: 775 mm  
(25.2" × 34.9" × 30.5")



**Full version**

H: 1980 mm × L: 887 mm × W: 775 mm  
(77.9" × 34.9" × 30.5")

	Base Module 1	Additional Module 2	Full System
Material	Stainless steel 304	Stainless steel 304	Stainless steel 304
Surface finish	Blast	Blast	Blast
Weight	30.5 kg	26 kg	82.5 kg
Flexboy® tray compartments	2 places for 20 L trays or 4 places for 5 or 10 L trays	4 places for 20 L trays or 8 places for 5 or 10 L trays	10 places for 20 L trays or 20 places for 5 or 10 L trays
Peristaltic pumps compartment	Yes - Dimension: 887 × 775 × 300 mm (34.9" × 30.5" × 11.8")	No	Yes - Dimension: 887 × 775 × 300 mm (34.9" × 30.5" × 11.8")
Wheels	Four swivel wheels (two with brakes)	-	Four swivel wheels (two with brakes)
Additional features	Built in wheels (non autoclavable)	-	Full = 1 base + 2 additional modules

Ordering Information

Order Code	Description
Flexboy® Racks	
FFA102707	Flexboy® Rack   Base Module 1
FFA102714	Flexboy® Rack   Module 2
Flexboy® Trays	
FFA102705	Flexboy® Tray and Lid 5 L
FFA102715	Flexboy® Tray and Lid 10 L
FFA102716	Flexboy® Tray and Lid 20 L


Please note: Custom versions are available upon request.

**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.: SPL2004-e | Order No.: 85030-533-50 | Status: 08 | 31 | 2021