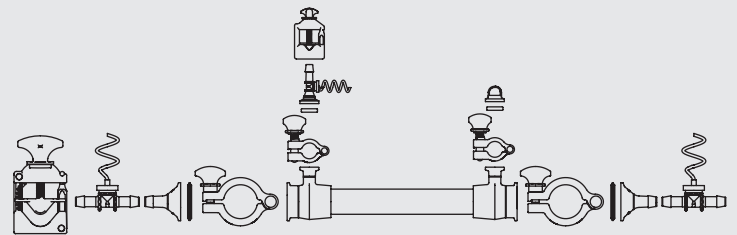


Flow Connector Kit Investigator

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



Executive Summary

Flow Connector Kit – Investigator connects your hollow fiber Investigator module to your assembly tubing. The kit comes with PendoTECH pressure sensors that enable monitoring of transmembrane pressure (TMP).

Unique Selling Points

Hardware selected to connect your Investigator 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

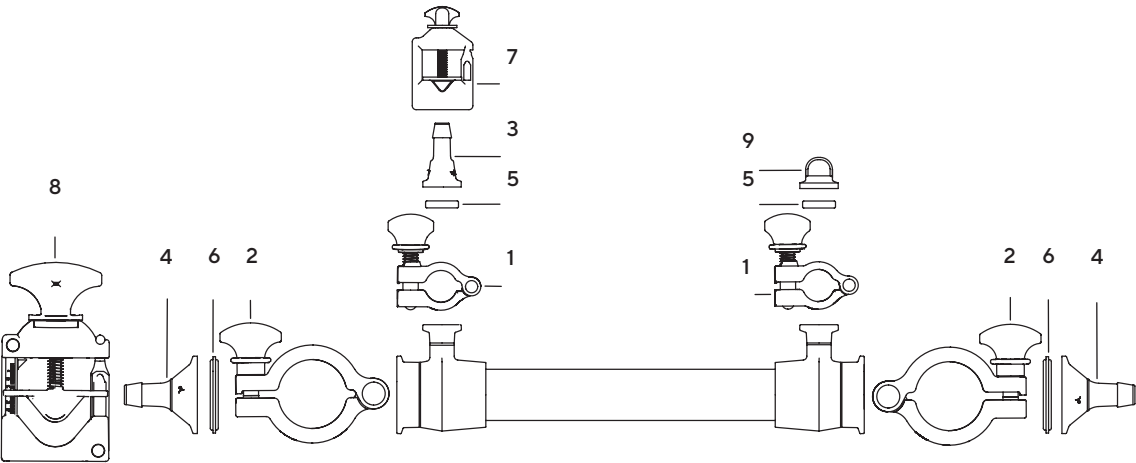
Sartorius Hollow Fiber Modules are designed for biopharmaceutical processing. They can be used for the following applications:

- Upstream**
 - Cell harvest
 - Cell or lysate clarification
- Downstream**
 - Concentration
 - Diafiltration

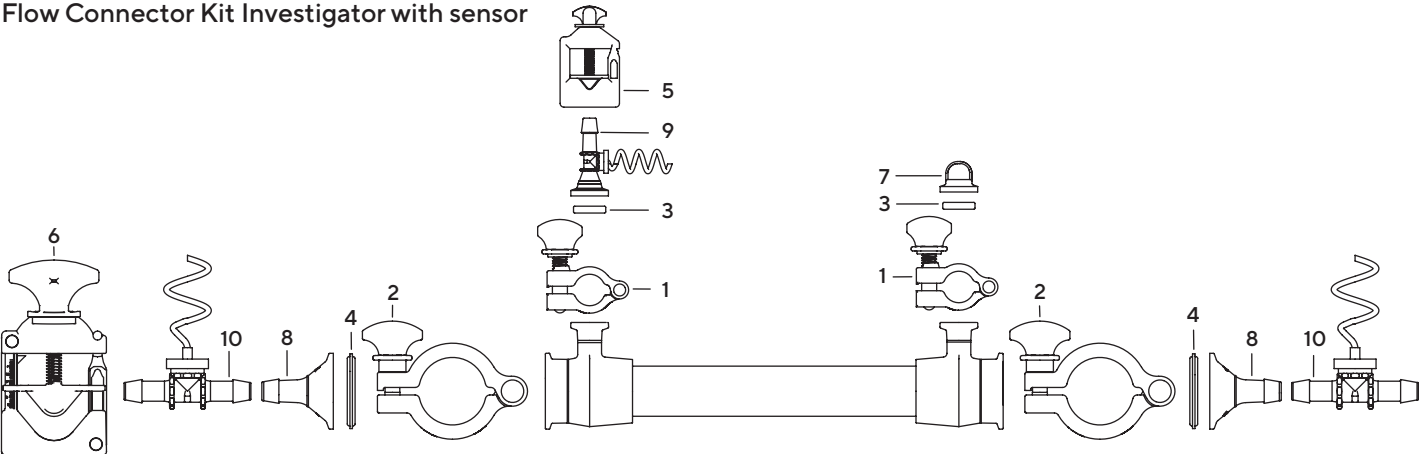
Technical Specifications

| 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 | Glass Reinforced White Nylon |
| Not in fluid contact | Clamps | Glass Reinforced White Nylon |

Flow Connector Kit Investigator no sensor



Flow Connector Kit Investigator with sensor



| Specification | Item Number | Description | Use | Quantity |
|-----------------|-------------|---------------------------------------------------|----------------------|-----------------------|
| without Sensors | 1 | 0.75 inch TC White Plastic Clamps | For Permeate | 2 |
| | 2 | 1.5 inch TC Clamps | For Retentate Feed | 2 |
| | 3 | 0.75 inch TC × 0.375 inch Barb (Reducer) | For Permeate | 1 |
| | 4 | 1.5 inch TC × 0.5 inch Barb (Reducer) | For Retentate Feed | 2 |
| | 5 | 0.75 inch TC Silicone Gaskets | For Permeate | 4 (2 extra not shown) |
| | 6 | 1.5 inch TC Silicone Gaskets | For Retentate Feed | 4 (2 extra not shown) |
| | 7 | 0.375 inch Pinch Valve | For Permeate | 1 |
| | 8 | 0.75 inch Pinch Valve | For Retentate | 1 |
| | 9 | 0.75 inch TC Cap | For Permeate | 1 |
| with Sensors | 1 | 0.75 inch TC White Plastic Clamps | For Permeate | 2 |
| | 2 | 1.5 inch TC Clamps | For Retentate Feed | 2 |
| | 3 | 0.75 inch TC Silicone Gaskets (2 extra not shown) | For Permeate | 4 (2 extra not shown) |
| | 4 | 1.5 inch TC Silicone Gaskets (2 extra not shown) | For Retentate Feed | 4 (2 extra not shown) |
| | 5 | 0.375 inch Pinch Valve | For Permeate | 1 |
| | 6 | 0.75 inch Pinch Valve | For Retentate | 1 |
| | 7 | 0.75 inch TC Cap | For Permeate | 1 |
| | 8 | 1.5 inch TC × 0.5 inch Barb (Reducer) | For Retentate Feed | 2 |
| | 9 | Single-use Pressure Sensor 0.75 inch TC | For Permeate | 1 |
| | 10 | Single-use Pressure Sensors, 0.5 inch Barb | For Retentate Feed | 2 |



Ordering Information


| | Green Line, Steamer Line, Reuse Line | |
|---------------|--------------------------------------|-----------------|
| Specification | with Sensors | without Sensors |
| Order Code | FCINVKIT-SEN01 | FCINVKIT-00001 |

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