

TakeOne®

Simplified Sampling Solutions for Stainless Steel Equipment



Product Information

TakeOne® aseptic sampling solutions have been designed to integrate seamlessly with stainless steel processes and collect perfectly representative samples with no risk of false-positives for each process step and scale. TakeOne® products are delivered fully assembled, in individual packs, gamma sterilized, integrity tested and ready for immediate use allowing you to bypass parts washing, equipment prep and assembly with confidence.

Features and Benefits

- Ensures representative sampling, avoiding dead-leg zones
- Cost-effective with a ready-to-use, single-use design
- Suitable for all stages, from R&D to GMP manufacturing
- Offers a diverse range of collection containers for different assays, including bags, bottles, tubes, and valves for syringe access
- Available in a variety of fluid contact materials and container sizes
- Designed for ease of use and ergonomic handling

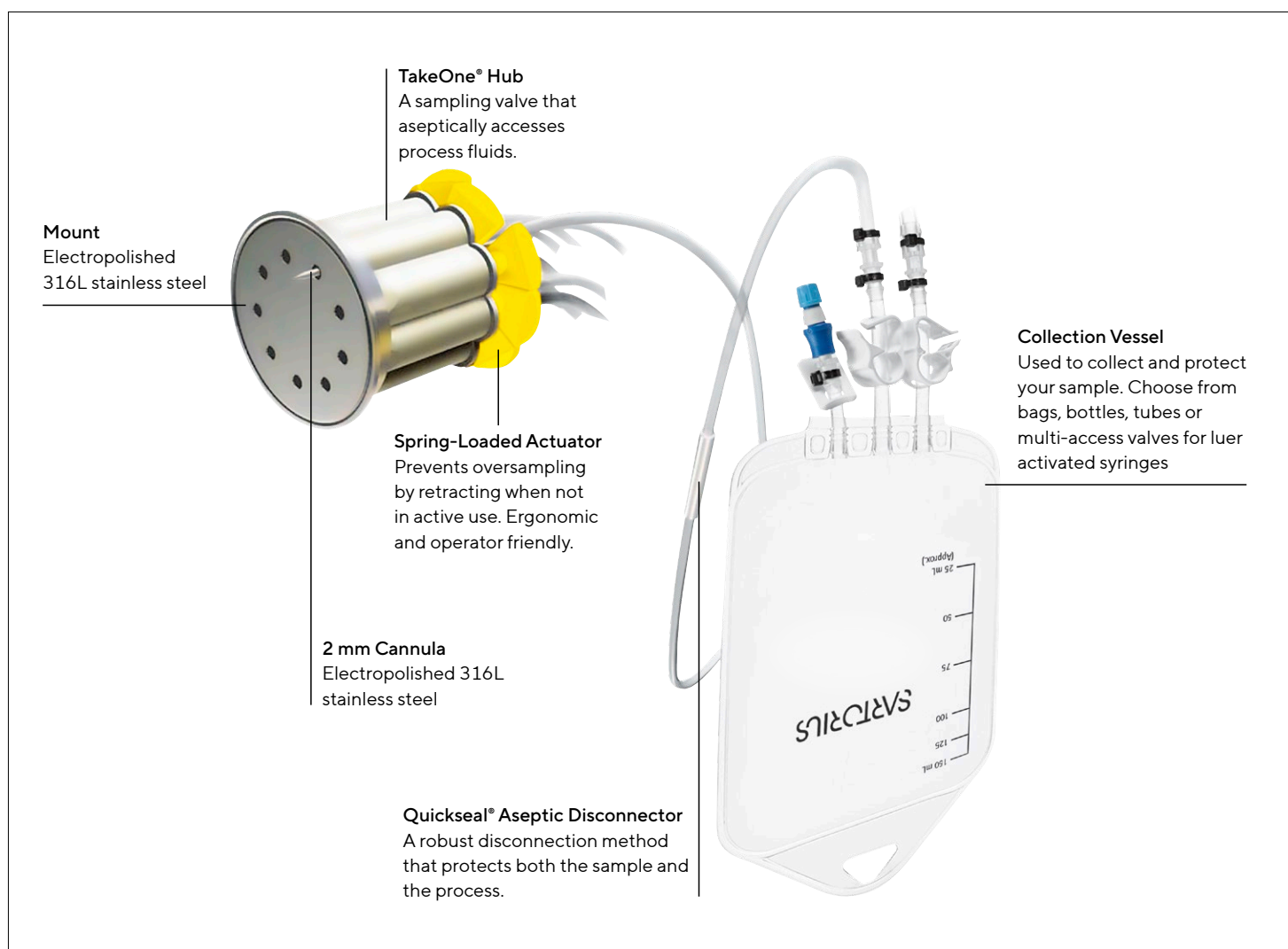
Introduction

TakeOne® Aseptic Sampling Solutions for Stainless Steel Equipment

Aseptic sampling is a principal component of an effective microbial control program. Not all samples are equal and we don't treat them that way. Whether it is sterility, bioburden, endotoxin testing or routine process monitoring, TakeOne® has been designed to meet your processes demands.

This aseptic sampling solution has been engineered to take representative samples, giving exact results with no risk of false-positives thanks to its pre-assembled, sterile and closed design. It collects samples for analysis while protecting the process fluid from adventitious agents.

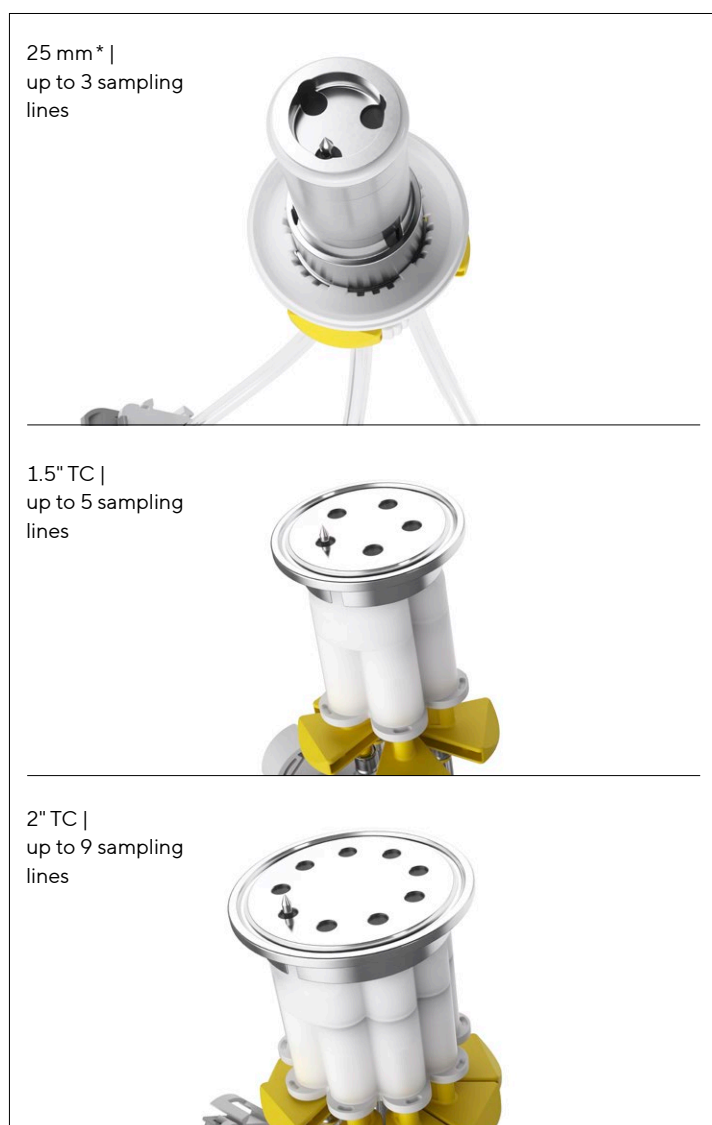
Figure 1: *Key Components*



Robust, Safe for Operators And Protective of Process Integrity

Careful selection of the materials of construction means TakeOne® sampling solutions are engineered to perform. TakeOne® connects to the sampling point with a 316L stainless steel tank mount that was selected to optimize safety and robustness. Steel-to-steel fittings are secure and trusted during thermal cycling and high pressure, especially steam-in-place. Any other material would be a compromise on safety, susceptible to leaking and be difficult to validate.

Figure 2: *Three Fitting Formats*

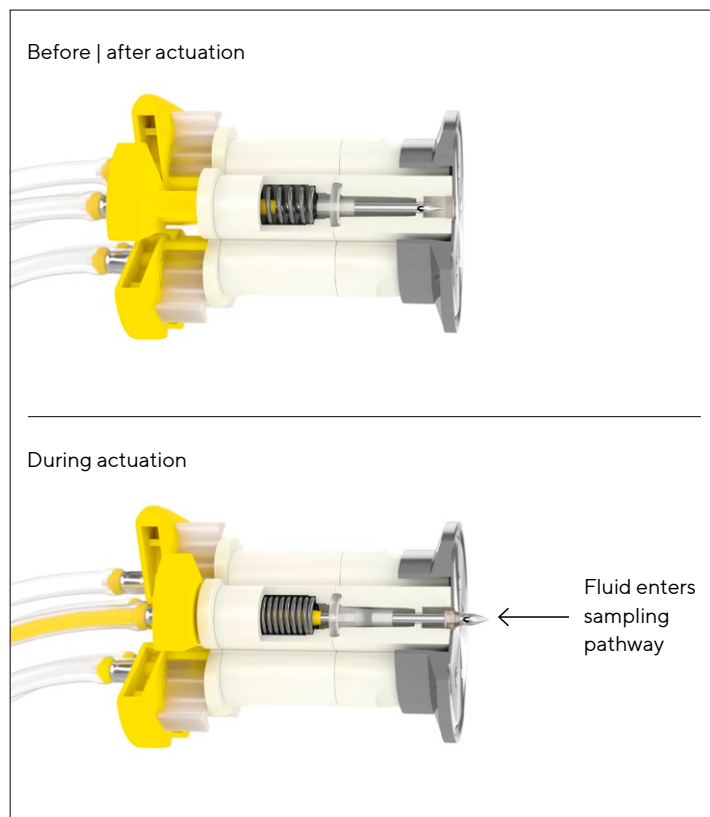


*25 mm TakeOne® fitting requires an accessory to adapt the vessel's native, threaded port to make it tri clamp compatible. Our adaptors are 316L stainless steel. Once installed, they can remain in place through use and cleaning | sanitization procedures.

Completely single-use which eliminate the risk of contamination from improper or ineffective cleaning and incorrect assembly of reusable parts.

- Ready-to-use and has no spare, reused parts and is easy to install, manage and use. Eliminate the risk of production delays or batch contaminations from lost or damaged parts.
- TakeOne® sampling solutions are assembled in ISO 7 certified cleanrooms. Each unit undergoes a variety of in-process quality inspections including 100% visual inspection and 100% leak testing (pressure decay) of each sampling line ensures proper assembly and integrity of the internal aseptic chamber.
- The cleverly designed, internal aseptic chamber houses the sampling needle and remains intact, preserving the aseptic sampling pathway, after dozens of actuations, protecting the production process from contamination.

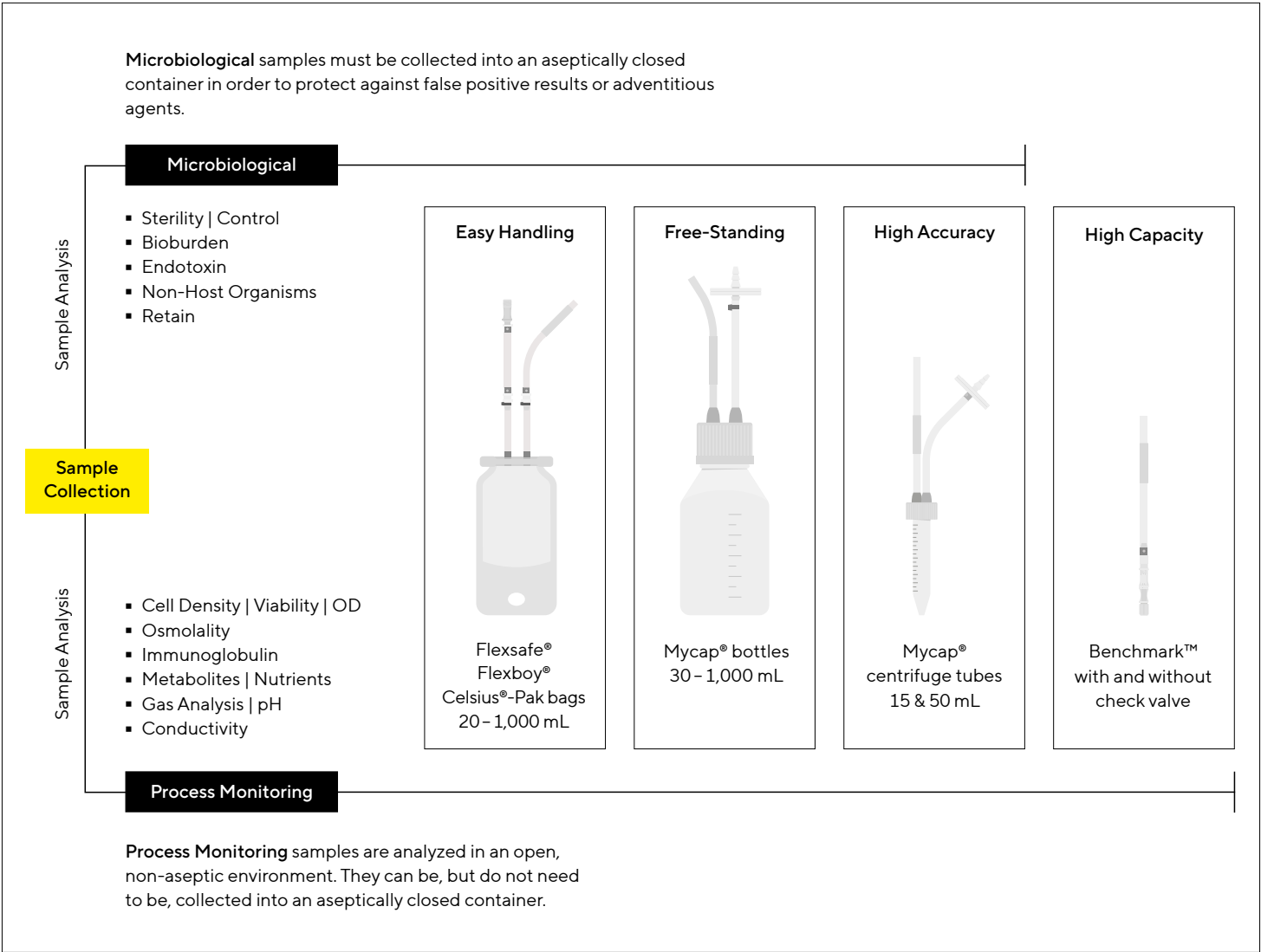
Figure 3: *Aseptic Chamber*



Simply Smarter Sampling

Categorization by sample type reveals the ideal balance between cost per sample and sampling capacity.

Figure 4: Sampling Container Selection Based on Assay Type

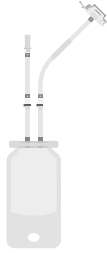



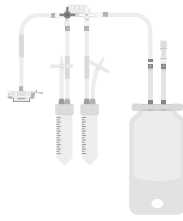
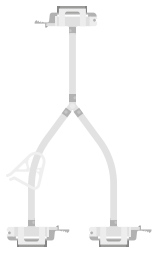


Collection Container Types

TakeOne® products are available with a diverse variety of container types in different fluid contact materials, volumes and formats to accommodate various process fluids, assay types and handling preferences.

Sartorius’ wide range includes bags in 3 different films, multi-access sampling valves, as well as, bottles and tubes that feature Mycap®, a robust aseptic closure suitable for all biopharmaceutical applications. Designs are available in both TPE and silicone tubing variations with aseptic connectors or sealed weldable tubing.

Table 1: TakeOne® Flex Is Available in a Wide Range of Types, Materials, and Volumes

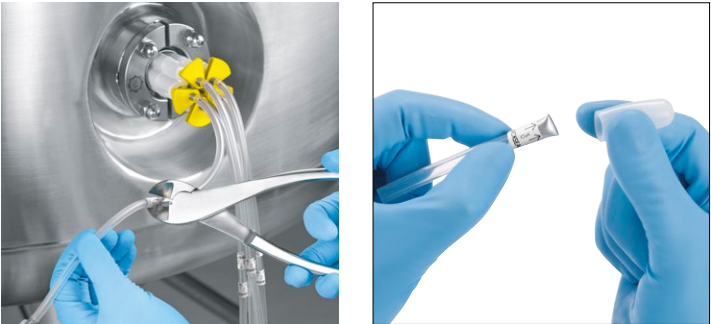
					
Flexsafe®, Flexboy®, Celsius® Bags*	Mycap® Centrifuge Tubes	Mycap® Bottles	Benchmark™ Needleless Access Site	Stopcock Manifolds, incl. Purge Bag	2-Way Splitter
20 30 50 100 150 250 500 1,000 mL	15 50 mL	30 60 125 250 500 1,000 mL	With without check valve	2 5 collection vessels	Aseptic connections
PE EVA	PS PP	PETG (other materials available upon request)	SIL PP SAN	Per your selection	TPE SIL

*Per your bag brand selection

Quick and Easy Disconnect

Each TakeOne® sampling line includes a Quickseal® aseptic disconnect. Once the sample has been collected, operators easily cut the aluminum collar with a light and portable hand-held cutting tool. Quickseal® patented technology protects the sample and process vessel from contamination. A Quickseal® silicone protective cap then shields the cut collar.

Figure 5: How to Use Quickseal®



Integrating TakeOne® Into Your Processes

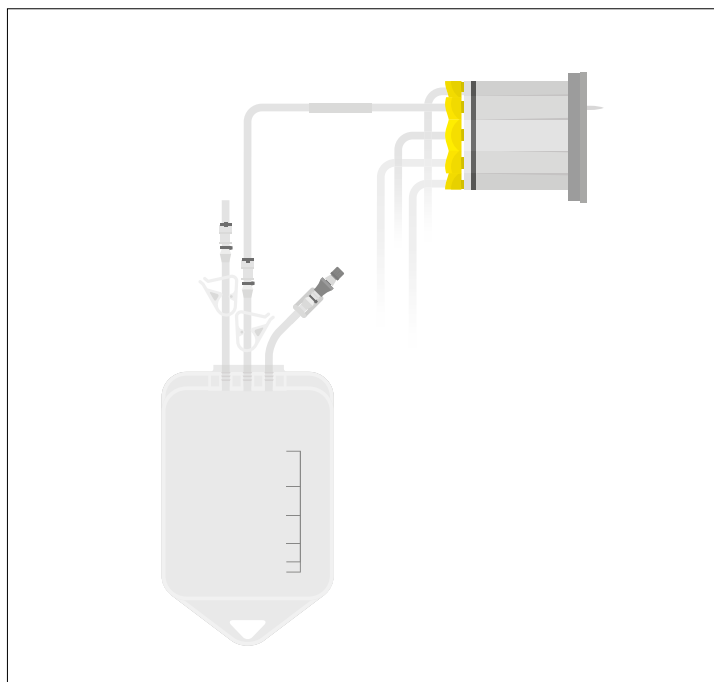
TakeOne® sampling solutions for stainless steel equipment are available in two flexible families to ensure all of your varied process needs are met from R&D to Process Development and full scale GMP production.

TakeOne® Standard to Order—Pre-Designed Configurations—A Complete Sampling Solution

TakeOne® Pre-Designed Configurations. Select and go, with aseptic sampling solutions that have sampling lines and collection containers preinstalled to the hub and ready for immediate installation on your vessel. You can save time, avoid handling needles and benefit from factory installed and pressure tested assemblies.

Our broad portfolio of individual and manifolded collection containers is available across a wide variety of standard Sartorius designs you can select from. We also offer custom configurations tailored to meet the exact needs of your sampling program. Please contact your local Sartorius representative for a custom configuration from our design services offerings.

Figure 6: *TakeOne® Standard to Order*



TakeOne® Flex—A modular Approach to Sampling

Sartorius recognizes that sampling can be complex and process-specific. TakeOne® Flex was created to meet these widely varied process demands and take the complexity out of sample planning by offering a modular range of container formats, volumes, and fluid contact materials and sampling hubs. Plan and adapt to any process situation by maintaining just a few part numbers (PN), and create tens or thousands of different sampling solutions while still optimizing for the most accurate and repeatable results using the precise volume the sample requires.

This plug and play approach allows you to quickly react on demand with point-of-use configuration that provide an ideal solution for unplanned or variable sampling needs. React with TakeOne® Flex by simply selecting and connecting the appropriate sample collection container at the moment the sample is required.

Sampling Hubs

Sartorius offers the TakeOne® aseptic sampling hubs with 2 types of aseptic connections options to provide configuration flexibility at your point-of-use, even mid-process.

Save time and avoid handling needles — sampling lines have already been pre-assembled on to the hub which is ready for immediate installation on your vessel. Weldable and plugged tubing or genderless AseptiQuik® connectors are used for fast and easy assembly of the sampling containers to the hubs

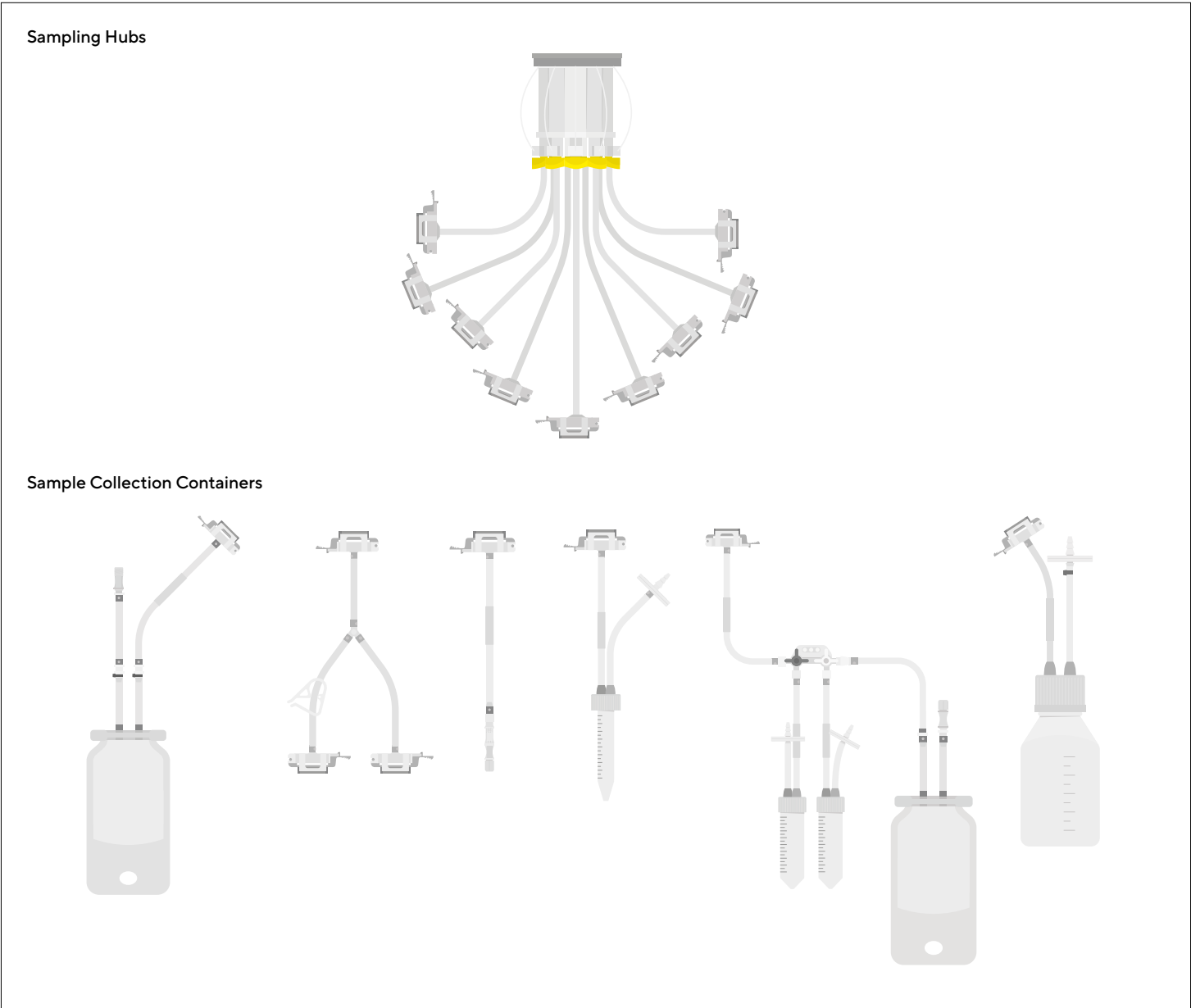
Sample Collection Containers

Our broad portfolio of individual or manifolded collection containers are available decoupled, so they can be aseptically connected to sampling hubs when and as needed, providing near endless flexibility to your sampling program. They can also be used as part of a comprehensive sampling platform, as they work equally well for sampling from single-use equipment and assemblies.

Easily Reduce and Expand

Avoid overwhelming your inventory management program. Maintain just a PN's and configure tens or thousands of different sampling solutions right at the point-of-use.

Figure 7: *Example of TakeOne® Flex Products*



TakeOne® Flex Unique Configuration Possibilities

With just a few sample container part numbers, you can create a multitude of different sampling configurations for each sampling hub used:

The number of different sampling container arrangements that can be configured with a sampling hub based on the number of sampling lines available, per fitting size, and the number of different sampling containers ordered.

Example: 2 different sampling container PN's and 1 × 3-line 25 mm hub allows you to configure up to 6 different sampling solutions.

10 different sampling containers PN's and 1 × 9-line 2" hub PN – allows you to configure up to 92,377 different sampling solutions.

Table 2: Possible Configuration Variations




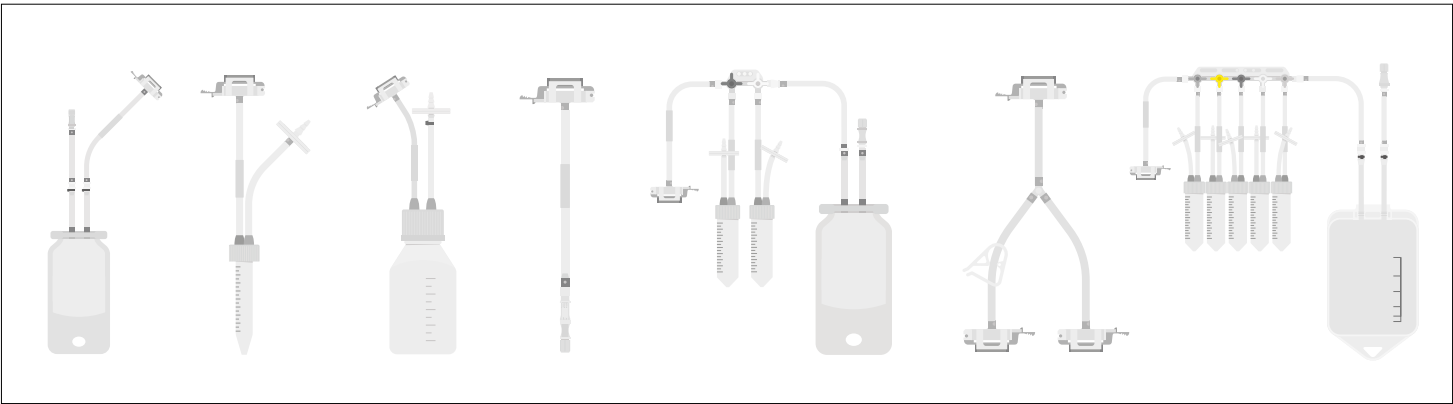
Sampling Hub Type	25 mm 3 Lines	1.5" TC 5 Lines	2" TC 9 Lines
Quantity of Unique Collection Container PN's On-Hand			
	Number of Possible Configuration Variations		
	2	20	54
	4	125	714
	6	461	5,004
	8	1,286	24,309
	10	3,002	92,377

Figure 10: Examples of Collection Container PN's



Design Services

When you need something fine-tuned to your exact needs, we also offer custom configurations tailored to meet the exact needs of your sampling program.

Configured to Order

Easily configure a sampling device with your Sartorius representative.

- Qualified design space
- Select from well characterized components
- Tubing materials
- Fitting types
- Collection containers

Engineered to Order

Fully engineered solutions tailored to your needs.

- Application-based expertise to help optimize your process
- Modify standard features or build completely customized designs
- A wider variety of components offered:
- Alternative tube materials
- Collection vessels
- Fittings
- Etc.

Technical Specifications

Equipment Interface | Mounting Mechanism

25 mm port Ingold® port	Up to 3 sampling lines
1.5" Sanitary TC close coupled	Up to 5 sampling lines
2" Sanitary TC close coupled	Up to 9 sampling lines

Operating Parameters for Sampling Hubs

Max SIP	1 cycle @60 mins at 43 psi 3 bar 143.7 °C
SIP cycles after actuation	10 cycles @60 mins at 29 psi 2 bar 134 °C
Actuations per line	40
Temperature range	-55 °C to 300 °C
Operating temperature range	-20 °C to 50 °C (configuration dependent)
Device burst pressure	129 psi 8.9 bar before actuation 105 psi 7.3 bar after actuation
Recommended max operating pressure	45 psi 3.1 bar at 25 °C (do not overfill or pressurize collection vessels)

Manufacturing

Manufacturing environment	ISO 7
Sterilization method	Gamma irradiation (25 kGy - 45 kGy)
Device release criteria	100% leak-tested (pressure decay)

Materials of Contact (Device Only)

See Collection Container Ordering Information for more details

Septum	Platinum-cured silicone
Equipment interface	316L stainless steel, electropolished (< 15 µin Ra)
2 mm cannula	316L stainless steel, electropolished (< 15 µin Ra)
Tubing material	Silicone (Pt), C-Flex® (374) or Tuflux® TPE

Ordering Information

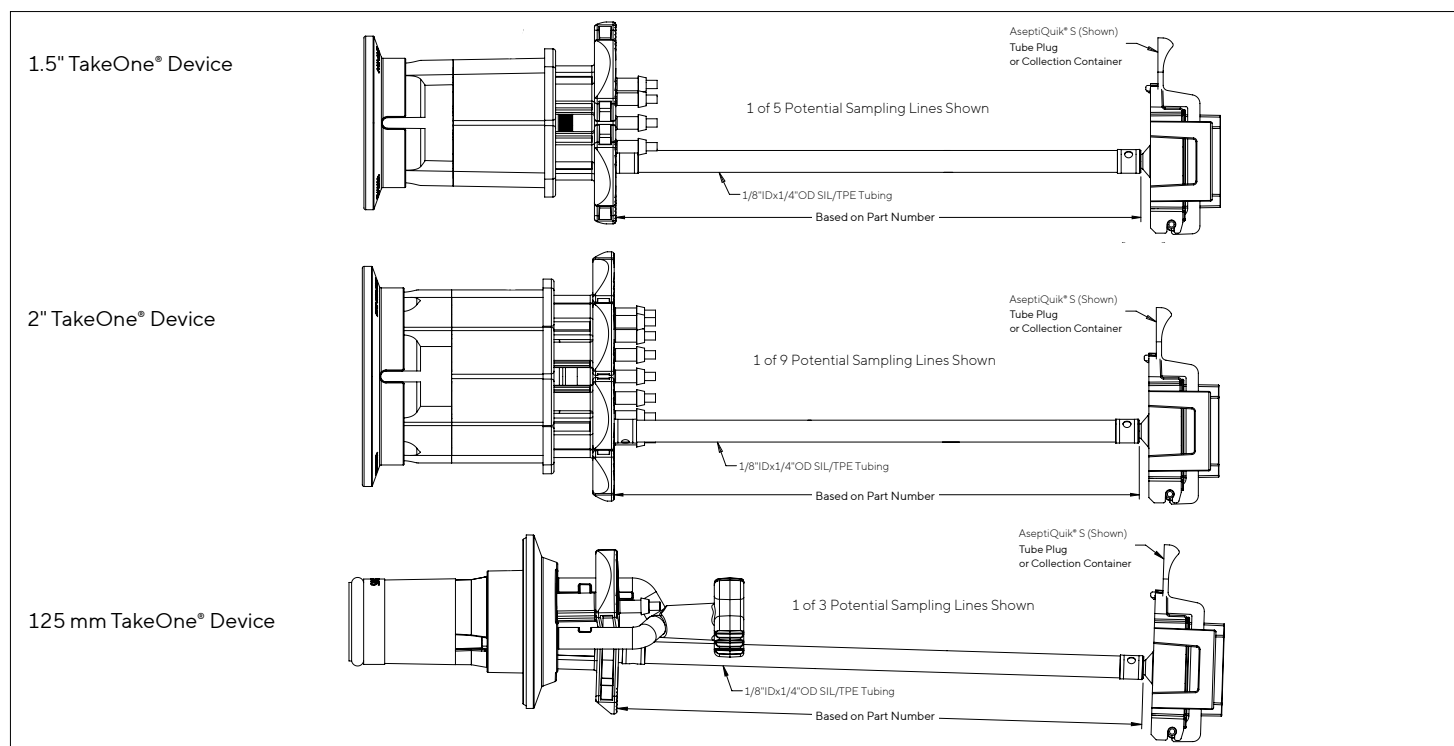
TakeOne® STO – Pre-Designed Configurations*

For nominal assembly dimensions please refer to the illustrations for TakeOne® Flex sampling hubs (omit tubing) and to the illustrations for the appropriate TakeOne® Flex Sample Collection Containers (omit aseptic connector). In combination you can assess the preconfigured assembly's approximate overall footprint.

Hub Size and Fitting	Sampling Lines	Catalog Number	Description	Fluid Contact Materials
25 mm port (Ingold®)	3	INGP03STOT0001	TakeOne® STO, 25 mm, 3 × 60 mL PETG bottle, Tuflux® TPE	316 L stainless steel, silicone (Pt), high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), Tuflux® TPE
		INGP03STOT0002	TakeOne® STO, 25 mm, 3 × 15 mL centrifuge tube, Tuflux® TPE	316 L stainless steel, silicone (Pt), polystyrene, high density polyethylene (HDPE), Tuflux® TPE
		INGP03STOT0003	TakeOne® STO, 25 mm, 3 × 250 mL Flexsafe® bag, Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, polyethylene (PE S80), Tuflux® TPE
		INGP03STOT0004	TakeOne® STO, 25 mm, 3 × (2 × 50 mL Flexsafe® bag), Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, polyethylene (PE S80), high density polyethylene (HDPE), Tuflux® TPE
		INGP03STOT0005	TakeOne® STO, 25 mm, 3 × (2 × 60 mL PETG bottle), Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), Tuflux® TPE
1.5" Sanitary TC	5	24TC05STOT0001	TakeOne® STO, 1.5", 5 × 50 mL Flexsafe® bag, Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, polyethylene (PE S80), Tuflux® TPE
		24TC05STOT0002	TakeOne® STO, 1.5", 5 × 250 mL Flexsafe® bag, Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, polyethylene (PE S80), Tuflux® TPE
		24TC05STOT0003	TakeOne® STO, 1.5", 5 × 60 mL PETG bottle, Tuflux® TPE	316 L stainless steel, silicone (Pt), high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), Tuflux® TPE
		24TC05STOT0004	TakeOne® STO, 1.5", 5 × 15 mL centrifuge tube, Tuflux® TPE	316 L stainless steel, silicone (Pt), polystyrene, high density polyethylene (HDPE), Tuflux® TPE
		24TC05STOT0005	TakeOne® STO, 1.5", 5 × (2 × 50 mL Flexsafe® bag), Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, polyethylene (PE S80), high density polyethylene (HDPE), Tuflux® TPE
		24TC05STOT0006	TakeOne® STO, 1.5", 5 × (2 × 60 mL PETG bottle), Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), Tuflux® TPE
2" Sanitary TC	9	32TC09STOT0001	TakeOne® STO, 2", 9 × 50 mL Flexsafe® bag, Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, polyethylene (PE S80), Tuflux® TPE
		32TC09STOT0002	TakeOne® STO, 2", 9 × 250 mL Flexsafe® bag, Tuflux® TPE	316 L Stainless steel, Silicone (Pt), polycarbonate, polyethylene (PE S80), Tuflux® TPE
		32TC09STOT0003	TakeOne® STO, 2", 9 × 15 mL centrifuge tube, Tuflux® TPE	316 L stainless steel, silicone (Pt), polystyrene, high density polyethylene (HDPE), Tuflux® TPE
		32TC09STOT0004	TakeOne® STO, 2", 9 × 60 mL PETG bottle, Tuflux® TPE	316 L stainless steel, silicone (Pt), high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), Tuflux® TPE
		32TC09STOT0005	TakeOne® STO, 2", 9 × (2 × 50 mL Flexsafe® bag), Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, high density polyethylene (HDPE), polyethylene (PE S80), Tuflux® TPE
		32TC09STOT0006	TakeOne® STO, 2", 9 × (2 × 60 mL PETG bottle), Tuflux® TPE	316 L stainless steel, silicone (Pt), polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), Tuflux® TPE

*Products come individually double bagged
Ingold® is a registered trademark of Mettler-Toledo

TakeOne® Flex – Point-of-Use Configuration



Sampling Hub Bodies | Silicone (Pt)* With Aseptic Connector

Hub Size and Fitting	Sampling Lines	Catalog Number	Tube Material	Sampling Line Connection Type	Fluid Contact Materials
1.5" Sanitary TC	3	24TC03PASQLT01	Silicone (Pt) 1/8" x 1/4" Length 6"	AseptiQuik® S	316L stainless steel, silicone (Pt), polycarbonate
1.5" Sanitary TC	5	24TC05PASQLT01			
2" Sanitary TC	5	32TC05PASQLT01			
2" Sanitary TC	9	32TC09PASQLT01			
25 mm port (Ingold®)	3	INGP03PASQLT01			

Sampling Hub Bodies | TPE With Aseptic Connector*

Hub Size and Fitting	Sampling Lines	Catalog Number	Tube Material	Sampling Line Connection Type	Fluid Contact Materials
1.5" Sanitary TC	3	24TC03CASQLT01	C-Flex® 1/8" x 1/4" Length 6"	AseptiQuik® S	316L stainless steel, silicone (Pt), polycarbonate, C-Flex® (374)
	5	24TC05CASQLT01			
2" Sanitary TC	5	32TC05CASQLT01			
	9	32TC09CASQLT01			
25 mm port (Ingold®)	3	INGP03CASQLT01			
1.5" Sanitary TC	5	24TC05TPLG01	Tuflux® TPE 1/8" x 1/4" Length 12"	Plug for welding	316L stainless steel, silicone (Pt), polypropylene
2" Sanitary TC	9	32TC09TPLG01			
25 mm port (Ingold®)	3	INGP03TPLG01			

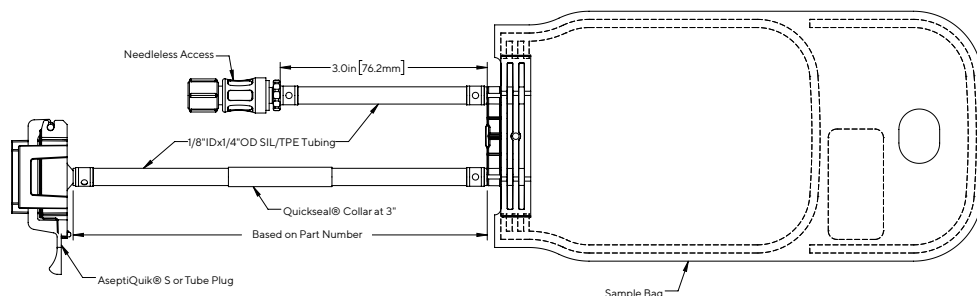
Sampling Hub Bodies | TPE for Sterile Welding*

Hub Size and Fitting	Sampling Lines	Catalog Number	Tube Material	Sampling Line Connection Type	Fluid Contact Materials
1.5" Sanitary TC	5	24TC05TPLG01	Tuflux® TPE 1/8" x 1/4" Length 12"	Plug for welding	316L stainless steel, silicone (Pt), polypropylene
2" Sanitary TC	9	32TC09TPLG01			
25 mm port (Ingold®)	3	INGP03TPLG01			

* Products come individually double bagged.
Ingold® is a registered trademark of Mettler-Toledo

TakeOne® Flex | Sample Containers

Single Collection Container



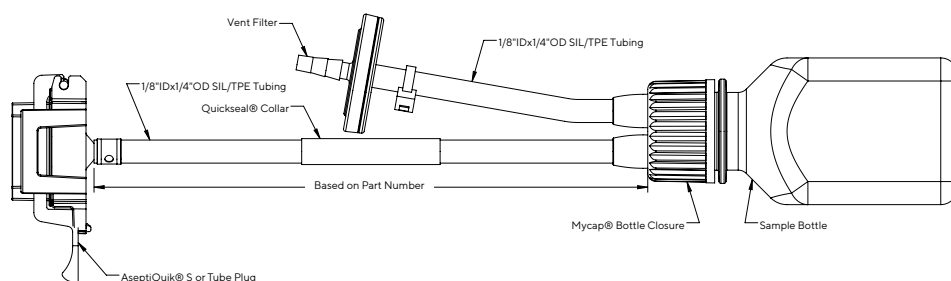
Bags With Aseptic Connector

Container Type	Volume	Catalog Number TPE Tube	Catalog Number Silicone (Pt) Tube	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Flexsafe® bag	20 mL	FLS0021CASQLT	FLS0021PASQLT	AseptiQuik® S ⅝" x ¼" Length 6" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Polyethylene (S80)	Polycarbonate, polypropylene (PP), silicone (Pt)
	50 mL	FLS0051CASQLT	FLS0051PASQLT				
	150 mL	FLS0151CASQLT	FLS0151PASQLT				
	250 mL	FLS0251CASQLT	FLS0251PASQLT				
	500 mL	FLS0501CASQLT	FLS0501PASQLT				
	1,000 mL	FLS1001CASQLT	FLS1001PASQLT				
Flexboy® bag	50 mL	FBY0051CASQLT	FBY0051PASQLT	AseptiQuik® S ⅝" x ¼" Length 6" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Ethylene-vinyl acetate (EVA S71)	Polycarbonate, polypropylene (PP), silicone (Pt)
	150 mL	FBY0151CASQLT	FBY0151PASQLT				
	250 mL	FBY0251CASQLT	FBY0251PASQLT				
	500 mL	FBY0501CASQLT	FBY0501PASQLT				
	1,000 mL	FBY1001CASQLT	FBY1001PASQLT				
Celsius® Pak bag	30 mL	CSP0031CASQLT	—	AseptiQuik® S ⅝" x ¼" Length 6" Quickseal® 3"	Luer fitting	Ethylene-vinyl acetate (EVA S71-360 µm)	Polycarbonate, polypropylene (PP), silicone (Pt)
	100 mL	CSP0101CASQLT	—				

Bags for Sterile Welding

Container Type	Volume	Catalog Number	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Flexsafe® bag	20 mL	FLS0021TPLG	Tuflux® TPE with Plug ⅝" x ¼" Length 12" Quickseal® 3"	Benchmark™ for needle-free syringe	Polyethylene (S80)	Polycarbonate, polypropylene, silicone (Pt)
	50 mL	FLS0051TPLG				
	150 mL	FLS0151TPLG				
	250 mL	FLS0251TPLG				
	500 mL	FLS0501TPLG				
	1,000 mL	FLS1001TPLG				
Celsius® Pak bag	30 mL	CSP0031TPLG	Tuflux® TPE with Plug ⅝" x ¼" Length 12" Quickseal® 3"	Luer fitting	Ethylene-vinyl acetate (EVA S71-360 µm)	Polypropylene
	100 mL	CSP0101TPLG				

*List does not include tubing type material



Bottles With Aseptic Connector

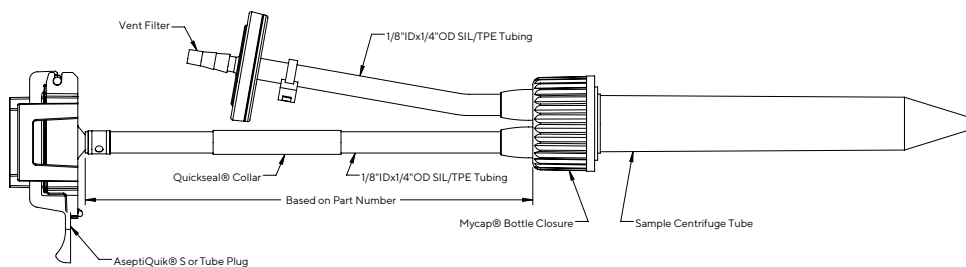
Container Type	Volume	Catalog Number TPE Tube	Catalog Number Silicone (Pt) Tube	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Square bottle	30 mL	PTG0031CASQLT	PTG0031PASQLT	AseptiQuik® S 1/8" x 1/4" Length 6" Quickseal® 3"	Minisart® vent filter	Polyethylene terephthalate glycol-modified (PETG)**	Polycarbonate, silicone (Pt)
	60 mL	PTG0061CASQLT	PTG0061PASQLT				
	125 mL	PTG0121CASQLT	PTG0121PASQLT				
	250 mL	PTG0251CASQLT	PTG0251PASQLT				
	500 mL	PTG0501CASQLT	PTG0501PASQLT				
	1,000 mL	PTG1001CASQLT	PTG1001PASQLT				

Bottles for Sterile Welding

Container Type	Volume	Catalog Number	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Square bottles	30 mL	MYCAPNGPG0030F	Tuflux® TPE with Plug 1/8" x 1/4" Length 15.5" Quickseal® 3" Diptube	25 mm Minisart® vent filter	Polyethylene terephthalate glycol-modified (PETG)**	HDPE, polycarbonate, polypropylene, silicone (Pt)
	60 mL	MYCAPNGPG0060F	Tuflux® TPE with Plug 1/8" x 1/4" Length 15.5" Quickseal® 3" Diptube	25 mm Minisart® vent filter		
	125 mL	MYCAPNGPG0125F	Tuflux® TPE with Plug 1/8" x 1/4" Length 14" Quickseal® 2" Diptube			
	250 mL	MYCAPNGPG0250F	Tuflux® TPE with Plug 1/8" x 1/4" Length 18" Quickseal® 2" Diptube		Polyethylene terephthalate glycol-modified (PETG)**	Polycarbonate, polypropylene, silicone (Pt)
	500 mL	MYCAPNGPG0500F	Tuflux® TPE with Plug 1/8" x 1/4" Length 17" Quickseal® 3" Diptube	50 mm Midisart® vent filter		
	1,000 mL	MYCAPNGPG1,000F	Tuflux® TPE with Plug 1/8" x 1/4" Length 21.75" Quickseal® 4.5" Diptube			

* List does not include tubing type material

** Other materials available upon request



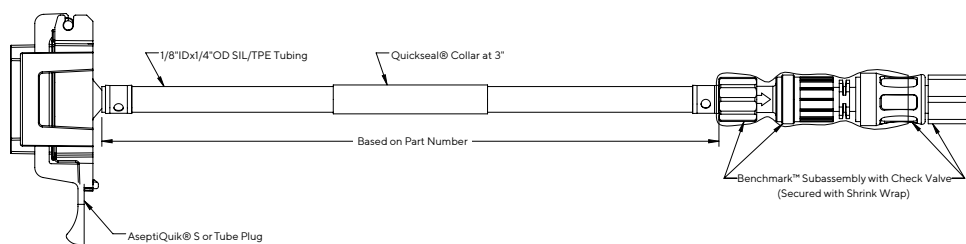
Centrifuge Tubes With Aseptic Connector

Container Type	Volume	Catalog Number TPE Tube	Catalog Number Silicone (Pt) Tube	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Centrifuge tube	15 mL	PSCT0151CASQLT	PSCT0151PASQLT	AseptiQuik® S ⅝" x ¼" Length 6" Quickseal® 3"	Minisart® vent filter	Polystyrene	Polycarbonate, high-density polyethylene (HDPE), silicone (Pt)
	50 mL	PPCT0051CASQLT	PPCT0051PASQLT	AseptiQuik® S ⅝" x ¼" Length 6" Quickseal® 3"	Minisart® vent filter	Polypropylene (PP)	Polycarbonate, silicone (Pt)

Centrifuge Tubes for Sterile Welding

Container Type	Volume	Catalog Number	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Centrifuge tube	15 mL	MYCAPCPSG0015F	Tuflux® TPE with Plug ⅝" x ¼" Length 13.5" Quickseal® 2.5" Diptube	25 mm Minisart® vent filter	Polystyrene	High density polyethylene (HDPE), silicone (Pt), polypropylene
	50 mL	MYCAPCPPG0050F	Tuflux® TPE with Plug ⅝" x ¼" Length 13.5" Quickseal® 2" Diptube	25 mm Minisart® vent filter	Polypropylene (PP)	Silicone (Pt), polypropylene

*List does not include tubing type material



Benchmark™ Needleless Access Site With Aseptic Connector

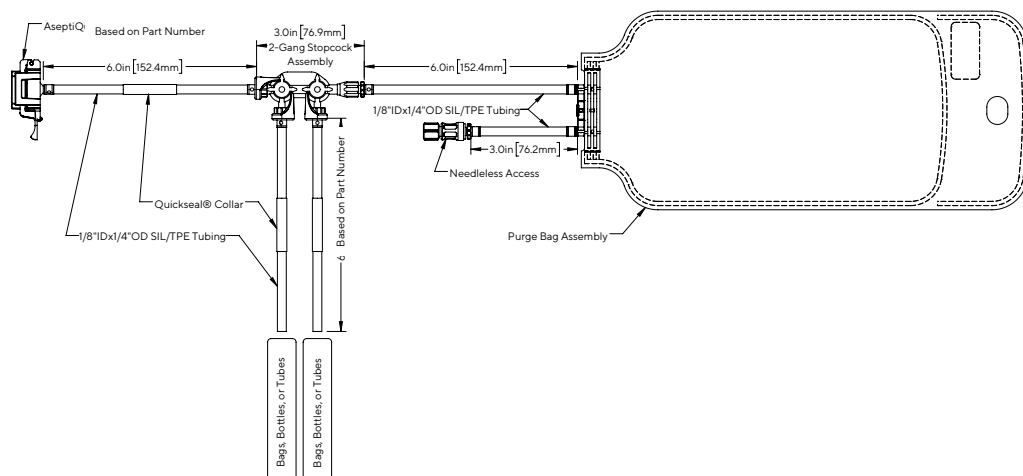
	Catalog Number TPE Tube	Catalog Number Silicone (Pt) Tube	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Benchmark™ with check valve	BNCHCV1CASQLT	BNCHCV1PASQLT	AseptiQuik® S ⅜" × ¼" Length 6" Quickseal® 3"	—	As selected	Polycarbonate, polypropylene (PP), styrene acrylonitrile (SAN), silicone (Pt)
w/out check valve	BNCH1CASQLT	BNCH1PASQLT	AseptiQuik® S ⅜" × ¼" Length 6" Quickseal® 3"	—	As selected	Polycarbonate, silicone (Pt)

Benchmark™ Needleless Access Site for Sterile Welding

Catalog Number	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Benchmark™ BNCHCV1TPLG	Tuflux® TPE with Plug ⅜" × ¼" Length 12" Quickseal® 3"	2 × Tuflux® TPE with Plug ⅜" × ¼" Length 12" Quickseal® 3", Benchmark™ with check valve	—	Polypropylene, silicon (Pt), polycarbonate, SAN
BNCH1TPLG	Tuflux® TPE with Plug ⅜" × ¼" Length 12" Quickseal® 3"	3 × Tuflux® TPE with Plug ⅜" × ¼" Length 12" Quickseal® 3", Benchmark™	—	Polypropylene, silicon (Pt), polycarbonate

*List does not include tubing type material

Manifolded Collection Containers



2-Gang Manifold With Aseptic Connector

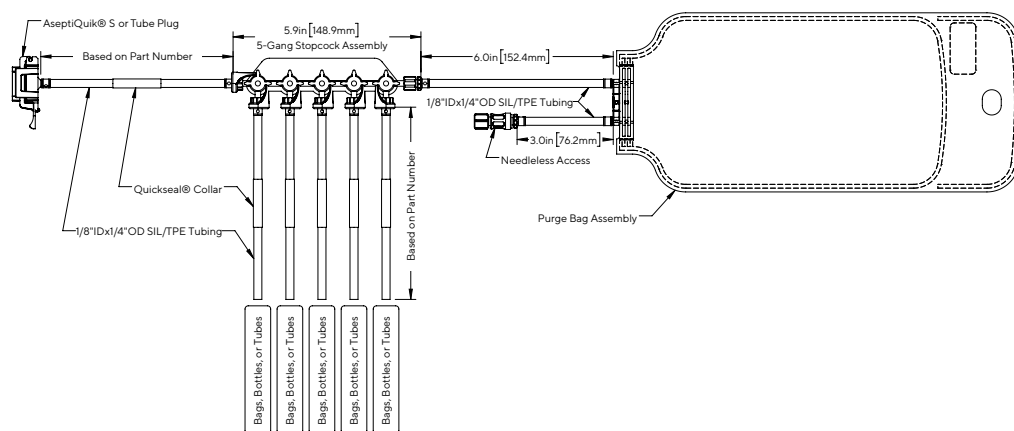
Container Type	Volume	Catalog Number TPE Tube	Catalog Number Silicone (Pt) Tube	Container Inlet	Container Outlet	Fluid Contact Materials*
Flexsafe® bags on stopcock manifold	2 × 20 mL	FLS0022CASQLT	FLS0022PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Polycarbonate, high-density polyethylene (HDPE), polyethylene (PE S80), polypropylene (PP), silicone (Pt)
	2 × 50 mL	FLS0052CASQLT	FLS0052PASQLT			
	2 × 150 mL	FLS0152CASQLT	FLS0152PASQLT			
	2 × 250 mL	FLS0252CASQLT	FLS0252PASQLT			
	2 × 500 mL	FLS0502CASQLT	FLS0502PASQLT			
	2 × 1,000 mL	FLS1002CASQLT	FLS1002PASQLT			
Flexboy® bags on stopcock manifold	2 × 50 mL	FBY0052CASQLT	FBY0052PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Polycarbonate, high-density polyethylene (HDPE), ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone (Pt)
	2 × 250 mL	FBY0252CASQLT	FBY0252PASQLT			
	2 × 1,000 mL	FBY1002CASQLT	FBY1002PASQLT			
Centrifuge tube on stopcock manifold	2 × 15 mL	PSCT0152CASQLT	PSCT0152PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Minisart® vent filter	Polycarbonate, polystyrene, high density polyethylene (HDPE), silicone (Pt)
	2 × 50 mL	PPCT0052CASQLT	PPCT0052PASQLT			
Bottle on stopcock manifold	2 × 60 mL	PTG0062CASQLT	PTG0062PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Minisart® vent filter	Polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), silicone (Pt)
2-way splitter		2WAY001CASQLT	2WAY001PASQLT	AseptiQuik® S ⅝" × ¼" Length 3"	AseptiQuik® S (× 2) ⅝" × ¼" Length 3"	Polycarbonate, high density polyethylene (HDPE), silicone (Pt)

*List does not include tubing type material

2-Gang Manifold for Sterile Welding

Container Type	Volume	Catalog Number	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Flexsafe® bags on stopcock manifold	2 × 20 mL	FLS0022TPLG	Tuflux® with Plug ⅝" × ¼" Length 12" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Polyethylene (S80)	Polycarbonate, high-density polyethylene (HDPE), polyethylene, polypropylene (PP), silicone (Pt)
	2 × 50 mL	FLS0052TPLG				
	2 × 150 mL	FLS0152TPLG				
	2 × 250 mL	FLS0252TPLG				
	2 × 500 mL	FLS0502TPLG				
	2 × 1,000 mL	FLS1002TPLG				
Centrifuge tube on stopcock manifold	2 × 15 mL	PSCT0152TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	25 mm Minisart® vent filter	Polystyrene	Polycarbonate, high-density polyethylene (HDPE), polyethylene (S80), polypropylene (PP), silicone (Pt)
	2 × 50 mL	PPCT0052TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	25 mm Minisart® vent filter	Polypropylene	Polycarbonate, high-density polyethylene (HDPE), polyethylene (S80), polyethylene, polypropylene (PP), silicone (Pt)
Bottle on stopcock manifold	2 × 60 mL	PTG0062TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	25 mm Minisart® vent filter	Polyethylene terephthalate glycol-modified (PETG)	Polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), silicone (Pt), polypropylene, polyethylene (S80), polyethylene
2-way splitter	—	2WAY001TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	Tuflux® TPE with Plug (× 2) ⅝" × ¼" Length 12" Quickseal® 3"	—	Polypropylene

*List does not include tubing type material



5-Gang Manifold With Purge Vessel With Aseptic Connector

Container Type	Volume	Catalog Number TPE Tube	Catalog Number Silicone (Pt) Tube	Container Inlet	Container Outlet	Fluid Contact Materials*
Flexsafe® bags on stopcock manifold	5 × 20 mL	FLS0025CASQLT	FLS0025PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), polyethylene (PE S80), polypropylene (PP), silicone (Pt)
	5 × 50 mL	FLS0055CASQLT	FLS0055PASQLT			
	5 × 150 mL	FLS0155CASQLT	FLS0155PASQLT			
	5 × 250 mL	FLS0255CASQLT	FLS0255PASQLT			
	5 × 500 mL	FLS0505CASQLT	FLS0505PASQLT			
	5 × 1,000 mL	FLS1005CASQLT	FLS1005PASQLT			
Flexboy® bags on stopcock manifold	5 × 50 mL	FBY0055CASQLT	FBY0055PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Benchmark™ (for needle-free syringes)	Polycarbonate, high-density polyethylene (HDPE), ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone (Pt)
	5 × 250 mL	FBY0255CASQLT	FBY0255PASQLT			
	5 × 1,000 mL	FBY1005CASQLT	FBY1005PASQLT			
Centrifuge tube on stopcock manifold	5 × 15 mL	PSCT0155CASQLT	PSCT0155PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Minisart® vent filter	Polycarbonate, polystyrene, high-density polyethylene (HDPE), silicone (Pt)
	5 × 60 mL	PTG0065CASQLT	PTG0065PASQLT			
Bottle on stopcock manifold	5 × 60 mL	PTG0065CASQLT	PTG0065PASQLT	AseptiQuik® S ⅝" × ¼" Length 6" Quickseal® 3"	Minisart® vent filter	Polycarbonate, high-density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), silicone (Pt)
	5 × 125 mL	PTG0125CASQLT	PTG0125PASQLT			
	5 × 250 mL	PTG0255CASQLT	PTG0255PASQLT			

*List does not include tubing type material

5-Gang Manifold With Purge Vessel for Sterile Welding

Container Type	Volume	Catalog Number	Container Inlet	Container Outlet	Container Material	Other Fluid Contact Materials*
Flexsafe® bags on stopcock manifold	5 × 20 mL	FLS0025TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	Benchmark™ for needle-free syringe	Polyethylene (S80)	Polycarbonate, high-density polyethylene (HDPE), polyethylene (S80), polyethylene, polypropylene (PP), silicone (Pt)
	5 × 50 mL	FLS0055TPLG				
	5 × 150 mL	FLS0155TPLG				
	5 × 250 mL	FLS0255TPLG				
	5 × 500 mL	FLS0505TPLG				
	5 × 1,000 mL	FLS1005TPLG				
Centrifuge tube on stopcock manifold	5 × 15 mL	PSCT0155TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	25 mm Minisart® vent filter	Polystyrene	Polycarbonate, high-density polyethylene (HDPE), polyethylene (S80), polypropylene (PP), silicone (Pt)
	5 × 50 mL	PPCT0055TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	25 mm Minisart® vent filter	Polypropylene	Polycarbonate, high-density polyethylene (HDPE), polyethylene (S80), polyethylene, polypropylene (PP), silicone (Pt)
Bottle on stopcock manifold	5 × 60 mL	PTG0065TPLG	Tuflux® TPE with Plug ⅝" × ¼" Length 12" Quickseal® 3"	25 mm Minisart® vent filter	Polyethylene terephthalate glycol-modified (PETG)**	Polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate, silicone (Pt), polypropylene, polyethylene (S80), polyethylene
	5 × 125 mL	PTG0125TPLG				
	5 × 250 mL	PTG0255TPLG				

*List does not include tubing type material

TakeOne® Accessories



	Catalog Number	Fluid Contact Materials*
Quickseal® cutter (small diameter)	QSCUTTERSD	—
Quickseal® silicone caps, ¼" OD (50/pack)	QS04CAPSILNT	—
TakeOne® tab actuator lock (5/pack)	TKOTCHD	—
TakeOne® 25 mm depth accessory 1 (30 mm Ingold® port length)	INGDEPTH01*	316L Stainless Steel
TakeOne® 25 mm depth accessory 2 (46 mm Ingold® port length)	INGDEPTH02*	316L Stainless Steel
TakeOne® 25 mm depth accessory 3 (52 mm Ingold® port length)	INGDEPTH03*	316L Stainless Steel
TakeOne® 25 mm depth accessory 4 (40 mm Ingold® port length)	INGDEPTH04*	316L Stainless Steel

*For use on threaded 25 mm ports. Please consult you Sartorius representative if you need assistance determining port length.

Process Interface Compatibility

Type	Size	Number of Sampling Lines	Requires Multi-use Accessory	Part Number
TC	1.5"	× 5	No	—
	2"	× 9	No	—
25 mm Port Ingold® Safety Port	30 mm	× 3*	Yes	INGDEPTH01
	40 mm	× 3*	Yes	INGDEPTH04
	46 mm	× 3*	Yes	INGDEPTH02
	52 mm	× 3*	Yes	INGDEPTH03

* Uses the 25 mm TakeOne® single-use sampling hub.

Fast and Efficient Delivery

Rapidly adapt your sampling. Our off-the-shelf products* mean that TakeOne® is an extension of your operation and long delivery schedules are no longer a bottleneck to your timely processes.

*Select items from our standard portfolio are maintained in stock.

Sartorius, Your Partner to Innovation

At Sartorius, we strive to provide you with innovative products and technologies, and help you successfully implement these in your complex and quality-critical biomanufacturing processes across the globe. Our customers' problems are the objectives we set to overcome and help pioneer the way, setting standards for single-use products in a reliable and economical way.

Source all your needs for process optimization, customized validation and regulatory support with Confidence® Services for TakeOne®. Confidence® provides the closest interpretation of current regulatory requirements and industry standards by working with you to define relevant test conditions for all your process components. Services include, but are not limited to:

- Extractables assessment | Leachable studies
- Toxicological assessment | Safety evaluation
- Physio-Chemical studies

To support you in your regulatory needs, specific documentation is available upon request:

- Validation guide
- Technical drawing
- Inspection guide
- Instructions for use

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
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

www.sartorius.com