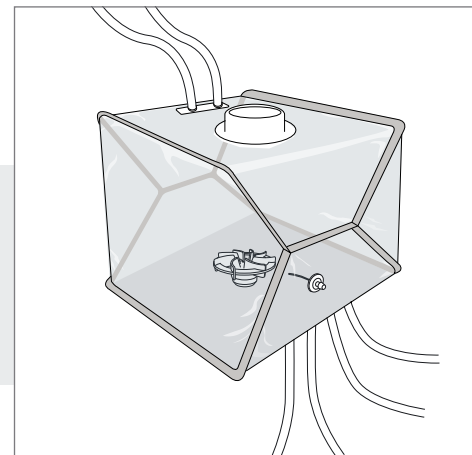


# Flexsafe® Pro Mixer Pre-designed Solutions

## Downstream Intermediates

Flexsafe® Pro Mixer with Thermowell for Post Filtration



Please visit  
[www.sartorius.com/flexsafe-pro-mixer](http://www.sartorius.com/flexsafe-pro-mixer)  
 or contact your local sales representative  
 for more information.

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	FMS500094	1	No	Closed 8" Top Port	1/2" ID TuFlux®	1/2" ID TuFlux®	1/4" ID TuFlux®	3/8" ID TuFlux®	3/8" ID TuFlux®	NA
100 L	FMS500095	1	No		Silicone 1 m (40")	Silicone 1 m (40")	Silicone 0.1 m (4")	Silicone 1.5 m (60")	Silicone 1.5 m (60")	
200 L	FMS500096	2	No		+ 1/2" ID TuFlux® TPE	+ 1/2" ID TuFlux® TPE	+ Clave needleless sampling port	+ MPC Female	+ MPC Female	
400 L	FMS500097	2	No		0.5 m (20")	0.5 m (20")		+ sealing cap	+ sealing cap	
650 L	FMS500098	1	No		+ 1/2" Opta® SFT female	+ 1/2" Tri-Clamp				
1,000 L	FMS500099	1	No			1 1/2" sanitary flange with gasket, cap & union				