

Chromatography
Systems Capabilities

Simplifying Progress

SARTORIUS

Chromatography Systems Head to Head

Sartorius System

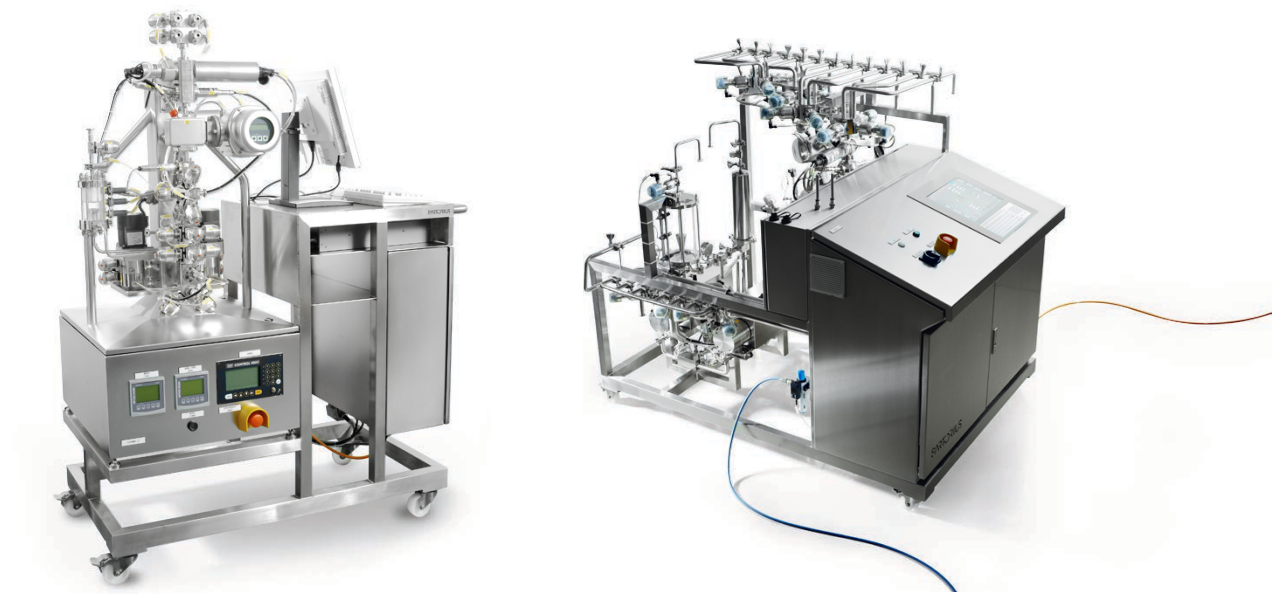
Cytiva System

Single-Use Systems

Flowrate Capabilities	System	System	Flowrate Capabilities
10 to 500 L h	Sartorius S.U.T. Chrom 3/8"	ÄKTA ready with Low Flow Kit	3 to 175 L h
20 to 900 L h	Sartorius S.U.T. Chrom 1/2"	ÄKTA ready with High Flow Kit	7.5 to 510 L h
		ÄKTA ready XL, 3/4" kit	45 to 1900 L h
		ÄKTA ready XL, 1" kit	90 to 3500 L h

Multi-Use Systems

Flowrate Capabilities	System	System	Flowrate Capabilities
1 to 100 L h	PKP2	Bioprocess Modular 1/4 in	1 to 60 L h
1 to 150 L h	PKP3	ÄKTA process, 6 mm or 3/8 in	4 to 180 L h
10 to 100 L h	PK10	ÄKTA process, 10 mm or 1/2 in	15 to 600 L h
25 to 400 L h	PK25	ÄKTA process, 1 in	45 to 1800 L h
50 to 1200 L h	PK50	Bioprocess Modular 1.5 in	200 to 5000 L h
100 to 2800 L h	PK300		
300 to 5000 L h	PK500		
600 to 6000 L h	PK600		
900 to 9000 L h	PK900	Bioprocess Modular 2 in	400 to 12000 L h
1200 to 12000 L h	PK1200		





Chromatography Skid Capabilities to Process Sartobind® Devices

Luer lock

Hose barb or ¼" sanitary flange
























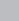








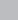

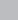

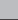

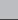


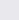
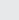

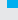
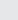
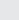


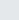
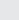
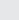


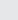
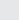
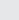
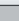
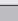
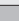
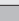
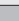
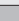
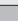
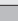
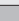
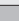

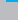
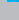
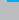
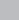

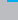
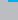
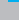
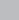










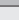
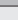
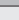
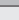
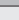
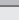
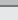
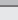
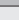
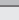

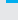
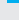
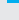
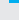

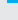
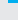
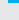
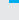










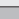
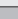
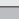
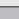
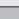
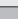
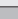
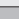
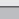

















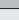
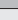
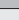
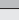
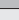
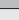
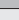
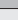
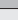
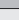
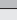
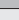












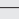
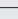
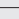
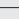
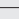
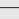
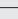
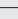
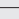
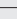
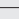

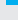
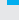
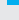
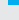
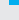
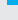
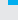
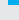
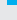
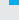















1 ½" sanitary flange

Sartorius System

Cytiva System

System	Single-Use Multi-Use	Flowrate	Detectors		
ÄKTA start	Multi-Use	0.5 to 5 mL min	Flow, pressure, UV, conductivity		
ÄKTA pure 25	Multi-Use	0.001 to 25 mL min	Flow, pressure, UV, conductivity, pH		
ÄKTA avant 25	Multi-Use	0.001 to 25 mL min	Flow, pressure, UV, conductivity, pH		
ÄKTA pure 150	Multi-Use	0.01 to 150 mL min	Flow, pressure, UV, conductivity, pH		
ÄKTA avant 150	Multi-Use	0.01 to 150 mL min	Flow, pressure, UV, conductivity, pH		
ÄKTA pilot	Multi-Use	4.00 to 800 mL min	Flow, pressure, UV, conductivity, pH		
BioProcess Modular ¼ in	Multi-Use	0.02 to 1 L min	Flow, pressure, UV, conductivity, pH		
PKP2	Multi-Use	0.05 to 1.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow		
PKP3	Multi-Use	0.05 to 2.50 L min	Air, pressure, conductivity, temperature, pH, UV, flow		
PK10	Multi-Use	0.17 to 1.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow		
ÄKTA ready with Low Flow Kit	Single-Use	0.05 to 2.92 L min	Flow, pressure, UV, conductivity, pH		
ÄKTA process, 6 mm polypropylene or ¾ in stainless steel	Multi-Use	0.07 to 3.00 L min	Flow, pressure, UV, conductivity, pH electrode (optional)		
PK25	Multi-Use	0.42 to 6.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow		
Sartorius S.U.T. Chrom ¾"	Single-Use	0.17 to 8.33 L min	Air, level, pressure, conductivity, temperature, pH, UV, flow		
ÄKTA ready with High Flow Kit	Single-Use	0.125 to 8.50 L min	Flow, pressure, UV, conductivity, pH		
ÄKTA process, 10 mm polypropylene or ½ in stainless steel	Multi-Use	0.25 to 10.00 L min	Flow, pressure, UV, conductivity, pH electrode (optional)		
Sartorius S.U.T. Chrom ½"	Single-Use	0.33 to 15.00 L min	Air, level, pressure, conductivity, temperature, pH, UV, flow		
PK50	Multi-Use	0.83 to 20.00 L min	Air, pressure, conductivity, temperature, pH, UV, flow		
ÄKTA process, 1 in polypropylene or stainless steel	Multi-Use	0.75 to 30.00 L min	Flow, pressure, UV, conductivity, pH electrode (optional)		
ÄKTA ready XL, ¾" kit	Single-Use	0.75 to 31.67 L min	Flow, pressure, UV, conductivity, pH electrode (optional)		
ÄKTA ready XL, 1" kit	Single-Use	1.50 to 58.33 L min	Flow, pressure, UV, conductivity, pH electrode (optional)		
PK300	Multi-Use	1.66 to 46.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow		
Bioprocess Modular 1.5 in	Multi-Use	3.33 to 83.33 L min	Flow, pressure, UV, conductivity, pH		
PK500	Multi-Use	5.00 to 83.33 L min	Air, pressure, conductivity, temperature, pH, UV, flow		

Size	Pico	Nano
Connection Type		
Membrane	Sartobind® Q Sartobind® S Sartobind® STIC Sartobind® Phenyl	
Bed Height		
Membrane Volume	0.08 mL	1 mL
Recommended Application	Flowthrough B&E	
Recommended Flowrate	10 - 30 MV min	
Minimum Flowrate	0.0008 L min	0.01 L min
Maximum Flowrate	0.0032 L min	0.03 L min

5"	10"	20"	30"	Jumbo	Cassette	Nano	5"	10"	20"	30"	Jumbo	Cassette
												
Sartobind® Q Sartobind® S Sartobind® STIC						Sartobind® Q Sartobind® S Sartobind® Phenyl						
4 mm						8 mm						
75 mL	200 mL	400 mL	600 mL	2500 mL	800 mL*	3 mL	150 mL	400 mL	800 mL	1200 mL	5000 mL	1600 mL*
Flowthrough						Bind and Elute						
10-30 MV min						5 MV min						
0.75 L min	2 L min	4 L min	6 L min	25 L min	8 L min*	0.015 L min	0.75 L min	2 L min	4 L min	6 L min	25 L min	8 L min*
2.25 L min	6 L min	12 L min	18 L min	75 L min	24 L min*	0.015 L min	0.75 L min	2 L min	4 L min	6 L min	25 L min	8 L min*
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												
												

Chromatography Skid Capabilities to Process Columns

				PRC Prepacked							
				10			50			100	
				100	300	600	100	300	600	100	300
<div><div></div>Sartorius System</div> <div><div></div>Cytiva System</div>				Column Diameter (mm)							
				Linear Velocity (cm hr)							
				Flowrate L min							
System	Single-Use Multi-Use	Flowrate	Detectors								
ÄKTA start	Multi-Use	0.5 to 5 mL min	Flow, pressure, UV, conductivity	■	■						
ÄKTA pure 25	Multi-Use	0.001 to 25 mL min	Flow, pressure, UV, conductivity, pH	■	■	■					
ÄKTA avant 25	Multi-Use	0.001 to 25 mL min	Flow, pressure, UV, conductivity, pH	■	■	■					
ÄKTA pure 150	Multi-Use	0.01 to 150 mL min	Flow, pressure, UV, conductivity, pH	■	■	■	■	■		■	
ÄKTA avant 150	Multi-Use	0.01 to 150 mL min	Flow, pressure, UV, conductivity, pH	■	■	■	■	■		■	
ÄKTA pilot	Multi-Use	4.00 to 800 mL min	Flow, pressure, UV, conductivity, pH		■	■	■	■	■	■	■
BioProcess Modular ¼ in	Multi-Use	0.02 to 1 L min	Flow, pressure, UV, conductivity, pH				■	■	■	■	■
PKP2	Multi-Use	0.05 to 1.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow				■	■	■	■	■
PKP3	Multi-Use	0.05 to 2.50 L min	Air, pressure, conductivity, temperature, pH, UV, flow				■	■	■	■	■
PK10	Multi-Use	0.17 to 1.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow						■	■	■
ÄKTA ready with Low Flow Kit	Single-Use	0.05 to 2.92 L min	Flow, pressure, UV, conductivity, pH					■	■	■	■
ÄKTA process, 6 mm polypropylene or ¾ in stainless steel	Multi-Use	0.07 to 3.00 L min	Flow, pressure, UV, conductivity, pH electrode (optional)					■	■	■	■
PK25	Multi-Use	0.42 to 6.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow								
Sartorius S.U.T. Chrom ⅜"	Single-Use	0.17 to 8.33 L min	Air, level, pressure, conductivity, temperature, pH, UV, flow						■		■
ÄKTA ready with High Flow Kit	Single-Use	0.125 to 8.50 L min	Flow, pressure, UV, conductivity, pH								■
ÄKTA process, 10 mm polypropylene or ½ in stainless steel	Multi-Use	0.25 to 10.00 L min	Flow, pressure, UV, conductivity, pH electrode (optional)								■
Sartorius S.U.T. Chrom ½"	Single-Use	0.33 to 15.00 L min	Air, level, pressure, conductivity, temperature, pH, UV, flow								■
PK50	Multi-Use	0.83 to 20.00 L min	Air, pressure, conductivity, temperature, pH, UV, flow								
ÄKTA process, 1 in polypropylene or stainless steel	Multi-Use	0.75 to 30.00 L min	Flow, pressure, UV, conductivity, pH electrode (optional)								
ÄKTA ready XL, ¾" kit	Single-Use	0.75 to 31.67 L min	Flow, pressure, UV, conductivity, pH electrode (optional)								
ÄKTA ready XL, 1" kit	Single-Use	1.50 to 58.33 L min	Flow, pressure, UV, conductivity, pH electrode (optional)								
PK300	Multi-Use	1.66 to 46.66 L min	Air, pressure, conductivity, temperature, pH, UV, flow								
Bioprocess Modular 1.5 in	Multi-Use	3.33 to 83.33 L min	Flow, pressure, UV, conductivity, pH								
PK500	Multi-Use	5.00 to 83.33 L min	Air, pressure, conductivity, temperature, pH, UV, flow								
PK600	Multi-Use	10.00 to 100.00 L min	Air, pressure, conductivity, temperature, pH, UV, flow								
PK900	Multi-Use	15.00 to 150.00 L min	Air, pressure, conductivity, temperature, pH, UV, flow								
Bioprocess Modular 2 in	Multi-Use	6.67 to 200.00 L min	Flow, pressure, UV, conductivity, pH								
PK1200	Multi-Use	20.00 to 200.00 L min	Air, pressure, conductivity, temperature, pH, UV, flow								

				Resolute DM																		
				Resolute Hydraulic DAP																		
Resolute FM				Resolute Linear																		
	200			400			600			800			1000			1500			2000			
600	100	300	600	100	300	600	100	300	600	100	300	600	100	300	600	100	300	600	100	300	600	
0.785	0.524	1.57	3.14	2.09	6.28	12.6	4.71	14.1	28.3	8.38	25.1	50.3	13.1	39.3	78.5	29.5	88.4	177	52.4	157	314	
■	■																					
■	■																					
■	■	■																				
■	■	■		■																		
■	■	■																				
■	■	■		■																		
■	■	■		■																		
■	■	■	■	■	■		■															
■	■	■	■	■	■		■															
■	■	■	■	■	■		■			■												
■	■	■	■	■	■		■			■												
■	■	■	■	■	■		■	■		■												
		■	■	■	■	■	■	■		■			■									
■	■	■	■	■	■		■	■	■	■	■		■			■						
■		■	■	■	■	■	■	■	■	■	■		■			■						
		■	■	■	■	■	■	■	■	■	■		■	■		■			■			
				■	■	■	■	■	■	■	■	■	■	■	■	■	■					
						■		■	■	■	■	■	■	■	■	■	■	■	■	■		
									■		■	■		■	■	■	■	■				
							</															

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