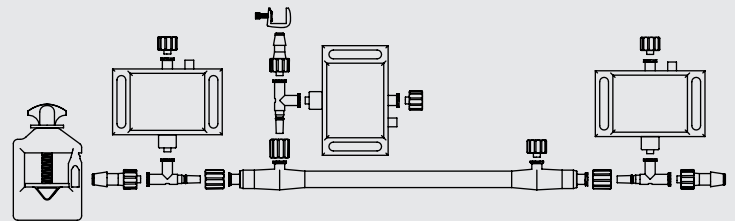


Flow Connector Kit Discover

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



Executive Summary

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 Discover hollow fiber module 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 Modules are designed for biopharmaceutical processing. They can be used for the following applications:

- Upstream**

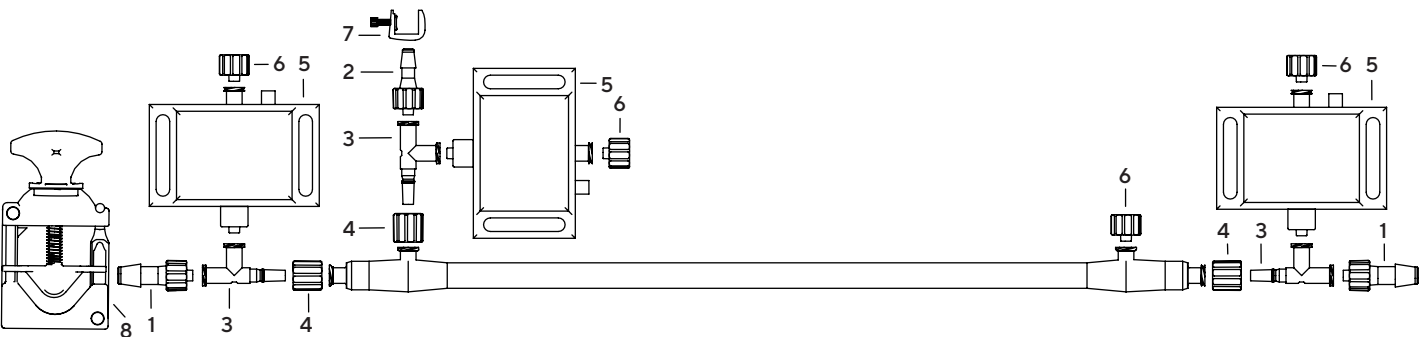
 - Cell harvest
 - Cell or lysate clarification
- Downstream**

 - 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

Flow Connector Kit Discover with sensor



Specification	Item Number	Description	Use	Quantity
without Sensors	1	Male Luer Integral Lock Ring to 0.25 inch Barb (Reducer)	For Retentate Feed	4 (2 extra not shown)
	2	Male Luer Integral Lock Ring to 0.15625 inch Barb (Reducer)	For Permeate	2 (1 extra not shown)
	3	Female Luer Thread Style to Male Luer Tee (with separate Luer lock ring)	For Permeate Retentate Feed	3
	4	Luer Lock Ring	For Permeate Retentate Feed	3
	5	Single-use Pressure Sensor with Luer	For Permeate Retentate Feed	3
	6	Male Luer Lock Ring Plugs	For Permeate Retentate Feed	8 (4 extra not shown)
	7	Pinch Valve for 0.09375 inch to 0.25 inch Tube OD	For Permeate	1
	8	0.375 inch Pinch Valve	For Retentate	1



Ordering Information


	Green Line, Steamer Line, Reuse Line
Specification	with Sensors
Order Code	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

Specifications subject to change without notice.

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

DIR: 2730230-000-00 | Status: 05 | 27 | 2021