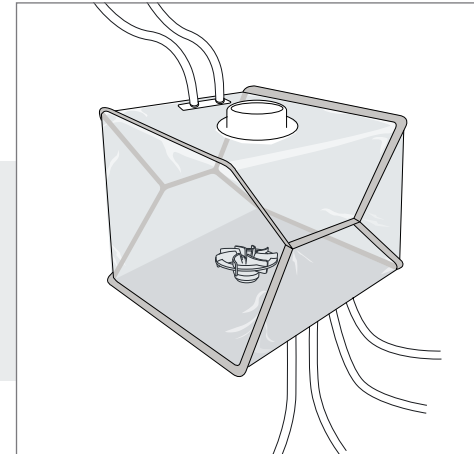


Flexsafe® Pro Mixer Pre-designed Solutions

Drug Product Formulation

Flexsafe® Pro Mixer Without Single-use Sensor and With Sterile Connections and Disconnections



Example of a 200 L mixing bag without single-use sensor and with sterile connections and disconnections

Ordering Information

Bag Volume	Part Number	Units per Box	Drain Valves	Top Port	Inlet	Outlet	Sampling	Addition 1	Addition 2	pH Adjustment
50 L	FMS500126	1	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 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + 3/8" Opta® SFT female	When drain valves: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + 3/8" Opta® SFT female	NA
	FMS500127	1	No							
100 L	FMS500128	1	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	
	FMS500129	1	No							

Flexsafe® Pro Mixer Pre-designed Solutions

Drug Product Formulation

Flexsafe® Pro Mixer Without Single-use Sensor and With Sterile Connections and Disconnections

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	FMS500130	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	NA
	FMS500131	2	No							
400 L	FMS500132	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	
	FMS500133	2	No							
650 L	FMS500134	1	Yes							
	FMS500135	1	No							
1,000 L	FMS500136	1	Yes							
	FMS500137	1	No							