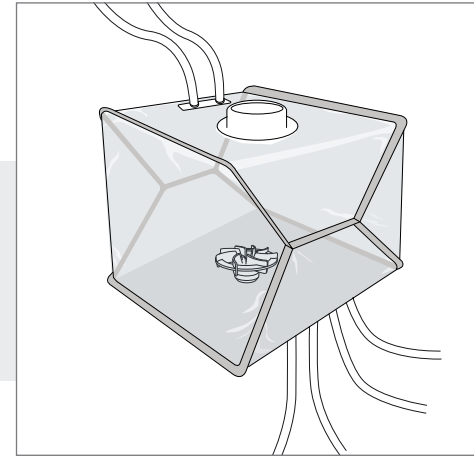


# Flexsafe® Pro Mixer Pre-designed Solutions

## Drug Product Formulation

Flexsafe® Pro Mixer Without Single-use Sensor and With Aseptic Connections and Disconnections under ISO Laminar Air Flow



Example of a 200 L mixing bag without single-use sensor and with aseptic connections and disconnections under ISO laminar air flow

### Ordering Information

Bag Volume	Part Number	Units per Box	Drain Valves	Top Port	Inlet	Outlet	Sampling	Addition 1	Addition 2	pH Adjustment
50 L	FMS500100	1	Yes	Closed 8" Top Port	1/2" ID TuFlux® Silicone 1,5 m (60")	1/2" ID TuFlux® Silicone 1,5 m (60")	1/4" ID TuFlux® Silicone 0.1 m (4")	When drain valves: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap	When drain valves: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap	NA
	FMS500101	1	No							
100 L	FMS500102	1	Yes		+ 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ Clave needleless sampling port	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	
	FMS500103	1	No							

# Flexsafe® Pro Mixer Pre-designed Solutions

## Drug Product Formulation

Flexsafe® Pro Mixer Without Single-use Sensor and With Aseptic Connections and Disconnections under ISO Laminar Air Flow

Please visit  
[www.sartorius.com/flexsafe-pro-mixer](http://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	FMS500104	2	Yes	Closed 8" Top Port	1/2" ID TuFlux® Silicone 1,5 m (60")	1/2" ID TuFlux® Silicone 1,5 m (60")	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
	FMS500105	2	No							
400 L	FMS500106	2	Yes		+ 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union				
	FMS500107	2	No							
650 L	FMS500108	1	Yes							
	FMS500109	1	No							
1,000 L	FMS500110	1	Yes					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	
	FMS500111	1	No							