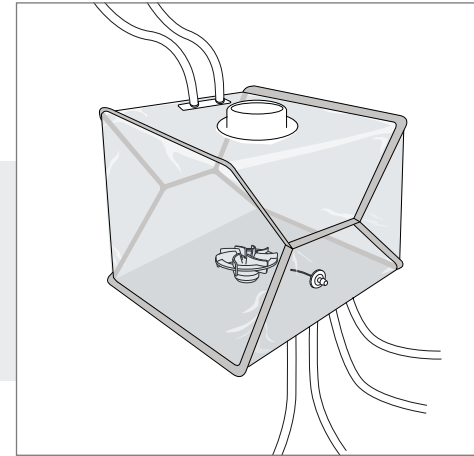


Flexsafe® Pro Mixer Pre-designed Solutions

Media Formulation

Intelligent Flexsafe® Pro Mixer Bag with Thermowell



Example of a 200 L mixing bag equipped with 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	FMS500044	1	Yes	Opened 8" Top Port with gasket, cap & union	1/2" ID TuFlux®	1/2" ID TuFlux®	1/4" ID TuFlux®	When drain valves: Top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap	When drain valves: Top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + MPC female + sealing cap
	FMS500045	1	No		Silicone 1,5 m (60")	Silicone 1,5 m (60")	Silicone 0.1 m (4")			
100 L	FMS500046	1	Yes		+ 1/2" Tri-Clamp	+ 1/2" Tri-Clamp	+ Clave needleless sampling port	When no drain valve: Bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap	When no drain valve: Bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap	
	FMS500047	1	No		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 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	FMS500048	2	Yes	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: Bottom line: 1/2" ID TuFlux® Silicone 1.5 m (60") + MPC Female + sealing cap	When drain valves: 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
	FMS500049	2	No							
400 L	FMS500050	2	Yes							
	FMS500051	2	No							
650 L	FMS500052	1	Yes							
	FMS500053	1	No							
1,000 L	FMS500054	1	Yes							
	FMS500055	1	No							
1,500 L	FMS500056	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" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	3/4" ID TuFlux® Silicone 1,5 m (60") + 3/4" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	
2,000 L	FMS500057	1	No							
2,500 L	FMS500058	1	No							
3,000 L	FMS500059	1	No							