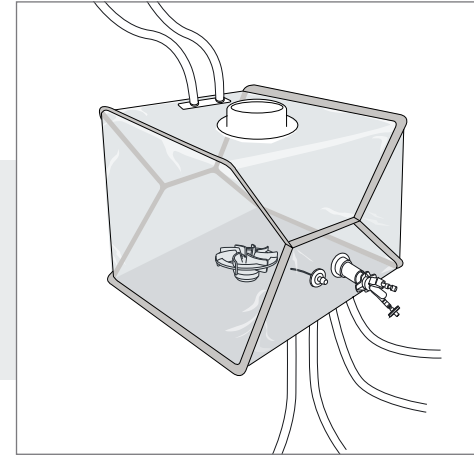


Flexsafe® Pro Mixer Pre-designed Solutions

Drug Substance Purification

Flexsafe® Pro Mixer with pH Sensor and Thermowell



Please visit
www.sartorius.com/flexsafe-pro-mixer
 or contact your local sales representative
 for more information.

Example of a 200 L mixing bag equipped
 with pH sensor & Thermowell

Ordering Information

Bag Volume	Part Number	Units per Box	Drain Valves	Top Port	Inlet	Outlet	Sampling	Addition 1	Addition 2	pH Adjustment
50 L	FMS500077	1	No	Closed 8" Top Port	1/2" ID TuFlux®	1/2" ID TuFlux®	1/4" ID TuFlux®	3/8" ID TuFlux®	NA	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + MPC female + sealing cap
100 L	FMS500079	1	No		Silicone 1,5 m (60")	Silicone 1,5 m (60")	Silicone 0.1 m (4")	Silicone 1.5 m (60")		
200 L	FMS500081	2	No		+ 1/2" Tri-Clamp	+ 1/2" Tri-Clamp	+ Clave needleless sampling port	+ MPC Female + sealing cap	3/8" ID TuFlux®	
400 L	FMS500083	2	No		1 1/2" sanitary flange with gasket, cap & union	1 1/2" sanitary flange with gasket, cap & union			Silicone 1.5 m (60")	
650 L	FMS500085	1	No						+ MPC Female + sealing cap	
1,000 L	FMS500087	1	No							