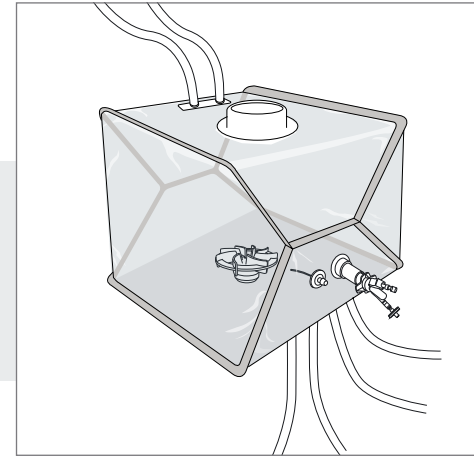


Flexsafe® Pro Mixer Pre-designed Solutions

Media Formulation

Intelligent Flexsafe® Pro Mixer Bag with pH Sensor & Thermowell



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	FMS500004	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	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + MPC female + sealing cap
	FMS500042	1	Yes		Silicone 1,5 m (60")	Silicone 1,5 m (60")	Silicone 0.1 m (4")			
100 L	FMS500005	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		
	FMS500043	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

Media Formulation

Intelligent Flexsafe® Pro Mixer Bag with pH 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	FMS500018	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	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + MPC female + sealing cap				
	FMS500038	2	Yes		Silicone 1,5 m (60")	Silicone 1,5 m (60")								
400 L	FMS500019	2	No		+ 1/2" Tri-Clamp	+ 1/2" Tri-Clamp								
	FMS500039	2	Yes		1 1/2" sanitary flange with gasket, cap & union	1 1/2" sanitary flange with gasket, cap & union								
650 L	FMS500020	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	
	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®	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	FMS500023	1	No		Silicone 3 m (119")	Silicone 3 m (119")								
2,500 L	FMS500024	1	No		+ 1" Tri-Clamp	+ 1" Tri-Clamp				+ 3/4" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ 3/4" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union			
3,000 L	FMS500025	1	No		1 1/2" sanitary flange with gasket, cap & union	1 1/2" sanitary flange with gasket, cap & union								