

Sartorius Hollow Fiber TFF Flow Connector Kit | Discover

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



Executive Summary

Hollow Fiber TFF Flow Connector Kit | Discover connects your hollow fiber Discover 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 Hollow Fiber TFF Module | Discover to a crossflow assembly. The kits include, as needed:

- Luer fittings
- Pinch valves
- Reducers
- Silicone gaskets
- Plugs
- Single-use pressure sensors to enable monitoring of transmembrane pressure (TMP)

Relevant Applications

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

Upstream

- Cell harvest
- Cell or lysate clarification

Downstream

- Concentration
- Diafiltration

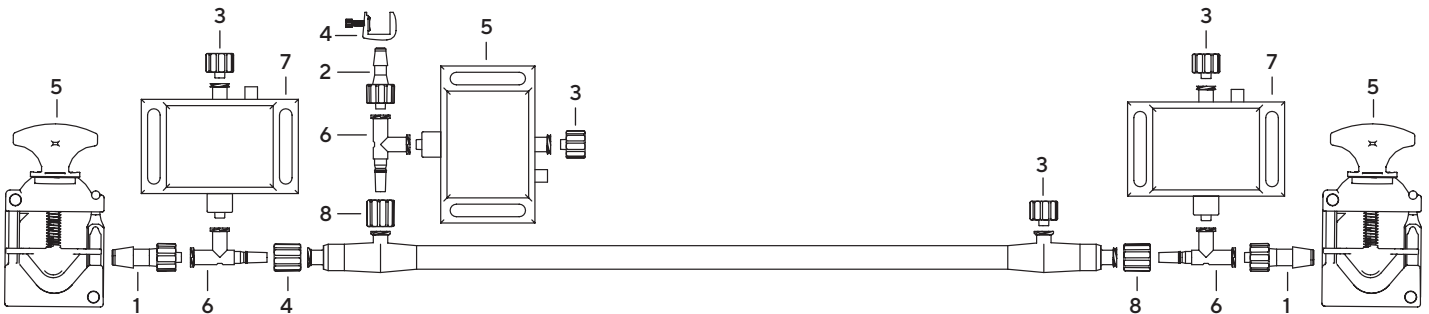
Technical Specifications

| Attribute | Item | Material |
|----------------------|--------------------|-------------------------|
| In fluid contact | Gaskets | Platinum-cured silicone |
| In fluid contact | Reducers and plugs | Polypropylene AF |
| In fluid contact | Sensors | Polysulfone |
| Not in fluid contact | Pinch Valves | Nylon |

Sartorius Hollow Fiber Flow Connector Kit | Discover Without Sensors



Sartorius Hollow Fiber Flow Connector Kit | Discover With Sensors



| Specification | Item Number | Description | Use | Quantity |
|-----------------|-------------|--|---------------------------------|----------|
| Without Sensors | 1 | Male Luer Integral Lock Ring to 0.25 inch Barb (Reducer) | For Retentate Feed | 2 |
| | 2 | Male Luer Integral Lock Ring to 0.15625 inch Barb (Reducer) | For Permeate | 1 |
| | 3 | Male Luer Lock Ring Plugs | For Permeate Retentate Feed | 1 |
| | 4 | Pinch Valve for 0.09375 inch to 0.25 inch Tube OD | For Permeate | 1 |
| | 5 | 0.375 inch Pinch Valve | For Retentate | 2 |
| With Sensors | 1 | Male Luer Integral Lock Ring to 0.25 inch Barb (Reducer) | For Retentate Feed | 2 |
| | 2 | Male Luer Integral Lock Ring to 0.15625 inch Barb (Reducer) | For Permeate | 1 |
| | 3 | Male Luer Lock Ring Plugs | For Permeate Retentate Feed | 4 |
| | 4 | Pinch Valve for 0.09375 inch to 0.25 inch Tube OD | For Permeate | 1 |
| | 5 | 0.375 inch Pinch Valve | For Retentate | 2 |
| | 6 | Female Luer Thread Style to Male Luer Tee (with separate Luer lock ring) | For Permeate Retentate Feed | 3 |
| | 7 | Single-use Pressure Sensor with Luer | For Permeate Retentate Feed | 3 |
| | 8 | Luer Lock Ring | For Permeate Retentate Feed | 3 |

Note

May use retentate Pinch Valve to generate back pressure during operation of Hollow Fiber TFF Module.

Caution

Do not operate Hollow Fiber TFF Module without pressure monitoring capabilities.

Recommendation

Contact Pressure Sensor manufacturer, PendoTECH, for pressure monitor requirements.

Ordering Information


| | Single-use Line, Steamer Line, Re-use Line | |
|---------------|--|----------------|
| Specification | Without Sensors | With Sensors |
| Order Code | FCDISKIT-00001 | FCDISKIT-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