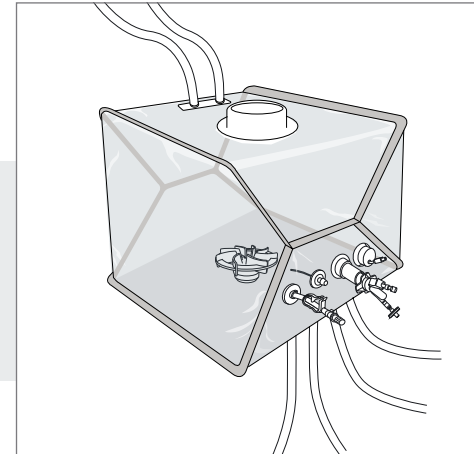


Flexsafe® Pro Mixer Pre-designed Solutions

Buffer Formulation

Intelligent Flexsafe® Pro Mixer Bag with pH Sensor, Conductivity sensor & Thermowell



Example of a 200 L mixing bag equipped with pH sensor, conductivity 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 | FMS500014 | 1 | No | Opened 8" Top Port with gasket, cap & union | 1/2" ID TuFlux® | 1/2" ID TuFlux® | 1/4" ID TuFlux® | When drain valves: NA | When drain valves: NA | 2 top lines: 1/4" ID TuFlux® |
| | FMS500015 | 1 | Yes | | Silicone 1,5 m (60") | Silicone 1,5 m (60") | Silicone 0.1 m (4") | | | |
| 100 L | FMS500017 | 1 | No | | + 1/2" Tri-Clamp | + 1/2" Tri-Clamp | + Clave needleless sampling port | When no drain valve: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap | When no drain valve: 1 top line: 3/8" ID TuFlux® Silicone 0.15 m (6") + MPC Female + sealing cap | + MPC female + sealing cap |
| | FMS500016 | 1 | Yes | | 1 1/2" sanitary flange with gasket, cap & union | 1 1/2" sanitary flange with gasket, cap & union | | | | |

Flexsafe® Pro Mixer Pre-designed Solutions

Buffer Formulation

Intelligent Flexsafe® Pro Mixer Bag with pH Sensor,
Conductivity sensor & Thermowell

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 | FMS500006 | 2 | No | Opened 8" Top Port with gasket, cap & union | 1/2" ID TuFlux® | 1/2" ID TuFlux® | 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 no drain valve: 1 bottom line: 3/8" 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 When no drain valve: 1 bottom line: 3/8" 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 |
| | FMS500007 | 2 | Yes | | 1/2" ID TuFlux® Silicone 1,5 m (60") | Silicone 1,5 m (60") | | | | |
| 400 L | FMS500008 | 2 | No | | + 1/2" Tri-Clamp | + 1/2" Tri-Clamp | | | | |
| | FMS500009 | 2 | Yes | | 1 1/2" sanitary flange with gasket, cap & union | 1 1/2" sanitary flange with gasket, cap & union | | | | |
| 650 L | FMS500010 | 1 | No | | | | | | | |
| | FMS500011 | 1 | Yes | | | | | | | |
| 1,000 L | FMS500012 | 1 | No | | | | | | | |
| | FMS500013 | 1 | Yes | | | | | | | |
| 1,500 L | FMS500000 | 1 | No | | Opened 4" Top Port with gasket, cap & union | 1" ID TuFlux® | | | | |
| 2,000 L | FMS500001 | 1 | No | 1" ID TuFlux® Silicone 3 m (119") | | Silicone 3 m (119") | | | | |
| 2,500 L | FMS500002 | 1 | No | + 1" Tri-Clamp | | + 1" Tri-Clamp | | | | |
| 3,000 L | FMS500003 | 1 | No | 1 1/2" sanitary flange with gasket, cap & union | | 1 1/2" sanitary flange with gasket, cap & union | | | | |