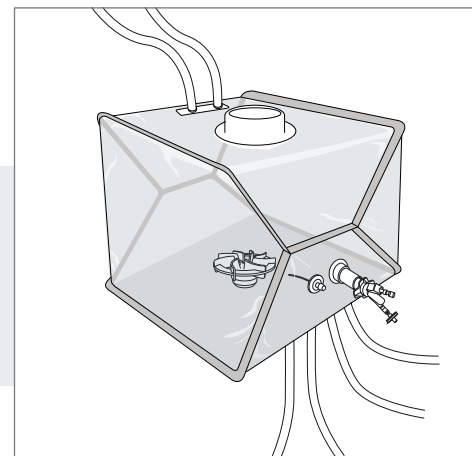


Flexsafe® Pro Mixer Pre-designed Solutions

Buffer Formulation

Intelligent Flexsafe® Pro Mixer with pH Sensor & 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
200 L	FMS500018	2	No	Opened 8" Top Port with gasket, cap & union	1/2" ID TuFlux® Silicone 1,5 m (60") + 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	1/2" ID TuFlux® Silicone 1,5 m (60") + 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	1/4" ID TuFlux® Silicone 0.1 m (4") + Clave needleless sampling port	When drain valves: 1 bottom line: 1/2" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap	When drain valves: 1 bottom line: 1/2" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + MPC female + sealing cap
	FMS500038	2	Yes							
400 L	FMS500019	2	No							
	FMS500039	2	Yes							
650 L	FMS500020	1	No							
	FMS500040	1	Yes							
1,000 L	FMS500021	1	No							
	FMS500041	1	Yes							
1,500 L	FMS500022	1	No	Opened 4" Top Port with gasket, cap & union	1" ID TuFlux® Silicone 3 m (119") + 1" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	1" ID TuFlux® Silicone 3 m (119") + 1" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	1/4" ID TuFlux® Silicone 0.1 m (4") + Clave needleless sampling port	3/4" ID TuFlux® Silicone 1,5 m (60")	3/4" ID TuFlux® Silicone 1,5 m (60")	
2,000 L	FMS500023	1	No							
2,500 L	FMS500024	1	No							
3,000 L	FMS500025	1	No							