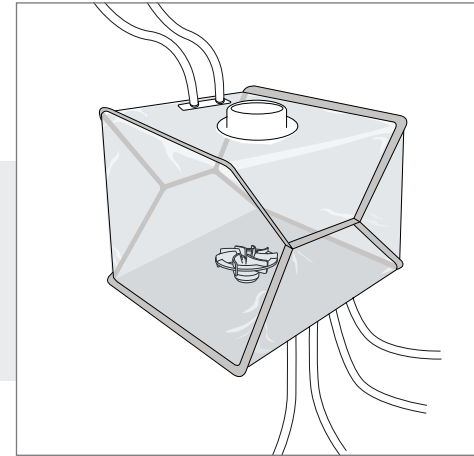


Flexsafe® Pro Mixer Pre-designed Solutions

Media Formulation

Flexsafe® Pro Mixer Without Single-use Sensor



Please visit
www.sartorius.com/flexsafe-pro-mixer
 or contact your local sales representative
 for more information.

Example of a 200 L mixing bag
 without single-use sensor

Ordering Information

Bag Volume	Part Number	Units per Box	Drain Valves	Top Port	Inlet	Outlet	Sampling	Addition 1	Addition 2	pH Adjustment
50 L	FMS500060	1	Yes	Opened 8" Top Port with gasket, cap & union	1/2" ID TuFlux® Silicone 1,5 m (60") + 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	1/2" ID TuFlux® Silicone 1,5 m (60") + 1/2" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	1/4" ID TuFlux® Silicone 0.1 m (4") + Clave needleless sampling port	NA	NA	NA
	FMS500061	1	No							
100 L	FMS500062	1	Yes							
	FMS500063	1	No							
200 L	FMS500064	2	Yes							
	FMS500065	2	No							
400 L	FMS500066	2	Yes							
	FMS500067	2	No							
650 L	FMS500068	1	Yes							
	FMS500069	1	No							
1,000 L	FMS500070	1	Yes							
	FMS500071	1	No							
1,500 L	FMS500072	1	No	NA	1" ID TuFlux® Silicone 3 m (119")	1" ID TuFlux® Silicone 3 m (119")	1/4" ID TuFlux® Silicone 0.1 m (4")	3/4" ID TuFlux® Silicone 1,5 m (60")	3/4" ID TuFlux® Silicone 1,5 m (60")	
2,000 L	FMS500073	1	No							
2,500 L	FMS500074	1	No							
3,000 L	FMS500075	1	No		+ 1" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ 1" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ Clave needleless sampling port	+ 3/4" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	+ 3/4" Tri-Clamp 1 1/2" sanitary flange with gasket, cap & union	