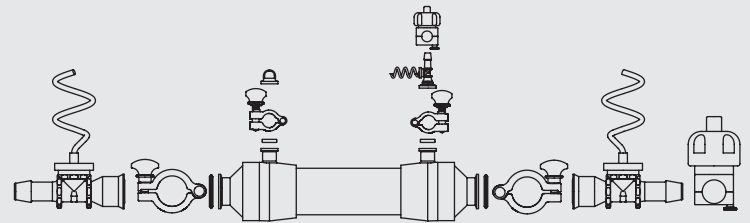


# Flow Connector Kit mini- Bioproducer

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



## Executive Summary

Flow Connector Kit – mini-Bioproducer connects your hollow fiber mini-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 mini-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**

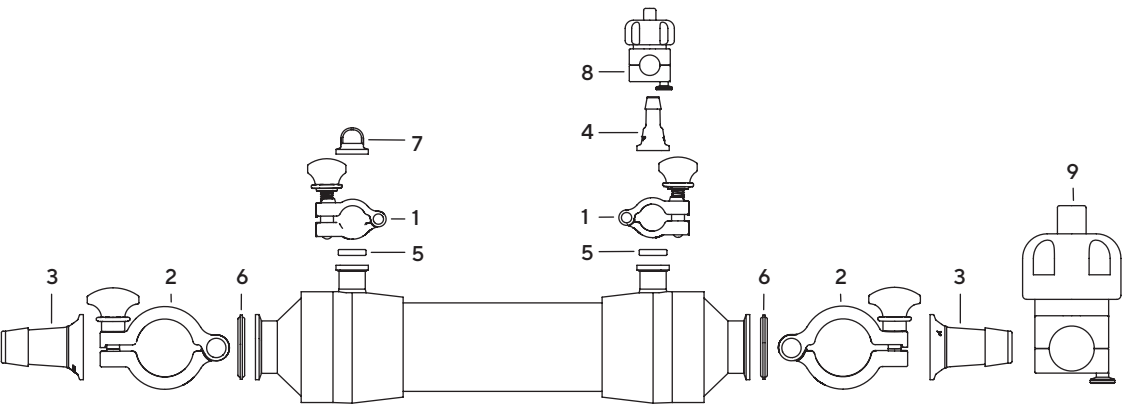
  - Cell harvest
  - Cell or lysate clarification
- Downstream**

  - Concentration
  - Diafiltration

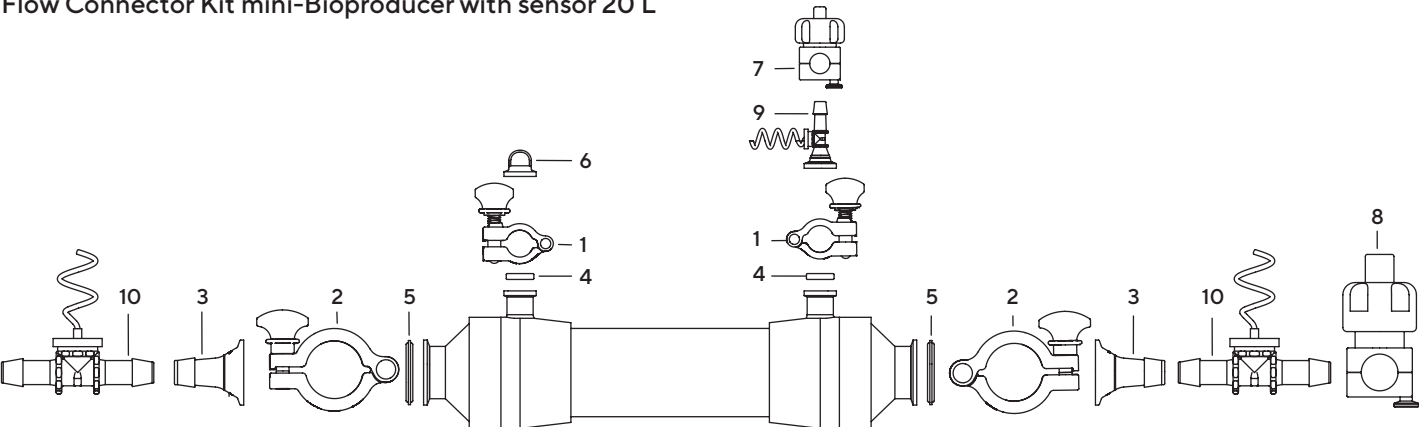
## Technical Specifications

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

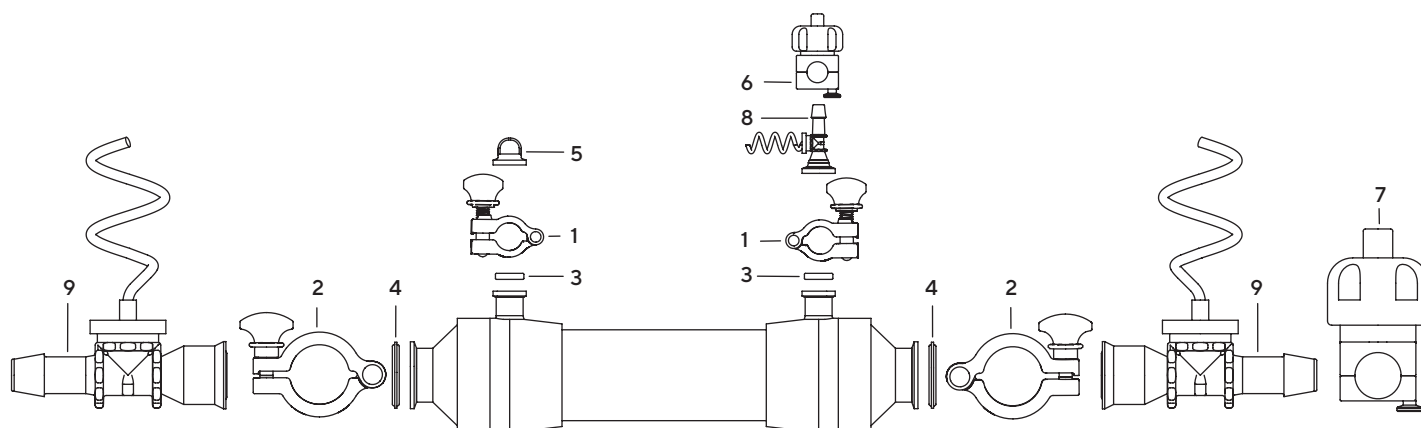
Flow Connector Kit mini-Bioproducer no sensor



Flow Connector Kit mini-Bioproducer with sensor 20 L



## Flow Connector Kit mini-Bioproducer with sensor 40 L



Specification	Item Number	Description	Use	Quantity
without Sensors	1	0.75 inch TC Clamps	For Permeate	2
	2	1.5 inch TC Clamps	For Retentate   Feed	2
	3	1.5 inch TC × 1 inch Barb (Reducers)	For Retentate   Feed	2
	4	0.75 inch TC × 0.375 inch Barb (Reducers)	For Permeate	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	0.75 inch TC Cap	For Permeate	1
	8	0.375 inch Pinch Valve	For Permeate	1
	9	1 inch Pinch Valve	For Retentate	1
with Sensors 20 L/min	1	0.75 inch TC Clamps	For Permeate	2
	2	1.5 inch TC Clamps	For Retentate   Feed	2
	3	1.5 inch TC × 0.75 inch Barb (Reducers)	For Retentate   Feed	2
	4	0.75 inch TC Silicone Gaskets	For Permeate	4 (2 extra not shown)
	5	1.5 inch TC Silicone Gaskets	For Permeate	4 (2 extra not shown)
	6	0.75 inch TC Cap	For Permeate	1
	7	0.375 inch Pinch Valve	For Permeate	1
	8	0.75 inch Pinch Valve	For Permeate	1
	9	Single-use Pressure Sensor, 0.75 inch TC to TC to 0.375 inch Barb	For Permeate	1
	10	Single-use Pressure Sensors, 0.75 inch Barb	For Retentate   Feed	2
with Sensors 40 L/min	1	0.75 inch TC Clamps	For Permeate	2
	2	1.5 inch TC Clamps	For Retentate   Feed	2
	3	0.75 inch TC Silicone Gaskets	For Permeate	4 (2 extra not shown)
	4	1.5 inch TC Silicone Gaskets	For Retentate   Feed	4 (2 extra not shown)
	5	0.75 inch TC Cap	For Permeate	1
	6	0.375 inch Pinch Valve	For Permeate	1
	7	1 inch Pinch Valve	For Retentate	1
	8	Single-use Pressure Sensors 0.75 inch TC × 0.375 inch Barb	For Permeate	2
	9	Single-use Pressure Sensors 1 inch TC × 1 inch Barb	For Retentate   Feed	2

# Ordering Information


Green Line, Steamer Line, Reuse Line			
Specification	with Sensors		without Sensors
	20 L/min	40 L/min	
Order Code	FCMPRK IT-S2001	FCMPRK IT-S4001	FCMPRK IT-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: 2738561-000-00 | Status: 05 | 27 | 2021