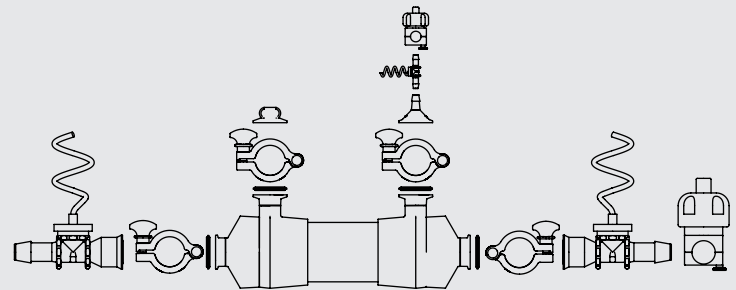


Flow Connector Kit Bioproducer

Hardware Selected
to Connect Your Specific
Hollow Fiber Module
to a Crossflow Assembly



Executive Summary

Flow Connector Kit Bioproducer connects your hollow fiber Bioproducer module to your assembly tubing. The kits come with PendoTECH pressure sensors that enable monitoring of transmembrane pressure (TMP).

Unique Selling Points

Hardware selected to connect your Bioproducer hollow fiber module to a crossflow assembly. The kits include, as needed:

- Clamps
- Pinch valves
- Reducers
- Silicone gaskets
- Caps
- Single-use pressure sensors to enable monitoring of transmembrane pressure (TMP)

Relevant Applications

Sartorius Hollow Fiber Modules are designed for biopharmaceutical processing. They can be used for the following applications:

- Upstream**

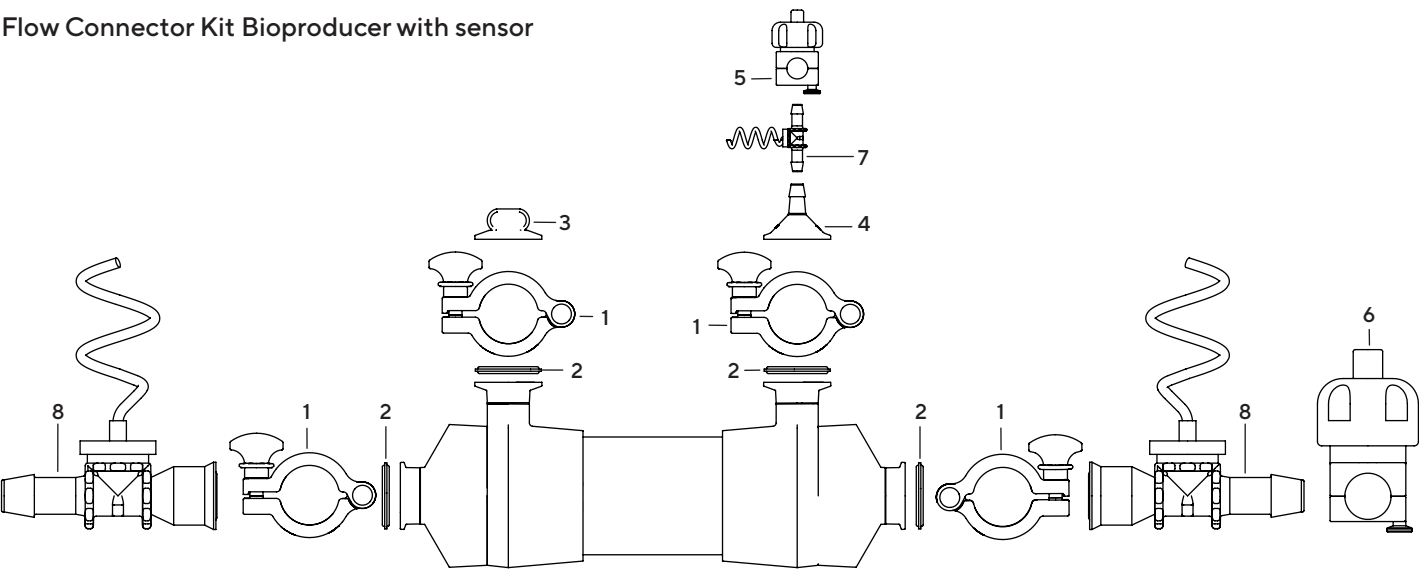
 - Cell harvest
 - Cell or lysate clarification
- Downstream**

 - Concentration
 - Diafiltration

Technical Specifications

| Attribute | Item | Material |
|----------------------|-------------------|------------------------------|
| In fluid contact | Gaskets | Platinum-cured silicone |
| In fluid contact | Reducers and caps | Polypropylene AF |
| In fluid contact | Sensors | Polysulfone |
| Not in fluid contact | Pinch Valves | Stainless Steel |
| Not in fluid contact | Clamps | Glass Reinforced White Nylon |

Flow Connector Kit Bioproducer with sensor



| Item Number | Description | Use | Quantity |
|-------------|--|---------------------------------|-----------------------|
| 1 | 1.5 inch TC Clamps | For Permeate Retentate Feed | 4 |
| 2 | 1.5 inch TC Silicone Gaskets | For Permeate Retentate Feed | 8 (4 extra not shown) |
| 3 | 1.5 inch TC Cap | For Permeate | 1 |
| 4 | 1.5 inch TC to 0.375 inch Barb (Reducer) | For Permeate | 1 |
| 5 | 0.375 inch Pinch Valve | For Permeate | 1 |
| 6 | 1 inch Pinch Valve | For Retentate | 1 |
| 7 | Single-use Pressure Sensors 0.375 inch Barb × 0.375 inch Barb | For Permeate | 1 |
| 8 | Single-use Pressure Sensors 1 inch TC × 1 inch Barb | For Retentate Feed | 1 |



Ordering Information


| | Green Line, Steamer Line, Reuse Line | |
|---------------|--------------------------------------|-----------------|
| Specification | with Sensors | without Sensors |
| Order Code | FCPROKIT-SEN01 | - |

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 Copyright Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

DIR: 2730229-000-00 | Status: 05 | 27 | 2021