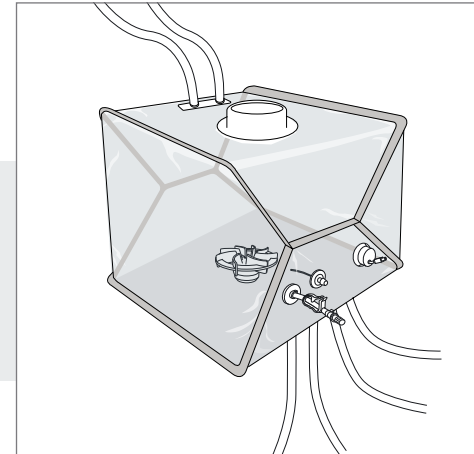


Flexsafe® Pro Mixer Pre-designed Solutions

Buffer Formulation

Intelligent Flexsafe® Pro Mixer with Conductivity sensor & Thermowell



Example of a 200 L mixing bag equipped with conductivity 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	FMS500112	1	No	Opened 8" Top Port with gasket, cap & union	1/2" ID TuFlux®	1/2" ID TuFlux®	1/4" ID TuFlux®	When drain valves: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap	NA	NA
	FMS500138	1	Yes		Silicone 1,5 m (60")	Silicone 1,5 m (60")	Silicone 0.1 m (4")			
100 L	FMS500113	1	No		+ 1/2" Tri-Clamp	+ 1/2" Tri-Clamp	+ Clave needleless sampling port	When no drain valve: 1 bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap		
	FMS500139	1	Yes		1 1/2" sanitary flange with gasket, cap & union	1 1/2" sanitary flange with gasket, cap & union				

Flexsafe® Pro Mixer Pre-designed Solutions

Buffer Formulation

Intelligent Flexsafe® Pro Mixer with Conductivity sensor & Thermowell

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

Ordering Information

Bag Volume	Part Number	Units per Box	Drain Valves	Top Port	Inlet	Outlet	Sampling	Addition 1	Addition 2	pH Adjustment					
200 L	FMS500026	2	No	Opened 8" Top Port with gasket, cap & union	1/2" ID TuFlux®	1/2" ID TuFlux®	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	NA					
	FMS500027	2	Yes		Silicone 1,5 m (60")	Silicone 1,5 m (60")									
400 L	FMS500028	2	No		+ 1/2" Tri-Clamp	+ 1/2" Tri-Clamp									
	FMS500029	2	Yes		1 1/2" sanitary flange with gasket, cap & union	1 1/2" sanitary flange with gasket, cap & union									
650 L	FMS500030	1	No									When no drain valve: 1 bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap	When no drain valve: 1 bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap		
	FMS500031	1	Yes												
1,000 L	FMS500032	1	No												
	FMS500033	1	Yes												
1,500 L	FMS500034	1	No		Opened 4" Top Port with gasket, cap & union	1" ID TuFlux®					1" ID TuFlux®	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	FMS500035	1	No			Silicone 3 m (119")					Silicone 3 m (119")				
2,500 L	FMS500036	1	No	+ 1" Tri-Clamp		+ 1" Tri-Clamp									
3,000 L	FMS500037	1	No	1 1/2" sanitary flange with gasket, cap & union		1 1/2" sanitary flange with gasket, cap & union									