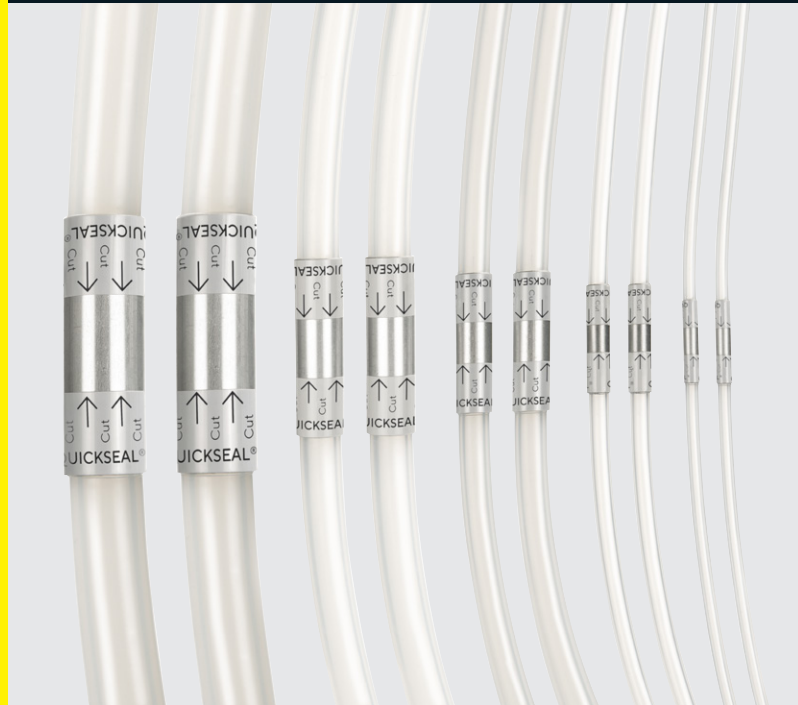


Quickseal[®]

Aseptic Disconnections



Product Information

Installed on an array of tube materials including platinum-cured and thermoplastic elastomers such as C-Flex[®] and Tuflux[®] TPE, Quickseal[®] is the quickest aseptic tube-separation technique available today. Extensively tested, Quickseal[®] is proven aseptic and the seal stronger than the tubing it is part of.

Applications

Quickseal[®] aseptic disconnections are used after fluid transfer to disconnect single-use transfer lines and bag or bottle assemblies used in biopharmaceutical applications. Quickseal[®] aseptic disconnections provide aseptic disconnection in non-classified and classified environments. It can be applied to multiple types and sizes of tubing.

Description

Quickseal® can be installed on a variety of tube materials across a range of tube sizes. The tubing runs continuously through the collar for an uninterrupted fluid path.

The Quickseal® collar is bonded to the outside of the tubing. The cut collar holds the walls of the tubing together to create the aseptic seal.

Quickseal® can be included as a component on single-use tube, bag or bottle assemblies, sampling manifolds, single-use bioreactors and more.

Quickseal® reduces operating time and improves reliability for biopharmaceutical and other high-purity fluid handling industries.



Collar is uncut

Collar has been cut and filter removed

Technical Data

Specifications

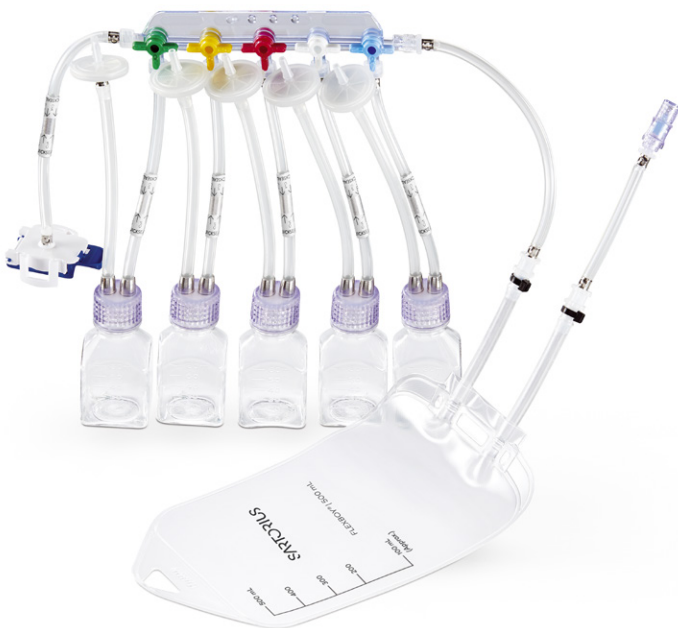
Quickseal® Aseptic Disconnections

Tubing Dimensions	
OD inch	1/8" 1/4" 3/8" 7/16" 1/2" 5/8" 3/4" 1" 1 1/8"
OD mm	3.2 6.4 9.6 11.1 12.7 15.9 19.0 25.4 28.6
Tubing Material	
	Silicone (Liveo PHARMA® 50 STHT-C®, Pumpsil®) TPE (C-Flex®, Tuflux® TPE) Ask about other tube materials
Quickseal® Position	Variable
Tubing Length	
inch	6" 12" 18" 24" 36"
mm	~150 305 460 610 915 Other tube lengths available on request
Sealing Time	< 4 seconds
Operating Steps	1
Seal Type	Mechanical
Fluid Contact Material	Tubing Material
Component Material	Aluminum (3003)
Validation Testing	
	Container Closure (Ambient) Container Closure (Freeze Thaw) Burst Test Tensile Pull Test



1/8" | 3.2 mm

1 1/8" | 28.6 mm



Quickseal® collars on inlet line to each Mycap® bottle assembly

Please note:

Mycap® and Tuflux® TPE are registered trademarks of Sartorius Stedim Biotech C-Flex® and STHT-C® is a registered trademark of St. Gobain Liveo. Pharma® is a registered trademark of Dupont Polymers Inc.

Cutting Tool Specifications: Large Diameter Cutter

Materials of Construction (cutting head)	304 Stainless Steel
Weight	4 lbs (2.2 kg)
Width	2" (60 mm)
Depth	4 ¼" (108 mm)
Length	14 ¾" (375 mm)
Sound	< 80 dB at 1 meter
Vibration	< 2.5 m/s ²
Cutting Time	4 seconds
Closing Speed	6 mm/s
Cuts per Charge	≈300
Battery Voltage	18V
Charging Time	20 min



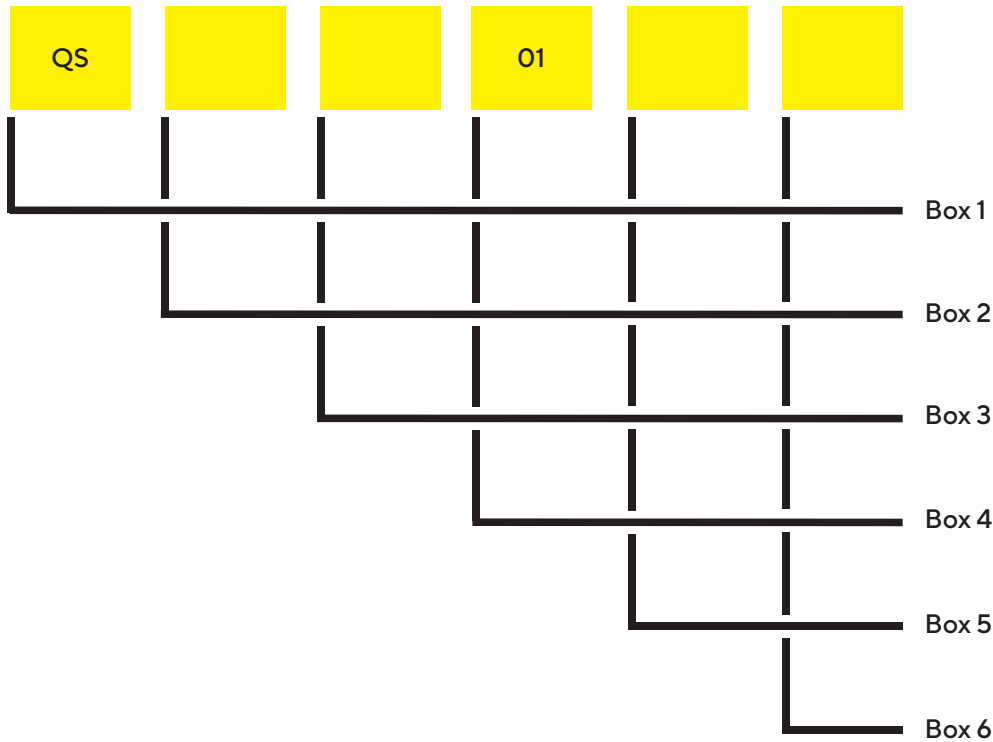
Cutting Tool Specifications: Small Diameter Cutter

Quickseal® Collar Size Range	< ¼" (6.4 mm)
Materials of Construction	420 Stainless Steel
Weight	¾ lbs (345g)
Width	½" (13 mm)
Depth	1 ½" (27 mm)
Length	8 ½" (206 mm)



Ordering Information

Quickseal® can be included on a Sartorius single-use assembly or can be ordered as standalone component. Build the part number according to the instructions below to order Quickseal® as a standalone component.



Explanation

Box 1

Fixed Value(QS). Indicates Quickseal® product family.

Box 2

Indicates tubing material

P Platinum-cured silicone. Pumpsil® when 1/8" ID x 1/4" OD, STHT-C otherwise

PG Platinum-cured silicone. STHT-C when 1/8" ID x 1/4" OD

PD Platinum-cured silicone. Liveo® Pharma 50

C Thermoplastic Elastomer. C-Flex®

T Thermoplastic Elastomer. Tuflux® TPE inquire for other tubing materials.

Box 3

Indicates internal diameter (ID) and outer diameter (OD) of tubing

0204 1/8" ID x 1/4" OD

0406 1/4" ID x 3/8" OD

0407 1/4" ID x 7/8" OD

0408 1/4" ID x 1/2" OD

0610 3/8" ID x 5/8" OD

0812 1/2" ID x 3/4" OD

1216 3/4" ID x 1" OD

1218 3/4" ID x 1-1/8" OD

Box 4

Fixed Value(01). Indicates number of collars on tube.

Box 5

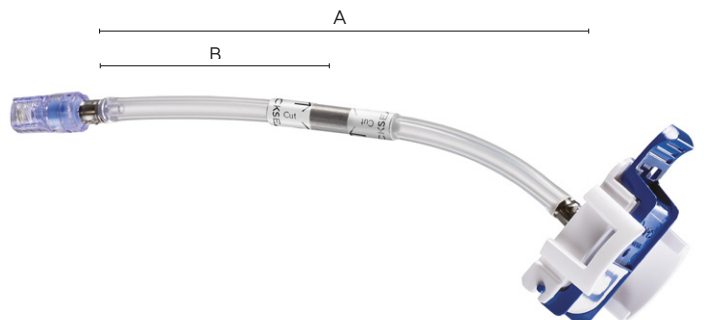
Length from edge of tubing to center of Quickseal® collar. Dimension B on illustration.

Box 6

Overall Length (inches). Dimension A on illustration. Selections available as 6 inch increments.

Example:

QST0204010612 is 1/8" ID x 1/4" OD Tuflux® TPE, 12" overall length with Quickseal® centered at 6"



Quickseal® Protective Caps

The Quickseal® protective caps can be installed after collars are cut. The caps are made from flexible and durable silicone. They are suitable for a wide range of temperatures.

Installation is simple:

1. Grasp the cap near the open end between index finger and thumb
2. Squeeze the cap until the opening is wide enough for the cut Quickseal® collar
3. Insert the Quickseal® collar and release



Order Code	Description
QS04CAPSILNT	Silicone protection caps for ¼" (6.4 mm) OD Quickseal® collars (pack of 50)
QS06CAPSILNT	Silicone protection caps for ⅜" (9.6 mm) OD Quickseal® collars (pack of 50)
QS07CAPSILNT	Silicone protection caps for ⅞" (11.1 mm) OD Quickseal® collars (pack of 50)
QS08CAPSILNT	Silicone protective caps for ½" (12.7 mm) OD Quickseal® collars (pack of 50)
QS10CAPSILNT	Silicone protection caps for ⅝" (16 mm) OD Quickseal® collars (pack of 50)
QS12CAPSILNT	Silicone protection caps for ¾" (19.0 mm) OD Quickseal® collars (pack of 50)
QS18CAPSILNT	Silicone protection caps for 1.125" (28.5 mm) OD Quickseal® collars (pack of 50)

Quickseal® Cutting Tools

The Quickseal® cutting tools were designed for single-person operation with safety, simplicity and portability in mind. Place the collar in the cutting and cut the Quickseal® collar.




Order Code	Description
QSCUTTERSD	Small diameter Quickseal® cutter for tubes up to 0.25" (6.4 mm) OD
QSCUTTERLD	Large diameter Quickseal® cutter for tubes up to 1.125" (28.6 mm) OD; 120V battery charger and two 18V batteries
QSCUTTERLD230V	Large diameter Quickseal® cutter for tubes up to 1.125" (28.6 mm) OD; 230V battery charger and two 18V batteries
QSCUTTERREADER	Electronic reader for battery-powered large diameter Quickseal® cutter

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

Publication No.: SP-1512-e | Order No.: 85037-556-32 | Status: 10 | 01 | 2021