

# SARTOFLOW<sup>®</sup> Systems

Crossflow Solutions for Lab and Process Development



# Scalable Systems with Multiple Options



	SARTOFLOW® Smart	SARTOFLOW® Study	SARTOFLOW® Advanced
<b>Filter surface area</b>	50 cm <sup>2</sup> – 0.14 m <sup>2</sup>	0.1 – 2.1 m <sup>2</sup>	0.1 – 2.1 m <sup>2</sup>
<b>Automation</b>	Manual	Manual	Semi-automatic
<b>Application area</b>	PD, R&D, cGMP	R&D, GLP	R&D, cGMP production
<b>Min. working volume*</b>	<20 mL	200 mL or 400 ml (PP tank)	200 mL
<b>Cleaning   sterilisation</b>	CIP	CIP	CIP
<b>Pump</b>	4-piston Diaphragm 1 – 100 L/h @ 3 bar	4-piston Diaphragm 700 L/h @ 4 bar	4-piston Diaphragm 700 L/h @ 4 bar
<b>Parameters automatically regulated by the system**</b>	P <sub>in</sub> , ΔP (Optional: level control)	P <sub>in</sub> , ΔP (Optional: Retentate flow, level control)	P <sub>in</sub> , P <sub>out</sub> , ΔP, TMP, level control (Optional: Retentate flow)

\* with 1 Sartocor® Slice cassette

\*\* P<sub>in</sub>: Inlet pressure | P<sub>out</sub>: Retentate pressure | P<sub>per</sub>: Permeate pressure | ΔP: Differential pressure | TMP: Trans-Membrane pressure

## Options

DCU-4 with touch screen	•	•	•
Sartocor® Slice filter holder (standard)	Sartocor® Slice 200	•	•
Holder for Sartocor® Slice 50	(•)	-	-
Sartocor® Slice single-use filter holder (manual or hydraulic)	-	-	-
Sartocor® filter holder for up to 3 Sartocor® ECO cassettes (2.1 m <sup>2</sup> )	-	(•)	(•)
Pressure indicator feed and retentate	•	•	•
Pressure sensor permeate	•	•	•
Load cell	•	•	•
Flow-meter retentate	-	(•)	(•)
Flow-meter permeate	Balance	•	•
Automatic valve retentate	-	-	•
Automatic valve permeate	-	-	-
Conductivity probe retentate	(Vessel)	-	(•)
Conductivity probe permeate	-	-	(•)
pH probe retentate	(Vessel)	-	- (ETO)
pH probe permeate	-	-	- (ETO)
UV probe permeate	-	-	-
Temperature probe	(Vessel)	(•)	•
Pump head heat exchanger	-	-	-
Dry run protection	-	-	-
1 <sup>st</sup> peristaltic pump (filling of the tank and   or diafiltration)	(•)	(•)	•
2 <sup>nd</sup> peristaltic pump (filling of the tank and   or diafiltration)	-	(•)	(•)
1 L Vessel	•	-	-
10 L polypropylene vessel	-	•	-
10 L stainless steel tank	-	(•)	(•)
10 L stainless steel tank with a sight glass	-	(•)	(•)
Double-jacket tank, 10 L, stainless steel, with a sight glass	-	(•)	(•)
Bag holder for single-use crossflow bags	(•)	(•)	(•)
Standard & cGMP documentation (Level 0 & 1)	Level 0	Level 0	•
Documentation package – Basic Qualification Test (Level 2)	(•)	-	(•)
BQT Execution package in FAT (Level 3)	(•)	-	(•)
BQT Execution package in SAT (Level 4) contact service	(•)	-	(•)
SCADA MFCS   win 3.0 software (data acquisition upgradeable with options)	(•)	(•)	(•)
Recipe control (S88) software module (optional for MFCS   win 3.0)	(•)	(•)	(•)
21 CFR Part 11 software module (optional for MFCS   win 3.0)	(•)	(•)	(•)

• Standard (•) Optional – (ETO) Not available or only customized solution

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

Phone +49.551.308.0  
Fax +49.551.308.3289



[www.sartorius.com/crossflow-systems](http://www.sartorius.com/crossflow-systems)