



TakeOne® Flex

Plan for adaptability with the TakeOne® Flex aseptic sampling solution

Benefits

Quickly React, on Demand

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.

Easily Reduce, and Expand

Avoid overwhelming your inventory management program. Maintain just a few stock keeping units (SKU) and build tens or thousands of designs.



Product Information

Sartorius Stedim Biotech now offers the TakeOne® aseptic sampling device with aseptic connectors to provide configuration flexibility at your point-of-use.

Efficient Sampling with TakeOne®

- Save time and avoid handling needles – Sampling lines have already been pre-assembled in the body of the device which is ready for immediate installation on your vessel. The genderless AseptiQuik® connectors ensure you a fast and easy sampling container assembly.
- Keep safe during SIP – TakeOne® face plate is made of 316L stainless steel to guarantee full compatibility with your tank mount. The device has been validated to pass 10 SIP cycles.
- Rest assured with reliable performance batch-to-batch – Sample and product integrity are preserved. The unique design of TakeOne® integrates an aseptic chamber to ensure a closed fluid pathway with no risks of leakage. All individual sampling lines have been 100% leak tested.

TakeOne® Flex Unique Configuration Possibilities:

TakeOne® Flex has a vast array of different sample containers; Single-use bags | bottles | centrifuge tubes | needleless syringes | manifolds.

With just a few sample container types, create a multitude of different sampling configurations:

Container SKU Qty	1.5" – 5 lines	2" – 9 lines
2	20	54
4	125	714
6	461	5,004
8	1,286	24,309
10	3,002	92,377

Simply Select, and Connect



Select sample container



Connect

Quick, and Easy Disconnect

Each TakeOne® Flex sampling line assembly 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.



Fast, and Efficient Delivery

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

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.

* On selected items only. Please contact your local sales specialist.

Sampling Containers

Tailored to your needs, choose from:
Bags, bottles, syringe access sites and tubes.



Flexboy® Bags

- Easy handling
- 50 | 150 | 250 | 500 | 1000 mL



Celsius®-Pak Bags

- Freeze applications
- 30 | 100 mL



PETG Bottles

- Free standing
- 30 | 60 | 125 | 250 | 500 | 1000 mL



Centrifuge Tubes

- High accuracy
- 15 | 50 mL



BENCHMARK®

- High capacity sampling
- Optional check valve (for needle-free syringes)



Manifolds with purge bag

- Increase Capacity

Technical Data

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	
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 (device only)
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)	
Septum	Platinum-cured Silicone
Mounting Mechanism	316L Stainless Steel, electropolished (<15 µin Ra)
2 mm Cannula	316L Stainless Steel, electropolished (<15 µin Ra)
Tubing Material	Silicone(Pt) or C-Flex® (374)
AseptiQuik® S Connector	Polycarbonate, USP Class VI, ADCF Platinum-cured Silicone (seal)

Regulatory Compliance	
Silicone Elastomer for Closures and Tubing	USP 381 EP 3.1.9 21CFR177.2600
Biocompatibility	USP <88> Class VI
Bacterial Endotoxin	USP <85> (<0.125 EU/mL)
Particulate Matter	USP <788> (≤3 particles/mL >25 microns; ≤10 particles/mL >10 microns)

Ingold® is a registered trademark of Mettler-Toledo

AseptiQuik® is a registered trademark of Colder Products Company

C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation

Ordering Information

A minimum order quantity of 5 applies for each item.

TakeOne® Flex | Silicone(Pt)

Catalog Number	Device Size & Fitting	Sampling Lines	Tube Material	Fitting	Fluid Contact Materials
24TC03PASQLT01	1.5" Sanitary TC	3	Sil(Pt) 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate
24TC05PASQLT01	1.5" Sanitary TC	5	Sil(Pt) 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate
32TC05PASQLT01	2" Sanitary TC	5	Sil(Pt) 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate
32TC09PASQLT01	2" Sanitary TC	9	Sil(Pt) 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate
INGP03PASQLT01	25 mm port (Ingold®)	3	Sil(Pt) 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate

TakeOne® Flex | TPE

Catalog Number	Device Size & Fitting	Sampling Lines	Tube Material	Fitting	Fluid Contact Materials
24TC03CASQLT01	1.5" Sanitary TC	3	C-Flex® 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate, C-Flex® (374)
24TC05CASQLT01	1.5" Sanitary TC	5	C-Flex® 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate, C-Flex® (374)
32TC05CASQLT01	2" Sanitary TC	5	C-Flex® 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate, C-Flex® (374)
32TC09CASQLT01	2" Sanitary TC	9	C-Flex® 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate, C-Flex® (374)
INGP03CASQLT01	25 mm port (Ingold®)	3	C-Flex® 1/8" x 1/4" Length 6"	AseptiQuik® S	316L Stainless Steel, Silicone (Pt), Polycarbonate, C-Flex® (374)

Ingold® is a registered trademark of Mettler-Toledo

TakeOne® Flex | Sample Containers

All items are individually packed and gamma sterilized for your convenience.
Contact your local sales specialist for custom-made sample containers.

Single-use Bags

	Catalog Number TPE tube	Catalog Number Silicone(Pt) tube	Container inlet	Container outlet	Fluid Contact Materials
Flexboy® bag 50 mL	FBY0051CASQLT	FBY0051PASQLT	AseptiQuik® S 1/8" x 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
Flexboy® bag 150 mL	FBY0151CASQLT	FBY0151PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
Flexboy® bag 250 mL	FBY0251CASQLT	FBY0251PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
Flexboy® bag 500 mL	FBY0501CASQLT	FBY0501PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
Flexboy® bag 1000 mL	FBY1001CASQLT	FBY1001PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, ethylene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
Celsius® -Pak bag 30 mL	CSP0031CASQLT	-	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	Luer fitting	Polycarbonate, ethylene-vinyl acetate (EVA S71 - 360 µm), polypropylene (PP), silicone(Pt)
Celsius® -Pak bag 100 mL	CSP0101CASQLT	-	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	Luer fitting	Polycarbonate, ethylene-vinyl acetate (EVA S71 - 360 µm), polypropylene (PP), silicone(Pt)

BENCHMARK® (for needle-free syringes)

	Catalog Number TPE tube	Catalog Number Silicone(Pt) tube	Container inlet	Container outlet	Fluid Contact Materials
BENCHMARK® with checkvalve	BNCHCV1CASQLT	BNCHCV1PASQLT	AseptiQuik® S 1/8" x 1/4" Length 6" QUICKSEAL® 3"	–	Polycarbonate, Polypropylene (PP), Styrene acrylonitrile (SAN), silicone(Pt)
BENCHMARK® w/out checkvalve	BNCH1CASQLT	BNCH1PASQLT	AseptiQuik® S 1/8" x 1/4" Length 6" QUICKSEAL® 3"	–	Polycarbonate, silicone(Pt)

Bottles

	Catalog Number TPE tube	Catalog Number Silicone(Pt) tube	Container inlet	Container outlet	Fluid Contact Materials
PETG Bottle 30 mL	PTG0031CASQLT	PTG0031PASQLT	AseptiQuik® S 1/8" x 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polyethylene terephthalate glycol-modified (PETG), silicone(Pt)
PETG Bottle 60 mL	PTG0061CASQLT	PTG0061PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polyethylene terephthalate glycol-modified (PETG), silicone(Pt)
PETG Bottle 125 mL	PTG0121CASQLT	PTG0121PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polyethylene terephthalate glycol-modified (PETG), silicone(Pt)
PETG Bottle 250 mL	PTG0251CASQLT	PTG0251PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polyethylene terephthalate glycol-modified (PETG), silicone(Pt)
PETG Bottle 500 mL	PTG0501CASQLT	PTG0501PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polyethylene terephthalate glycol-modified (PETG), silicone(Pt)
PETG Bottle 1000 mL	PTG1001CASQLT	PTG1001PASQLT	AseptiQuik® S 8" x 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polyethylene terephthalate glycol-modified (PETG), silicone(Pt)

Centrifuge tubes

	Catalog Number TPE tube	Catalog Number Silicone(Pt) tube	Container inlet	Container outlet	Fluid Contact Materials
Polystyrene Centrifuge Tube 15 mL	PSCT0151CASQLT	PSCT0151PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polystyrene, high density polyethylene (HDPE), silicone(Pt)
Polypropylene Centrifuge Tube 50 mL	PPCT0051CASQLT	PPCT0051PASQLT	AseptiQuik® S 8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polypropylene (PP), silicone(Pt)

Standard 2-gang manifolds (with purge vessel)

	Catalog Number TPE tube	Catalog Number Silicone(Pt) tube	Container inlet	Container outlet	Fluid Contact Materials
2-way splitter	2WAY001CASQLT	2WAY001PASQLT	AseptiQuik® S (× 3) 1/8" × 1/4" Length 6"	–	Polycarbonate, high density polyethylene (HDPE), silicone(Pt)
2 × 50 mL Flexboy® bags on stopcock manifold	FBY0052CASQLT	FBY0052PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), ethyl- ene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
2 × 250 mL Flexboy® bags on stopcock manifold	FBY0252CASQLT	FBY0252PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), ethyl- ene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
2 × 1000 mL Flexboy® bags on stopcock manifold	FBY1002CASQLT	FBY1002PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), ethyl- ene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
2 × 15 mL centrifuge tube on stopcock manifold	PSCT0152CASQLT	PSCT0152PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polystyrene, high density polyethylene (HDPE), silicone(Pt)
2 × 50 mL centrifuge tube on stopcock manifold	PPCT0052CASQLT	PPCT0052PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, high density polyethylene (HDPE), poly- propylene (PP), silicone(Pt)
2 × 60 mL PETG Bottle on stopcock manifold	PTG0062CASQLT	PTG0062PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), silicone(Pt)

Standard 5-gang manifolds (with purge vessel)

	Catalog Number TPE tube	Catalog Number Silicone(Pt) tube	Container inlet	Container outlet	Fluid Contact Materials
5 × 50 mL Flexboy® bags on stopcock manifold	FBY0055CASQLT	FBY0055PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), ethyl- ene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
5 × 250 mL Flexboy® bags on stopcock manifold	FBY0255CASQLT	FBY0255PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), ethyl- ene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
5 × 1000 mL Flexboy® bags on stopcock manifold	FBY1005CASQLT	FBY1005PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	BENCHMARK® (for needle-free syringes)	Polycarbonate, high density polyethylene (HDPE), ethyl- ene-vinyl acetate (EVA S71), polypropylene (PP), silicone(Pt)
5 × 15 mL centrifuge tube on stopcock manifold	PSCT0155CASQLT	PSCT0155PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, polystyrene, high density polyethylene (HDPE), silicone(Pt)
5 × 50 mL centrifuge tube on stopcock manifold	PPCT0055CASQLT	PPCT0055PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, high density polyethylene (HDPE), poly- propylene (PP), silicone(Pt)
5 × 60 mL PETG Bottle on stopcock manifold	PTG0065CASQLT	PTG0065PASQLT	AseptiQuik® S 1/8" × 1/4" Length 6" QUICKSEAL® 3"	MiniSart® Vent Filter	Polycarbonate, high density polyethylene (HDPE), polyethylene terephthalate glycol-modified (PETG), silicone(Pt)

TakeOne® Flex | Accessories

Catalog Number	Description
QSCUTTERSD	QUICKSEAL® Cutter (small diameter)
QS04CAPSILNT	QUICKSEAL® Silicone caps, 1/4" OD (50/pack)
TKOTCHD	TakeOne® tab actuator lock (5/pack)
INGDEPTH01	TakeOne® 25 mm depth accessory 1 (30 mm Ingold® port length)
INGDEPTH02	TakeOne® 25 mm depth accessory 2 (46 mm Ingold® port length)
INGDEPTH03	TakeOne® 25 mm depth accessory 3 (52 mm Ingold® port length)
INGDEPTH04	TakeOne® 25 mm depth accessory 3 (40 mm Ingold® port length)

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
Phone +49.551.308.0
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
Argentina +54.11.4721.0505
Brazil +55.11.4362.8900
Mexico +52.55.5562.1102
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.913.586.098
Russian Federation +7.812.327.53.27
Japan +81.3.4331.4300
China +86.21.6878.2300

Specifications subject to change without notice.
Copyright Sartorius Stedim Biotech GmbH. Printed
in the EU on paper bleached without chlorine.
Order No.: 1000055879
Version 03
Date: 2019/03/06