

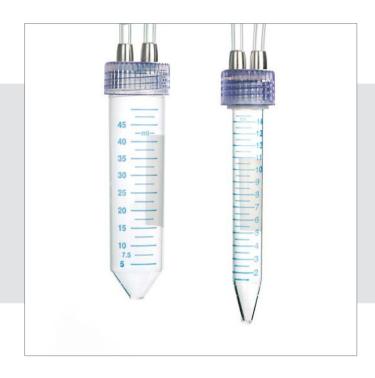
# MYCAP™ Centrifuge Tubes

# **Product Information**

MYCAP™ bottle closures are a one-piece, ready-to-use aseptic fluid transfer system.

The single-use design features integral tubing and a robust platinum-cured silicone seal.

MYCAP™ bottle assemblies include Falcon® 15mL and 50mL centrifuge tubes, are gamma irradiated and ready for immediate use.



Features	Benefits
– Ready-to-use	<ul><li>Save time</li><li>Eliminate errors</li></ul>
<ul><li>Single-use</li></ul>	<ul> <li>No contamination risk</li> </ul>
- One-For-All silicone closure	<ul> <li>Universal validation</li> <li>Robust &amp; reliable</li> <li>performance</li> </ul>
<ul> <li>One-piece construction</li> </ul>	<ul> <li>Integral closure</li> <li>Easy to handle</li> </ul>
<ul> <li>Easy customization</li> </ul>	<ul> <li>Configure MYCAP™ to your process needs</li> </ul>
<ul><li>One-For-All silicone closure</li><li>One-piece construction</li></ul>	<ul> <li>Universal validation Robust &amp; reliable performance</li> <li>Integral closure Easy to handle</li> <li>Configure MYCAP™</li> </ul>

#### **Applications**

MYCAP™ is ideal for critical fluid-handling applications.

 $\mathsf{MYCAP}^\mathsf{TM}$  bottle assemblies are gamma irradiated and ready for immediate use for:

- Buffer and Media Storage
- Buffer and Media Addition
- Bioreactor Feeds
- Sample Collection
- Bulk Harvest
- Product Pooling
- Bulk Intermediate Hold

Reduce Risk & Improve Performance

With  $\mathsf{MYCAP}^\mathsf{TM}$  bottle closures, every use is its first use.

There is no risk of batch-to-batch or product-to-product cross contamination.

MYCAP™ bottle assemblies pass studies of container closure by bacterial immersion.

#### One-For-All Closure System

At last - a truly universal bottle closure system.

All MYCAP™ closure systems are constructed using the same high performance platinum-cured silicone seal regardless of the tubing material, tubing size, port configuration, bottle type or bottle size.

Ease validation and reduce regulatory work by implementing MYCAP™ throughout the production process.

#### **Easy Implementation**

Alternative bottle closure systems include a number of reusable components which are costly to implement and manage.

Regulatory agencies demand a scientifically sound validation strategy for parts cleaning to avoid non-host or batch-to-batch contamination. The physical condition and life cycle of reused parts must be closely monitored to ensure cleaning efficacy.

Select pre-assembled and single-use MYCAP™ bottle assemblies to bypass the burden of validation and parts management.

#### Reliable Performance

The platinum-cured silicone seal of MYCAP $^{\text{m}}$  is suitable for a range of applications.

The closure to the bottle is secure so long as the seal maintains its physical properties. Silicone maintains its elastomeric properties from -70° to 260° Celcius.

Validation demonstrates a leak-free seal when MYCAP™ are installed with torqueforce from 4in-lbs to 60in-lbs.

#### Technical Data

#### **Specifications**

Elastomer Seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex® formulation 374
Cap Material	Polycarbonate (not fluid contact)
Bottle Style	Falcon® Polystyrene and Polypropylene Centrifuge Tube
Tube Ports	Inlet   Outlet, Air vent - 25 mm PES membrane MiniSart®
Fittings	Luer® Lock or tube plug
Sterilization	Gamma Irradiated

#### Manufacturing Environment

MYCAP™ bottle closures are constructed and packaged in an ISO 7 clean room.

Monthly cleaning with a schedule of LpH®, Vesphene® and Spor Klenz® reduce and control viable organisms. Viable organisms are measured quarterly to monitor the effectiveness of the clean room management and cleaning procedures and to be compliant to EU GMPs and ISO14698.

Line clearance, weekly cleaning of equipment and work surfaces and monthly cleaning of the cleanroom reduce and control non-viable particles. Non-viable readings are taken weekly to ensure  $0.5 \ \mu m/m^3$  and  $5.0 \ \mu m/m^3$  particles are within the ISO Class 7 acceptance criteria, per ISO 14644-1.

#### **Quality Assurance**

MYCAP™ bottle assemblies are manufactured in accordance with the key principles of cGMP.

Each MYCAP™ bottle closure is issued with a unique lot number on its' Certificate of Compliance. The lot number provides full traceability to all components used and operations performed.

The Certificate of Conformance for MYCAP™ bottle assemblies reports compliance to:

- USP <87>: Biological reactivity tests, in Vitro
- USP <88>: Biological reactivity tests, in Vivo
- USP <788>: Particulate matter
- USP <85>: Endotoxin
- Sterility Assurance (SAL) 10<sup>-6</sup> per ISO11137

#### **Shelf Life**

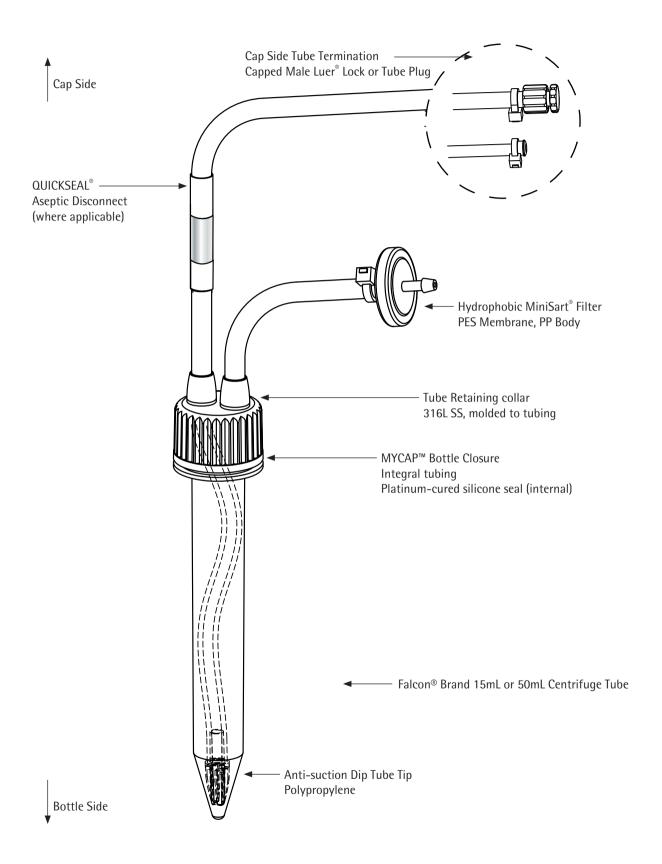
MYCAP™ bottle closures, components and component resins have shelf-life that exceeds 2 years.

#### TSE BSE Risk

The components and materials used in construction of MYCAP™ bottles assemblies are treated at high temperatures and undergo chemical reactions which far exceed the stringent requirements laid down in the documents CPMP|BWP|1230|98 and WHO|CPS|VPH|95.145 and or EMA|410|01, rev 3. These processing conditions are considered sufficient to inactivate TSE ft BSE transmitters.

C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation Luer® is a registered trademark of Becton Dickinson Dow Corning® is a registered trademark of Dow Corning Corporation LpH®, Vesphene® Spor Klenz® are registered trademarks of STERIS Corporation Nalgene® is a registered trademark of ThermoFisher Scientific

# MYCAP™ Bottle Assembly Anatomy



# **Ordering Information**

# Series A MYCAP™ Bottle Assemblies

- Optimized Tube Lengths including dip tube for fluid transfer
- Includes Falcon® Brand 15mL Polystyrene or 50mL Polypropylene Centrifuge Tube
- Platinum-Cured Silicone Tubing (Dow-Corning® PHARMA-50)
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	<b>Bottle Size</b>	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015A	2 port MYCAP™ bottle closure with 15mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 7.25"; capped male  Luer® fitting <b>Bottle Side:</b> To Bottom;  dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart <sup>®</sup> hydrophobic PES filter Bottle Side: None	1
MYCAPCPPG0050A	2 port MYCAP™ bottle closure with 50mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 7.5"; capped male Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart <sup>®</sup> hydrophobic PES filter Bottle Side: None	1

# **Series B MYCAP™ Bottle Closures**

- Optimized Tube Lengths including dip tube for fluid transfer
- Includes Falcon® Brand 15mL Polystyrene or 50mL Polypropylene Centrifuge Tube
- C-Flex® tubing (formulation 374)
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	<b>Bottle Size</b>	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015B	2 port MYCAP™ bottle closure with 15mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 13.25"; capped male  Luer® fitting <b>Bottle Side:</b> To Bottom;  dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart <sup>®</sup> hydrophobic PES filter Bottle Side: None	1
MYCAPCPPG0050B	2 port MYCAP™ bottle closure with 50mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 13.5"; capped male Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart <sup>®</sup> hydrophobic PES filter Bottle Side: None	1

# Series C MYCAP™ Bottle Assemblies

- Optimized Tube Lengths for including dip tube for fluid transfer
- Includes Falcon® Brand 15mL Polystyrene or 50mL Polypropylene Centrifuge Tube
- Platinum-Cured Silicone Tubing (Dow-Corning® PHARMA-50)
- QUICKSEAL® aseptic disconnect
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	<b>Bottle Size</b>	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015C	2 port MYCAP™ bottle closure with 15mL centrifuge tube	5 mL	Tube Size: $\frac{1}{8}$ " ID $\times \frac{1}{4}$ " OD Cap Side: 7.25"; QUICKSEAL® at 2.25"; capped male Luer® fitting Bottle Side: To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPCPPG0050C	2 port MYCAP™ bottle closure with 50mL centrifuge tube	50 mL	Tube Size: $\frac{1}{8}$ " ID $\times$ $\frac{1}{4}$ " OD Cap Side: 7.5"; QUICKSEAL® at 2.5"; capped male Luer® fitting Bottle Side: To Bottom; dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart® hydrophobic PES filter Bottle Side: None	1

# Series D MYCAP™ Bottle Assemblies

- Optimized Tube Lengths for including dip tube for fluid transfer
- Includes Falcon® Brand 15mL Polystyrene or 50mL Polypropylene Centrifuge Tube
- C-Flex tubing (formulation 374)
- QUICKSEAL® aseptic disconnect
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	<b>Bottle Size</b>	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015D	2 port MYCAP™ bottle closure with 15 mL centrifuge tube	15 mL	Tube Size: 1/8" ID × 1/4" OD Cap Side: 13.25"; QUICKSEAL® at 2.25"; plug at end Bottle Side: To Bottom; dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart <sup>®</sup> hydrophobic PES filter Bottle Side: None	1
MYCAPCPPG0050D	2 port MYCAP™ bottle closure with 50mL centrifuge tube	50 mL	Tube Size: $\frac{1}{8}$ " ID $\times$ $\frac{1}{4}$ " OD Cap Side: 13.5"; QUICKSEAL® at 2.5"; plug at end Bottle Side: To Bottom; dip tube tip	Tube Size: 1/8" ID × 1/4" OD Cap Side: 3"; Minisart <sup>®</sup> hydrophobic PES filter Bottle Side: None	1

#### **Sales and Service Contacts**

For further contacts, visit www.sartorius.com

#### Europe

**Germany** Sartorius Lab Instruments GmbH & Co. KG Weender Landstrasse 94–108 37075 Goettingen

Phone +49.551.308.0 Fax +49.551.308.3289

France & Suisse Romande

Sartorius France 2, rue Antoine Laurent de Lavoisier 7A de la Gaudrée 91410 Dourdan

Phone +33.1.70.62.50.00 Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH Modecenterstrasse 22 1030 Vienna

Phone +43.1.7965760.0 Fax +43.1.7965760.24

Sartorius Belgium N.V. Rue Colonel Bourg 105 1030 Bruxelles

Phone +32.2.756.06.90 Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy 00880 Helsinki

Phone +358.9.755.951 Fax +358.9.755.95.200

**Hungary** Sartorius Hungária Kft. Kagyló u. 5. 2092 Budakeszi

Phone +3623.457.227 Fax +3623.457.147

Ireland

Sartorius Ireland Ltd. Unit 41, The Business Centre Stadium Business Park Ballycoolin Road

Phone +353.1.8089050 Fax +353.1.8089388

Italy

Sartorius Italy S.r.l. Viale A. Casati, 4 20835 Muggiò (MB)

Phone +39.039.4659.1 Fax +39.039.4659.88

Netherlands

Sartorius Netherlands B.V. Phone +31 30 60 53 001

Fax +31.30.60.52.917

info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o. ul. Wrzesinska 70 62-025 Kostrzyn

Phone +48.61.6473830 Fax +48.61.6473839

Russian Federation

LLC "Sartorius RUS" Uralskava str. 4. Lit. B 199155 St. Petersburg

Phone +7.812.327.53.27 Fax +7.812.327.53.23

Spain & Portugal

Sartorius Spain, S.A. Avda. de la Industria, 32 Edificio PAYMA 28108 Alcobendas (Madrid)

Phone Spain +34.902.123.367 Phone Portugal +351.800.855.800 Fax Spain +34.91.358.96.23 Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG Ringstrasse 24a 8317 Tagelswangen (ZH)

Phone +41.44.746.50.00 Fax +41.44.746.50.50

II.K.

Sartorius UK Ltd. Longmead Business Centre Blenheim Road, Epsom Surrey KT19 9QQ

Phone +44.1372.737159 Fax +44.1372.726171

Ukraine

LLC "Biohit" Post Box 440 "B" 01001 Kiev, Ukraine

Phone +380.44.411.4918 Fax +380.50.623.3162

**Americas** 

IISA

Sartorius Corporation 5 Orville Drive, Suite 200 Bohemia, NY 11716

Phone +1.631.254.4249 Toll-free +1.800.635.2906 Fax +1.631.254.4253

Argentina Sartorius Argentina S.A. Int. A. Ávalos 4251 B1605ECS Munro Buenos Aires

Phone +54.11.4721.0505 Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda Avenida Senador Verqueiro 2962 São Bernardo do Campo CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900 Fax +55.11.4362.8901

Canada

Sartorius Canada Inc. 2179 Dunwin Drive #4 Mississauga, ON L5L 1X2

Phone +1.905.569.7977 Toll-Free +1.800.668.4234 Fax +1.905.569.7021

Sartorius de México S A de C V Circuito Circunvalación Poniente No. 149 Ciudad Satélite 53100, Estado de México México

Phone +52.5555.62.1102 Fax +52.5555.62.2942

Peru

Sartorius Peru S.A.C. Av. Emilio Cavenecia 264 San Isidro 15073 Lima, Perú

Phone +51.1.441 0158 Fax +51 1 422 6100

# Asia | Pacific

Sartorius Australia Pty. Ltd. Unit 5, 7-11 Rodeo Drive Dandenong South Vic 3175

Phone +61.3.8762.1800 Fax +61.3.8762.1828

Sartorius (Shanghai) Trading Co., Ltd. 3rd Floor, North Wing, Tower 1 No. 4560 Jinke Road Zhangjiang Hi-Tech Park Pudong District Shanghai 201210, P.R. China

Phone +86.21.6878.2300 Fax +86.21.6878.2882

**Hong Kong** Sartorius Hong Kong Ltd. Unit 1012, Lu Plaza 2 Wing Yip Street Kwun Tong Kowloon, Hong Kong

Phone +852.2774.2678 Fax +852.2766.3526

Sartorius Weighing India Pvt. Ltd. #69/2-69/3, NH 48, Jakkasandra, Nelamangala Tq 562 123 Bangalore, India

Phone +91.80.4350.5250 Fax +91.80.4350.5253

**Japan** Sartorius Japan K.K. 4th Fl., Daiwa Shinagawa North Bldg. 8–11, Kita-Shinagawa 1-chome Shinagawa-ku, Tokyo, 140–0001 Japan

Phone +81,3,3740.5408 Fax +81.3.3740.5406

Malaysia Sartorius Malaysia Sdn. Bhd Lot L3-E-3B, Enterprise 4 Technology Park Malaysia 57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622 Fax +60.3.8996.0755

Singapore Sartorius Singapore Pte. Ltd 1 Science Park Road, The Capricorn, #05-08A, Singapore Science Park II Singapore 117528

Phone +65 6872 3966 Fax +65.6778.2494

South Korea

Sartorius Korea Ltd. 8th Floor, Solid Space B/D, PanGyoYeok-Ro 220, BunDang-Gu SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700 Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd. 129 Rama 9 Road, Huaykwang Bangkok 10310

Phone +66.2643.8361-6 Fax +66.2643.8367