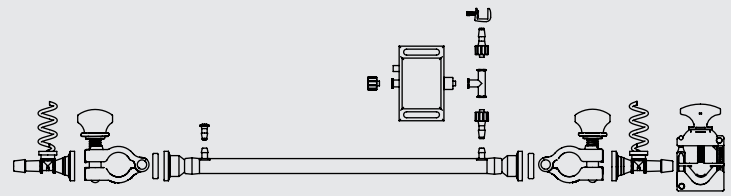


# Flow Connector Kit Explorer

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



## Executive Summary

Flow Connector Kit Explorer connects your hollow fiber Explorer 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 Explorer hollow fiber module to a crossflow assembly. The kits include, as needed:

- Clamps
- Pinch valves
- Reducers
- Silicone gaskets
- Plugs
- 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

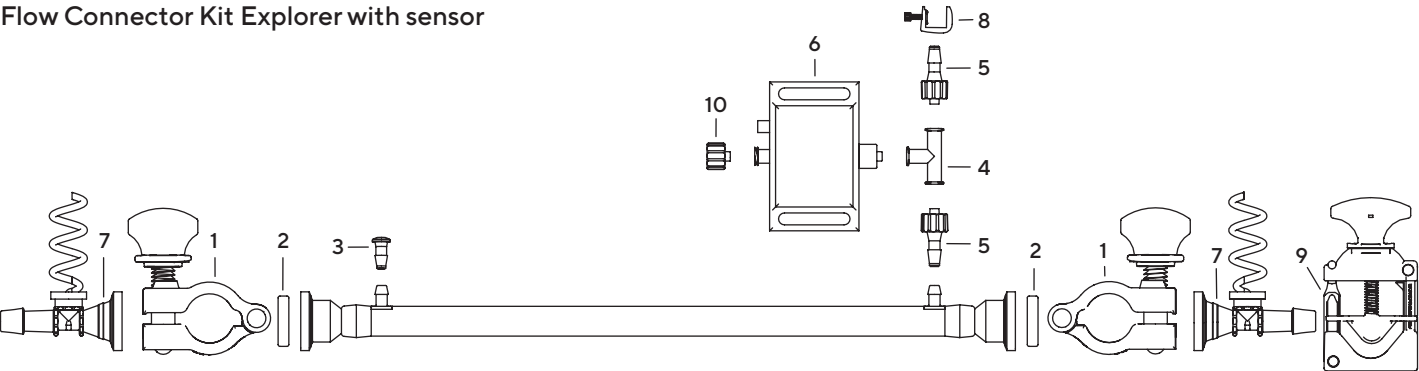
**Downstream**

- Concentration
- Diafiltration

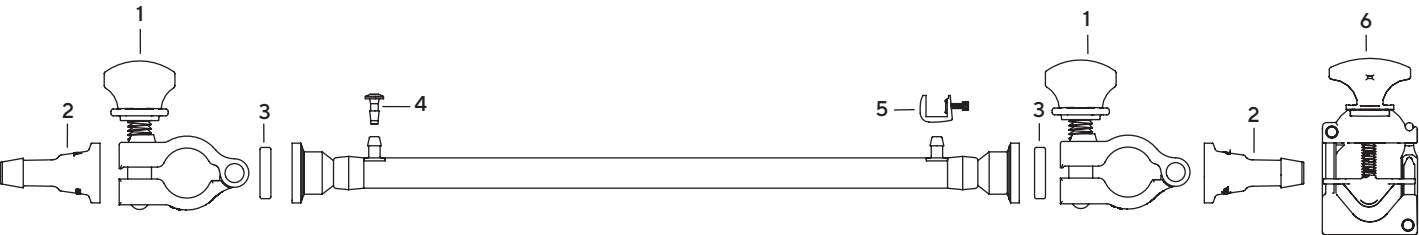
## Technical Specifications

Attribute	Item	Material
In fluid contact	Gaskets	Platinum-cured silicone
In fluid contact	Reducers	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 Explorer with sensor



Flow Connector Kit Explorer no sensor



Specification	Item Number	Description	Use	Quantity
with sensor	1	0.75 inch TC Clamps	For Retentate   Feed	2
	2	0.75 inch Silicone Gaskets	For Retentate   Feed	4 (2 extra not shown)
	3	Press in Plug, Barb	For Permeate	2 (1 extra not shown)
	4	TEE-Female Luer	For Permeate	1
	5	Luer × 0.1563 Barb (Reducers)	For Permeate	2 (1 extra not shown)
	6	Single Use Pressure Sensor Luer	For Permeate	1
	7	Single Use Pressure Sensor 0.75 inch TC to 0.375 inch Barb	For Retentate   Feed	1
	8	Pinch Valve for 0.375 inch to 0.25 inch Tube OD	For Permeate	1
	9	0.375 inch Pinch Valve	For Retentate	1
	10	Male Luer Lock Ring Plugs	For Permeate	3 (2 extra not shown)
no sensor	1	0.75 inch TC White Plastic Clamps	For Retentate   Feed	2
	2	0.75 inch TC × 0.375 inch Barb (Reducers)	For Retentate   Feed	2
	3	0.75 inch TC Silicone Gaskets	For Retentate   Feed	4 (2 extra not shown)
	4	Press in Plugs, 0.1875 inch Barb	For Retentate   Feed	2 (1 extra not shown)
	5	Pinch Valve for 0.09375 inch to 0.75 inch Tube OD	For Permeate	1
	6	0.375 inch Pinch Valve	For Retentate	1

## Ordering Information


	Green Line, Steamer Line, Reuse Line	
Specification	with Sensors	without Sensors
Order Code	FCEXPKIT-SEN01	FCEXPKIT-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](http://www.sartorius.com)

Specifications subject to change without notice.

© 2021 Copyright Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

DIR: 2765208-000-00 | Status: 05 | 27 | 2021