

Flexsafe® Bags for Liquid Storage, Shipping, and Mixing

Made-to-stock for rapid,
flexible integration across
workflows



Product Information

Our Flexsafe® bags for storage and mixing are available in selected made-to-stock configurations with short lead times. These simple, fully interconnectable designs are offered in a broad range of volumes to support scalability and easy integration into most process development and manufacturing steps.

Features and Benefits

- Reliable material performance: Flexsafe® S80 film provides excellent biocompatibility, chemical resistance, and robustness.
- Sustainable workflows: S80 film is ISCC PLUS certified and contains 71% biocircular content.
- End-to-end scalability: 20 mL – 3,000 L formats with versatile designs for process integration across modalities.
- Process-ready solutions: Full validation and extractables packages, ASTM D4169 shipping qualification, and mixing performance supported by CFD studies.

Relevant applications

- Research and development
- GMP and non-GMP applications
- Biopharmaceutical manufacturing

Relevant process steps

Applicable across upstream and downstream operations, from small-scale research and development to clinical and commercial manufacturing:

- Buffer | media preparation
 - Liquid–liquid or powder–liquid mixing
 - Storage and intermediate hold
 - Drug substance handling
 - Final formulation
 - Shipping
-

Overview

Selected Flexsafe® 2D, 3D, and mixing bags are available in made-to-stock configurations to help standardize and accelerate single-use workflows. Designed for routine storage, mixing, and shipping needs, the offering supports flexible integration across applications, with robust film performance and production aligned with sustainable manufacturing initiatives.

Flexsafe® S80 film performance

Flexsafe® S80 is a 400 µm co-extruded PE | EVOH | PE film designed to balance strength, flexibility, and barrier performance across storage, mixing, and shipping applications.

- Robust film construction combines high puncture resistance, elongation at break, and a large sealing window to support bag integrity across storage, mixing, and shipping workflows.
- Excellent biocompatibility helps maintain media quality during storage and supports consistent cell culture performance.
- Superior flexibility supports bag handling, installation, deployment in containers, and use in shipping and mixing applications.
- High gas and water barrier properties help preserve media, buffers, and drug substances during long-term storage.
- Inert PE contact layer supports use with non-buffered solutions.
- Low extractables profile supports critical drug substance and drug product process steps.
- ISCC PLUS-certified sites and mass-balance traceability support more sustainable production with 71% biocircular content.



Technical Specifications

Film validation and quality assurance

Feature	Test method
Film and connection robustness	- ASTM D882 Standard Test Method for Tensile Properties of Thin Plastic Sheeting - ISO 527-3 Plastics – Determination of Tensile Properties – Part 3: Test Conditions for Films and Sheets - ASTM D3787 Standard Test Method for Bursting Strength of Textiles – Constant-Rate-of-Traverse (CRT) Ball Burst Test (resistance to puncture) - ASTM F392 Standard Practice for Conditioning Flexible Barrier Materials for Flex Durability Internal standardized methods are used for connection testing and functional qualification under representative process conditions
Film gas transmission	- ASTM D3985 Standard Test Method for Oxygen Gas Transmission Rate Through Plastic Film and Sheeting Using a Coulometric Sensor - ISO 15106-3 Plastics – Film and Sheeting – Determination of Water Vapour Transmission Rate – Part 3: Coulometric Sensor Method - ASTM F2476 Standard Test Method for the Determination of Carbon Dioxide Gas Transmission Rate (CO₂TR) Through Barrier Materials Using an Infrared Detector
Biocompatibility	- USP <87> Biological Reactivity Tests, In Vitro - ASTM E3231 Standard Guide for Cell Culture Growth Assessment of Single-Use Material - USP <88> Biological Reactivity Tests, In Vivo - ISO 10993-5 Biological Evaluation of Medical Devices – Part 5: Tests for In Vitro Cytotoxicity
Chemical compatibility	ASTM D543-06 Standard Practices for Evaluating the Resistance of Plastics to Chemical Reagents
Material characterization: PE contact layer	EP 3.1.5 Polyethylene with Additives for Containers for Preparations for Parenteral Use and for Ophthalmic Preparations
Material characterization: PE bag port	EP 3.1.5 Polyethylene with Additives for Containers for Preparations for Parenteral Use and for Ophthalmic Preparations
Purity, extractables, and leachables	- USP <665> Plastic Components and Systems Used to Manufacture Pharmaceutical Drug Products and Biopharmaceutical Drug Substances and Products (extractable data) - EP 5.2.8 Minimising the Risk of Transmitting Animal Spongiform Encephalopathy Agents via Human and Veterinary Medicinal Products (TSE BSE control)
Cleanliness, particles, and sterilization	- USP <788> Particulate Matter in Injections - EP 2.9.19 Particulate Contamination: Sub-Visible Particles Visible particles: Representative products have been analyzed for particles ≥ 100 µm according to internally developed and validated methods. Results are monitored. - USP <85> Bacterial Endotoxins Test - EP 2.6.14 Bacterial Endotoxins - ISO 11737 Sterilization of Health Care Products – Microbiological Methods - ISO 11137 Sterilization of Health Care Products – Radiation - ISO 14644 Cleanrooms and Associated Controlled Environments - Gamma radiation dose mapping

Flexsafe® 2D Bags

Ordering number	Product	Description		
		Inlet	Outlet	Sampling line
FLS130122	Flexsafe® 2D 50 mL	TPE: ¼" × ⅜", 500 mm MPC female + cap	TPE: ¼" × ⅜", 500 mm MPC male + cap	N/A
FLS130017	Flexsafe® 2D 500 mL	TPE: ¼" × ⅜", 500 mm MPC female + cap	TPE: ¼" × ⅜", 500 mm MPC male + cap	N/A
FLS130129	Flexsafe® 2D 1 L	TPE: ¼" × ⅜", 500 mm MPC female + cap	TPE: ¼" × ⅜", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅜", 100 mm Luer + Clave® needleless sampling port
FLS130145	Flexsafe® 2D 3 L	TPE: ⅜" × ⅝", 500 mm MPC female + cap Clipster® in separate pouch	TPE: ⅜" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅜", 100 mm Luer + Clave® needleless sampling port
FLS130039	Flexsafe® 2D 5 L	TPE: ⅜" × ⅝", 500 mm MPC female + cap Clipster® in separate pouch	TPE: ⅜" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅜", 100 mm Luer + Clave® needleless sampling port
FLS130040	Flexsafe® 2D 10 L	TPE: ⅜" × ⅝", 500 mm MPC female + cap Clipster® in separate pouch	TPE: ⅜" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅜", 100 mm Luer + Clave® needleless sampling port
FLS130041	Flexsafe® 2D 20 L	TPE: ⅜" × ⅝", 500 mm MPC female + cap Clipster® in separate pouch	TPE: ⅜" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅜", 100 mm Luer + Clave® needleless sampling port
FLS130146	Flexsafe® 2D 50 L	TPE: ⅜" × ⅝", 500 mm MPC female + cap Clipster® in separate pouch	TPE: ⅜" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅜", 100 mm Luer + Clave® needleless sampling port



Flexsafe® 3D Bags

Ordering number	Product	Description		
		Inlet	Outlet	Sampling line
FIS320893	Flexsafe® 3D for Palletank® 50 L	Tuflux® SIL (Pt): ½" x ¾", 2,000 mm TPE: ½" x ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ½" x ¾", 2,000 mm TPE: ½" x ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS320894	Flexsafe® 3D for Palletank® 50 L	Tuflux® SIL (Pt): ½" x ¾", 2,000 mm TPE: ½" x ¾", 500 mm AseptiQuik® G	Tuflux® SIL (Pt): ½" x ¾", 2,000 mm TPE: ½" x ¾", 500 mm AseptiQuik® G	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS320895	Flexsafe® 3D for Palletank® 50 L	Tuflux® SIL (Pt): ½" x ¾", 2,000 mm TPE: ½" x ¾", 500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 2,000 mm TPE: ½" x ¾", 500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS501317	Flexsafe® 3D for Palletank® 100 L	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm, with Clipster® L AseptiQuik® G	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm, with Clipster® L AseptiQuik® G	N/A
FIS126489	Flexsafe® 3D for Palletank® 100 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS126730	Flexsafe® 3D for Palletank® 100 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm Sartopore® 2 0.2 µm, Size 8	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS500482	Flexsafe® 3D for Palletank® 200 L	Tuflux® SIL (Pt): ⅜" x ⅝", 2,500 mm TPE: ⅜" x ⅝", 500 mm AseptiQuik® G	Tuflux® SIL (Pt): ⅜" x ⅝", 2,500 mm TPE: ⅜" x ⅝", 500 mm AseptiQuik® G	N/A
FIS126252	Flexsafe® 3D for Palletank® 200 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm Sartopore® 2 0.2 µm, Size 9	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS126256	Flexsafe® 3D for Palletank® 200 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS126490	Flexsafe® 3D for Palletank® 200 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS500128	Flexsafe® 3D for Palletank® 500 L	Tuflux® SIL (Pt): ¾" x 1⅝", 300 mm AseptiQuik® G	Tuflux® SIL (Pt): ¾" x 1⅝", 300 mm AseptiQuik® G	N/A
FIS126253	Flexsafe® 3D for Palletank® 500 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm Sartopore® 2 0.2 µm, Size 0	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS126257	Flexsafe® 3D for Palletank® 500 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS126491	Flexsafe® 3D for Palletank® 500 L	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port
FIS126258	Flexsafe® 3D for Palletank® 1,000 L (Flexel® format)	Tuflux® SIL (Pt): ½" x ¾", 1,000 mm TPE: ½" x ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅝", 100 mm Luer + Clave® needleless sampling port

FIS126246	Flexsafe® 3D for Palletank® 1,000 L (Flexel® format)	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ½" × ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ½" × ¾", 500 mm MPX male + cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS302338	Flexsafe® 3D for Palletank® 1,000 L (LV format)	TPE: ½" × ¾", 1,500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 2,500 mm TPE: ½" × ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS501833	Flexsafe® 3D for Palletank® 1,000 L (LV format)	Inlet 1 Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", with union, gasket, and cap Inlet 2 Tuflux® SIL (Pt): ½" × ¾", 1,500 mm AsepticQuik® G	Outlet 1 Tuflux® SIL (Pt): ½" × ¾", 1,500 mm AsepticQuik® G Outlet 2 Tuflux® SIL (Pt): ½" × ¾", 1,500 mm AsepticQuik® G	N/A
FIS302339	Flexsafe® 3D for Palletank® 1,500 L (LV format)	TPE: ½" × ¾", 1,500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 2,500 mm TPE: ½" × ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS500947	Flexsafe® 3D for Palletank® 1,500 L (LV format)	Tuflux® SIL (Pt): ¾" × 1⅝", 2,500 mm TPE: ¾" × 1", 500 mm AsepticQuik® G	Tuflux® SIL (Pt): ¾" × 1⅝", 2,500 mm TPE: ¾" × 1", 500 mm AsepticQuik® G	N/A
FIS302340	Flexsafe® 3D for Palletank® 2,000 L (LV format)	TPE: ½" × ¾", 1,500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 2,500 mm TPE: ½" × ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS302341	Flexsafe® 3D for Palletank® 2,500 L (LV format)	TPE: ½" × ¾", 1,500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 2,500 mm TPE: ½" × ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS302342	Flexsafe® 3D for Palletank® 3,000 L (LV format)	TPE: ½" × ¾", 1,500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 2,500 mm TPE: ½" × ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port

Flexsafe® 3D Bags for Drums

Ordering number	Product	Description		
		Inlet	Outlet	Sampling line
FIS127903	Flexsafe® Drum 50 L (top drain)	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ½" × ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ½" × ¾", 500 mm MPX male + double female MPX cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS127876	Flexsafe® Drum 100 L (top drain)	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ½" × ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ⅝" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port
FIS127877	Flexsafe® Drum 200 L (top drain)	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ½" × ¾", 500 mm MPX female + cap	Tuflux® SIL (Pt): ½" × ¾", 1,000 mm TPE: ⅝" × ⅝", 500 mm MPC male + cap	Tuflux® SIL (Pt): ¼" × ⅝", 100 mm Luer + Clave® needleless sampling port

Flexsafe® Bags for Pro Mixer

Ordering number	Product	Description			
		Top port	Inlet	Outlet	Sampling line
FMS319774	Flexsafe® for Pro Mixer 5 L (powder)	Open 4" top port with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS319775	Flexsafe® for Pro Mixer 10 L (powder)	Open 4" top port with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS319776	Flexsafe® for Pro Mixer 20 L (powder)	Open 4" top port with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" × 5/8", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" × 5/8", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS319777	Flexsafe® for Pro Mixer 5 L	No top port	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS319778	Flexsafe® for Pro Mixer 10 L	No top port	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS319779	Flexsafe® for Pro Mixer 20 L	No top port	Tuflux® SIL (Pt): ⅜" × 5/8", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" × 5/8", 1,000 mm Tri-Clamp, 1½", with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS500061	Flexsafe® for Pro Mixer 50 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS500063	Flexsafe® for Pro Mixer 100 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS500065	Flexsafe® for Pro Mixer 200 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port
FMS500067	Flexsafe® for Pro Mixer 400 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" × ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¼" × 7/16", 100 mm Luer + Clave® needleless sampling port

FMS500069	Flexsafe® for Pro Mixer 650 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅞", 100 mm Luer + Clave® needleless sampling port
FMS500071	Flexsafe® for Pro Mixer 1,000 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¼" x ⅞", 100 mm Luer + Clave® needleless sampling port

Flexsafe® Bags for Magnetic Mixer

Ordering number	Product	Description				
		Top port	Inlet	Outlet	Addition	Sampling line
FMS301053	Flexsafe® for Magnetic Mixer 50 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301054	Flexsafe® for Magnetic Mixer 100 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301055	Flexsafe® for Magnetic Mixer 200 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301056	Flexsafe® for Magnetic Mixer 400 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301057	Flexsafe® for Magnetic Mixer 650 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301058	Flexsafe® for Magnetic Mixer 1,000 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301059	Flexsafe® for Magnetic Mixer 1,500 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ⅜" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port

Flexsafe® Bags for LevMixer®

Ordering number	Product	Description				
		Top port	Inlet	Outlet	Addition	Sampling line
FMS301047	Flexsafe® for LevMixer® 50 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¾" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301048	Flexsafe® for LevMixer® 100 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¾" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301049	Flexsafe® for LevMixer® 200 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¾" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301050	Flexsafe® for LevMixer® 400 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¾" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301051	Flexsafe® for LevMixer® 650 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¾" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port
FMS301052	Flexsafe® for LevMixer® 1,000 L	Open 8" top port with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ½" x ¾", 1,500 mm Tri-Clamp, 1½", sanitary flange, with union, gasket, and cap	Tuflux® SIL (Pt): ¾" x ⅝", 1,500 mm MPC female + cap	Tuflux® SIL (Pt): ¼" x ⅜", 100 mm Luer + Clave® needleless sampling port

Ordering Information

Ordering number	Product	Units per box	Minimum order quantity (units)
FLS130122	Flexsafe® 2D 50 mL	10	10
FLS130017	Flexsafe® 2D 500 mL	10	10
FLS130129	Flexsafe® 2D 1 L	10	10
FLS130145	Flexsafe® 2D 3 L	16	16
FLS130039	Flexsafe® 2D 5 L	16	16
FLS130040	Flexsafe® 2D 10 L	8	8
FLS130041	Flexsafe® 2D 20 L	8	8
FLS130146	Flexsafe® 2D 50 L	8	8
FIS320893	Flexsafe® 3D for Palletank® 50 L	6	6
FIS320894	Flexsafe® 3D for Palletank® 50 L	6	6
FIS320895	Flexsafe® 3D for Palletank® 50 L	6	6
FIS501317	Flexsafe® 3D for Palletank® 100 L	2	2
FIS126489	Flexsafe® 3D for Palletank® 100 L	3	3
FIS126730	Flexsafe® 3D for Palletank® 100 L	2	2
FIS500482	Flexsafe® 3D for Palletank® 200 L	2	2
FIS126252	Flexsafe® 3D for Palletank® 200 L	2	2
FIS126256	Flexsafe® 3D for Palletank® 200 L	3	3
FIS126490	Flexsafe® 3D for Palletank® 200 L	3	3
FIS500128	Flexsafe® 3D for Palletank® 500 L	2	2
FIS126253	Flexsafe® 3D for Palletank® 500 L	2	2
FIS126257	Flexsafe® 3D for Palletank® 500 L	3	3
FIS126491	Flexsafe® 3D for Palletank® 500 L	3	3
FIS126258	Flexsafe® 3D for Palletank® 1,000 L (Flexel® format)	3	3
FIS126246	Flexsafe® 3D for Palletank® 1,000 L (Flexel® format)	3	3
FIS302338	Flexsafe® 3D for Palletank® 1,000 L (LV format)	2	2
FIS501833	Flexsafe® 3D for Palletank® 1,000 L (LV format)	1	1
FIS302339	Flexsafe® 3D for Palletank® 1,500 L (LV format)	2	2
FIS500947	Flexsafe® 3D for Palletank® 1,500 L (LV format)	1	1
FIS302340	Flexsafe® 3D for Palletank® 2,000 L (LV format)	2	2
FIS302341	Flexsafe® 3D for Palletank® 2,500 L (LV format)	1	1
FIS302342	Flexsafe® 3D for Palletank® 3,000 L (LV format)	1	1
FIS127903	Flexsafe® Drum 50 L (top drain)	4	4
FIS127876	Flexsafe® Drum 100 L (top drain)	4	4
FIS127877	Flexsafe® Drum 200 L (top drain)	4	4

Ordering number	Product	Units per box	Minimum order quantity (units)
FMS319774	Flexsafe® for Pro Mixer 5 L (powder)	1	1
FMS319775	Flexsafe® for Pro Mixer 10 L (powder)	1	1
FMS319776	Flexsafe® for Pro Mixer 20 L (powder)	1	1
FMS319777	Flexsafe® for Pro Mixer 5 L	1	1
FMS319778	Flexsafe® for Pro Mixer 10 L	1	1
FMS319779	Flexsafe® for Pro Mixer 20 L	1	1
FMS500061	Flexsafe® for Pro Mixer 50 L	1	1
FMS500063	Flexsafe® for Pro Mixer 100 L	1	1
FMS500065	Flexsafe® for Pro Mixer 200 L	2	2
FMS500067	Flexsafe® for Pro Mixer 400 L	2	2
FMS500069	Flexsafe® for Pro Mixer 650 L	1	1
FMS500071	Flexsafe® for Pro Mixer 1,000 L	1	1
FMS301053	Flexsafe® for Magnetic Mixer 50 L	1	1
FMS301054	Flexsafe® for Magnetic Mixer 100 L	1	1
FMS301055	Flexsafe® for Magnetic Mixer 200 L	2	2
FMS301056	Flexsafe® for Magnetic Mixer 400 L	2	2
FMS301057	Flexsafe® for Magnetic Mixer 650 L	1	1
FMS301058	Flexsafe® for Magnetic Mixer 1,000 L	1	1
FMS301059	Flexsafe® for Magnetic Mixer 1,500 L	1	1
FMS301047	Flexsafe® for LevMixer® 50 L	1	1
FMS301048	Flexsafe® for LevMixer® 100 L	1	1
FMS301049	Flexsafe® for LevMixer® 200 L	2	2
FMS301050	Flexsafe® for LevMixer® 400 L	2	2
FMS301051	Flexsafe® for LevMixer® 650 L	1	1
FMS301052	Flexsafe® for LevMixer® 1,000 L	1	1

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178



For more information, visit
[sartorius.com](https://www.sartorius.com)

©2026 Sartorius. All rights reserved. Clipster®, Flexel®, Flexsafe®, Palletank®, and Tuflux® are registered trademarks of Sartorius or its subsidiaries.

AseptiQuik is a registered or unregistered trademark of Colder Products Company. Clave is a registered or unregistered trademark of ICU Medical, Inc. LevMixer is a registered or unregistered trademark of Pall Corporation. All other third-party trademarks are the property of their respective owners.

For details on the registrations please refer to <https://www.sartorius.com/en/patents-and-trademarks>

Last modified: 06 | 2026