

Takeone®

Design Services for Aseptic Sampling Solutions

Product Information

Have a particular need and can't find a standard solution? Sartorius offers different design services to fine-tune your sampling plan to your exact requirements.

Configure-to-Order

Standard features adjusted to your needs

Takeone® is configured from a database of pre-qualified components with proven functionalities. This provides the flexibility needed for an application-based single-use configuration whilst improving design and quotation turnaround times, manufacturing lead times and security of supply when compared with fully unique engineered custom designs.



Engineer-to-Order

Customize Takeone® design to match your unique requirements




This service allows the selection of a larger variety of solutions with added features, such as alternative tube materials or collection vessels, autoclavable options and more. If needed, entirely new components can be investigated with the potential to establish special supply terms with the supplier. In addition, you will be allocated a unique part number, so that the design can evolve together with your process without the hassle of changing references.



The robustness of such custom solutions relies on a complete chain of engineering, qualification, application development and validation services. Contact one of our specialists to initiate a project with a dedicated team of expert engineers to help you design your Takeone® aseptic sampling device.

Designing Takeone®

Configure your Takeone® from our library of common components and solutions in today's market:

| Tank fitting | 25 mm port (Ingold®) | 1.5" Sanitary TC | 2" Sanitary TC |
|----------------------------|---|---|--|
| |  |  |  |
| Quantity of sampling lines | 1-3 | 1-5 | 1-9 |
| Tubing material | Silicone (Pt) C-Flex® | | |
| Tubing length | 6" 12" with Quickseal® aseptic disconnect | | |
| Samples | Individual containers Flexboy® Bags: 50 150 250 500 1000 mL Celsius® Pak Bags: 30 100 mL PETG Bottles: 30 60 125 250 500 1000 mL Centrifuge Tubes: 15 mL (polystyrene) 50 mL (polypropylene) Benchmark® (access for needle-free syringes) | | Manifolds All of our individual containers can be manifolded together in 2-6 gang with a 500 mL purge bag. |

Our Experts Are Available to Help You

Whether it's questions you may have or the development of a completely new design, Sartorius is committed to helping you successfully achieve your sampling objectives. Contact your local sales specialist to start your personalized sampling plan solution.

Ingold® is a registered trademark of Mettler-Toledo.
C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation.

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 |
| 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 |
| Equipment interface | 316L Stainless Steel, electropolished (<15 µm Ra) |
| 2 mm cannula | 316L Stainless Steel, electropolished (<15 µm Ra) |
| Tubing material | Silicone(Pt) or C-Flex® (374) |

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) |

Ordering Information

All items are individually packed and gamma sterilized for your convenience.

| | Lead Time | Minimum Order Quantity |
|--------------------|-----------|------------------------|
| Configure-to-Order | 10 weeks | 5 |
| Engineer-to-Order | 12 weeks | 10 |




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 further contacts, visit
www.sartorius.com

Specifications subject to change without notice.

© 2021 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France

Order No.: 1000055880-e | Status: 09 | 30 | 2021