

# Biosafe® 110 Ports

Contained Aseptic Transfer  
of Components and Fluids



## Product Information

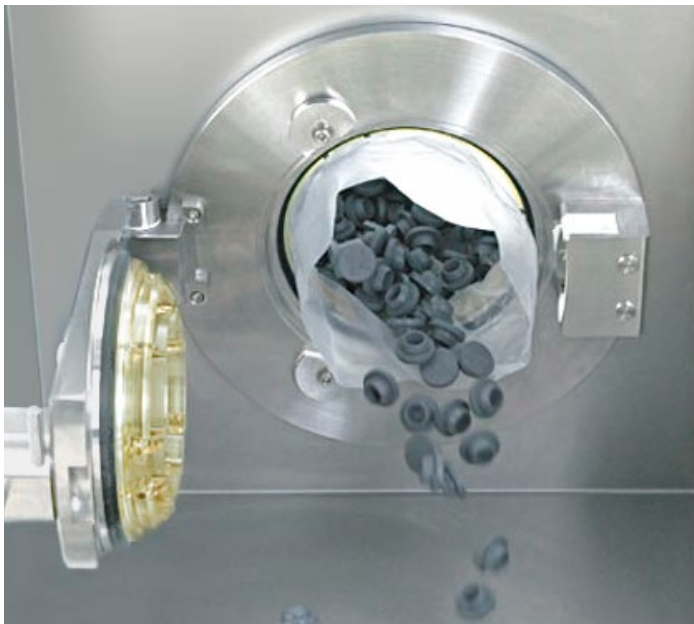
The Biosafe® 110 range of aseptic transfer ports offers reliable and easy-to-use solutions for the secure transfer of components and fluids while maintaining the integrity of the critical area – isolators, RABS and cleanrooms.

## Benefits

- Enhanced sterility assurance and viral segregation in aseptic processing
- Easy to use
- Reduced footprint
- Process safety
- Versatile technology
- Simplified maintenance and sterilization
- Cost effective


Applications	Upstream and Downstream Processing	Aseptic Processing	
Applications	Transfer of large volume support solutions held in lower classified environments to higher classification process zones. Examples of applications: <ul style="list-style-type: none"> <li>Media feed to N-2   N-1   bioreactor</li> <li>Buffer feed to chromatography columns</li> <li>Transfer out of higher classified zones: <ul style="list-style-type: none"> <li>fraction collection</li> <li>bulk intermediates</li> <li>bulk final product transferring out of purification</li> </ul> </li> </ul>	Discharge from autoclave <ul style="list-style-type: none"> <li>Stoppers</li> </ul>	Transfer into isolators or RABS <ul style="list-style-type: none"> <li>Drug products</li> <li>Entry of stoppers</li> <li>Entry and removal of QC test devices, tools and pumps</li> <li>Waste removal</li> </ul>
Biosafe® 110 port	<input checked="" type="checkbox"/>	-	<input checked="" type="checkbox"/> <sup>1</sup>
Biosafe® 110 bags (All Biosafe® bags are designed for connection to any Biosafe® 110 port.)	Rapid Aseptic Fluid Transfer (RAFT) systems preassembled with other Sartorius technologies such as Flexsafe® 3D and Flexboy® system. Complete assembly is gamma sterile.	Double connector gamma sterile	<ul style="list-style-type: none"> <li>Open autoclavable to be filled by end-users prior to autoclave sterilization and aseptic transfer.</li> <li>Prefilled autoclavable delivered ready-to-sterilize via component suppliers.</li> <li>Prefilled delivered gamma sterile by component suppliers.</li> <li>Closed gamma sterile for the removal of waste, tools, pumps and QC test devices.</li> </ul>

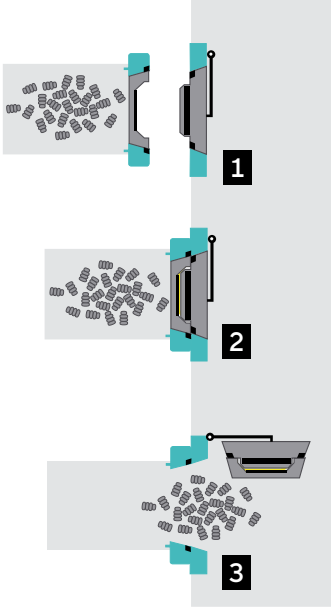
<sup>1</sup> Biosafe® 110 port with outside opening is the best choice to prevent air turbulence in RABS and glove usage in isolator.



# Operating Sequences

## Connecting a Biosafe® 110 Bag





**Approach**

- Wipe down the Biosafe® 110 port
- Open package of the Biosafe® 110 bag and remove the protecting pouch

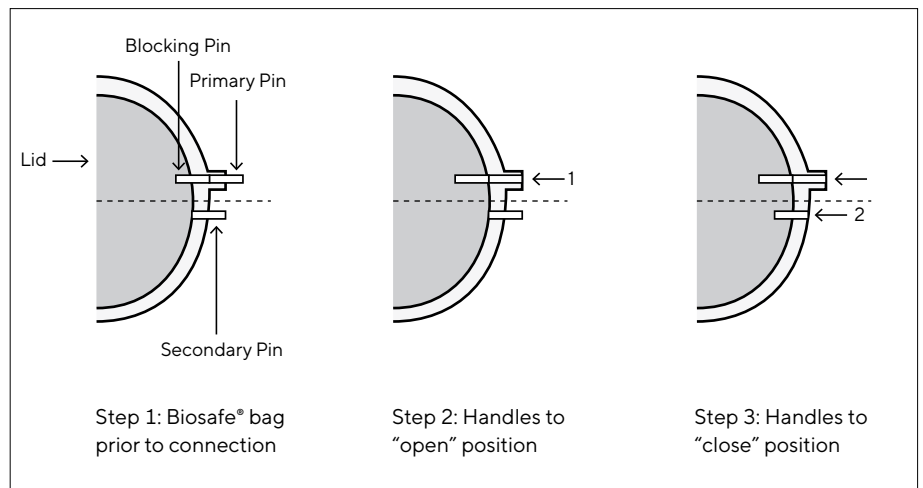
**Docking**

- Docking is secured by magnetic guidance on the Biosafe® 110 port
- The magnetic connection is further secured by mechanical locks.

**Opening and transfer**

- Open the double-door either from inside or outside the critical area.
- Aseptic transfer of components, fluids or powders.

Prior to the connection, the pins on the Biosafe® connector are in the “out” position (see beside). This is a proof that the Biosafe® 110 bag is ready to be connected to the Biosafe® 110 port and that it has not been used before.



## Maintenance, Decontamination and Sterilization of the Biosafe® 110 Port

When connected to the Biosafe® 110 port, the dummy service connector allows the door to be opened for the sterilization of the critical area and the inner side of the Biosafe® 110 port as well as for maintenance operations such as gasket replacement.

Outside View



Inside View



# Specifications

<b>Biosafe® 110 Port</b>	
Installation Requirements	<p>Wall thickness: 3–8 mm   0.08–0.31 in.</p> <ul style="list-style-type: none"> <li>▪ If the wall thickness exceeds 8 mm   0.31 in., the Biosafe® 110 port must be installed on a Biosafe® support which is then integrated into the wall.</li> <li>▪ If outside opening is chosen, the Biosafe® 110 port is systematically supplied on a Biosafe® support which is then integrated into the wall.</li> <li>▪ We highly recommend setting the port below or on a window so that the operator can see the other side.</li> <li>▪ Height of port for accessibility: for good access, the port axis must be 1.1 m to 1.4 m (43.3 in. to 52.12 in.) high from the operator standing reference.</li> </ul>
Weight (approx.)	15 kg   33.07 lb.
Operating Temperature Range	5 °C to 30 °C   41°F to 86°F
Maximum Temperature During Autoclave Cycle	N   A
Pressure Range During Autoclave Cycle	N   A
Materials of Construction <sup>1</sup>	Stainless Steel 316L; PETP; Silicone   EPDM for gaskets
Passage Diameter	110 mm   4.3 in.
Quality Standards	<ul style="list-style-type: none"> <li>▪ All materials are compliant with 21 CFR Part 177.2600 (EPDM Silicone), 21 CFR Part 177.1630 (PETP) and 21 CFR Part 177.2470 (PEEK)</li> </ul>
FAT   SAT	<p>Factory Acceptance Tests (FAT) and Site Acceptance Tests (SAT) are performed on each Biosafe® 110 port.</p> <ul style="list-style-type: none"> <li>▪ Only during FAT, air tightness is performed at several points of control: gasket, locking screws, handles positioning</li> <li>▪ Functional: positioning of gasket, positioning and manipulation of external and internal handle(s), lockers, mechanical securities</li> </ul>
Cleaning and Decontamination Agents	<ul style="list-style-type: none"> <li>▪ Purified water (WFI) or any neutral pH detergent</li> <li>▪ Ethanol or isopropyl alcohol (70% v v)</li> <li>▪ Peracetic acid solution (2% v v)</li> <li>▪ Hydrogen peroxide solution (2% v v)</li> </ul>

<sup>1</sup>The list provides construction materials in contact with the critical area (isolators, RABS, cleanrooms).

# Ordering information

## Biosafe® Ports

FAA109722	Biosafe® 110 port <sup>1</sup>
FAA109860	Biosafe® 110 port <sup>2</sup>
FAA109861	Biosafe® 110 port (w/ support) <sup>3</sup>
FAA109862	Biosafe® 110 port (w/ support) <sup>4</sup>
FAA308155	Biosafe® 110 port - 90° lever <sup>1</sup>
FAA308156	Biosafe® 110 port - 90° lever <sup>2</sup>
FAA308157	Biosafe® 110 port - 90° lever (w/o support) <sup>3</sup>
FAA308158	Biosafe® 110 port - 90° lever (w/o support) <sup>4</sup>

<sup>1</sup> Inside right opening

<sup>3</sup> Outside right opening

<sup>2</sup> Inside left opening

<sup>4</sup> Outside left opening

## Accessories, Options & Support for Biosafe® Ports

FAA109724	Biosafe® dummy service connector for Biosafe® 110 port
FAA114793	Biosafe® vertical support device for Biosafe® 110 port
FAA112077	Biosafe® inclined support device for Biosafe® 110 port

With each port ordered, please indicate the wall thickness on which the port will be installed, as this information is mandatory to proceed with manufacturing.

For further information or details on Site Acceptance Tests (SAT) and training at final customers, please contact your local Service representative.

## Spare Parts for Biosafe® 110 Ports

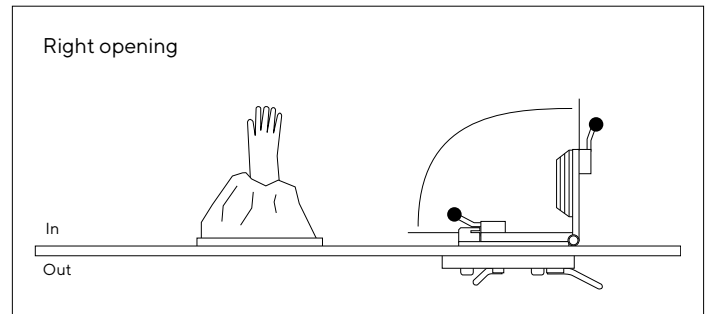
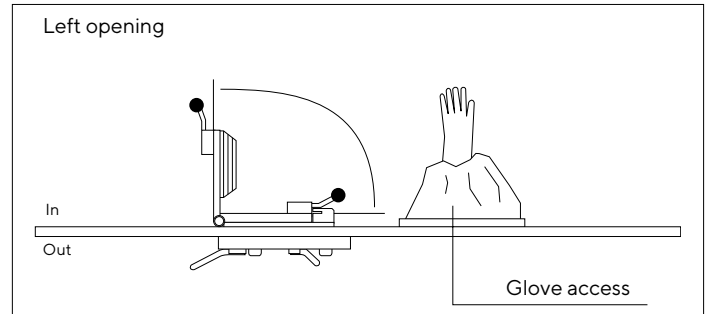
Upon order of Biosafe® 110 ports, you will receive a complete technical package including the list and prices of spare parts.

### Trademarks

Biosafe® and Biosteam® are Sartorius Stedim Aseptics registered trademarks.

## Left or Right-Hand Side Opening

You can choose to have an opening on the left or on the right-hand side as shown below:



## Lever on the Right or 90° Lever

You can choose to have the monolever on the right of the port or at 90° compared to the port as shown in the picture below:




**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)