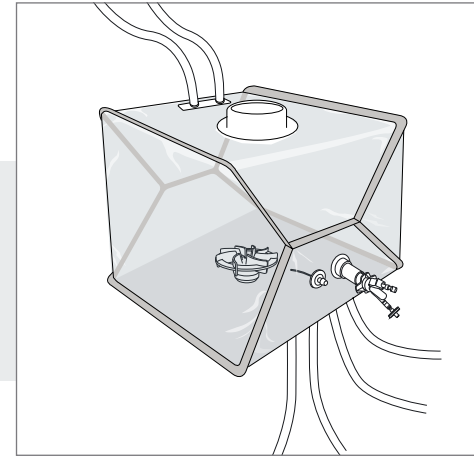


Flexsafe® Pro Mixer Pre-designed Solutions

Drug Product Formulation

Flexsafe® Pro Mixer with pH Sensor and 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	FMS500114	1	Yes	Closed 8" Top Port	1/2" TuFlux®	1/2" ID TuFlux®	1/4" ID TuFlux®	When drain valves: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + 3/8" Opta® SFT female	NA	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + 1/4" Opta® SFT female
	FMS500115	1	No		Silicone 1 m (40")	Silicone 1 m (40")	Silicone 0.1 m (4")			
100 L	FMS500116	1	Yes		+ 1/2" ID TuFlux®	+ 1/2" ID TuFlux®	+ Clave needleless sampling port	When no drain valve: 1 bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + 3/8" Opta® SFT female		
	FMS500117	1	No		TPE 0.5 m (20") + 1/2" Opta® SFT female	TPE 0.5 m (20") + 1/2" Opta® SFT male				

Flexsafe® Pro Mixer Pre-designed Solutions

Drug Product Formulation

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.

Ordering Information

Bag Volume	Part Number	Units per Box	Drain Valves	Top Port	Inlet	Outlet	Sampling	Addition 1	Addition 2	pH Adjustment
200 L	FMS500118	2	Yes	Closed 8" Top Port	1/2" TuFlux® Silicone 1 m (40")	1/2" ID TuFlux® Silicone 1 m (40")	1/4" ID TuFlux® Silicone 0.1 m (4")	When drain valves: 1 bottom line: 1/2" ID TuFlux® Silicone 1.5 m (60") + 1/2" Opta® SFT female	When drain valves: 1 bottom line: 1/2" ID TuFlux® Silicone 1.5 m (60") + 1/2" Opta® SFT female	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + 1/4" Opta® SFT female
	FMS500119	2	No							
400 L	FMS500120	2	Yes		+ 1/2" ID TuFlux® TPE 0.5 m (20") + 1/2" Opta® SFT female	+ 1/2" ID TuFlux® TPE 0.5 m (20") + 1/2" Opta® SFT male	+ Clave needleless sampling port	When no drain valve: 1 bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + 3/8" Opta® SFT female	When no drain valve: 1 bottom line: 3/8" ID TuFlux® Silicone 1.5 m (60") + 3/8" Opta® SFT female	
	FMS500121	2	No							
650 L	FMS500122	1	Yes							
	FMS500123	1	No							
1,000 L	FMS500124	1	Yes							
	FMS500125	1	No							