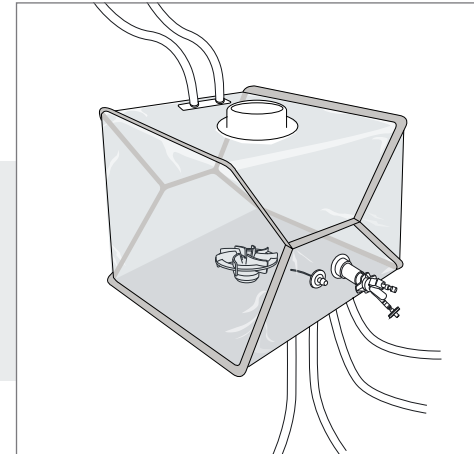


Flexsafe® Pro Mixer Pre-designed Solutions

Downstream Intermediates

Flexsafe® Pro Mixer with pH Sensor and Thermowell for Low pH Viral Inactivation and Neutralisation



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	FMS500076	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	NA	2 top lines: 1/4" ID TuFlux® Silicone 0.1 m (4") + MPC female + sealing cap
	FMS500077	1	No							
100 L	FMS500078	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		
	FMS500079	1	No							

Flexsafe® Pro Mixer Pre-designed Solutions

Downstream Intermediates

Flexsafe® Pro Mixer with pH Sensor and Thermowell for
Low pH Viral Inactivation and Neutralisation

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	FMS500080	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")	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
	FMS500081	2	No							
400 L	FMS500082	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	+ 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	
	FMS500083	2	No							
650 L	FMS500084	1	Yes							
	FMS500085	1	No							
1,000 L	FMS500086	1	Yes							
	FMS500087	1	No							