

# Mycap<sup>®</sup> Bottle Closures



## Product Information

Mycap<sup>®</sup> 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.

Available non-sterile for installation on autoclavable glass bottles, or fully integrated to the container, presterilized and ready for use.

# Table of Contents

<b>Mycap® GL45 Bottle Closures</b>	<b>4</b>
Applications	4
Technical Data	4
Image of Anatomy	5
Order Information	6
<b>Mycap® Bottle Assemblies</b>	<b>12</b>
Applications	12
Technical Data	12
Image of Anatomy	13
Order Information	14
<b>Mycap® Centrifuge Tubes</b>	<b>18</b>
Applications	18
Technical Data	18
Image of Anatomy	19
Order Information	20
<b>Mycap® Single-Bottle Assemblies for Media</b>	<b>23</b>
Applications	23
Technical Data	23
Image of Anatomy	24
Order Information	25
<b>Mycap® Manifold Bottle Assemblies for Media</b>	<b>29</b>
Applications	29
Technical Data	29
Image of Anatomy	30
Order Information	31

# Introduction

Features	Benefits
Ready-to-use	Save time Eliminate errors
Single-use	No contamination risk
One-for-all silicone closure	Universal validation Robust and reliable performance
One-piece construction	Integral closure Easy to handle
Easy customization	Configure Mycap® to your process needs

## Reduce Risk and Improve Performance

With Mycap® bottle closures, every use is its first use. There is no risk of batch-to-batch or product-to-product cross contamination. Mycap® bottle closures 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 closures to bypass the burden of validation and parts management.

## Reliable Performance

The platinum-cured silicone seal of Mycap® 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° Celsius.

GL45 Mycap® bottle closures can be field installed without use of any instruments or tools. A survey of random operators found the average torque applied by hand yields a leak-free closure. Validation demonstrates a leak-free seal when Mycap® are installed with torque force from 4in-lbs to 60in-lbs.

## 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 ISO 14698.

Line clearance, weekly cleaning of equipment and work surfaces and monthly cleaning of the clean-room reduce and control non-viable particles. Non-viable readings are taken weekly to ensure 0.5 µm/m<sup>3</sup> and 5.0 µm/m<sup>3</sup> particles are within the ISO Class 7 acceptance criteria, per ISO 14644-1.

## Quality Assurance

Mycap® bottle closures 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 GL45 Mycap® bottle closures reports compliance to:

- USP <87>: Biological reactivity tests, in Vitro
- USP <88>: Biological reactivity tests, in Vivo
- USP <788>: Particulate matter
- USP <85>: Endotoxin

## Shelf Life

Mycap® bottle closures, components and component resins have shelf-life that exceeds two years.

## TSE | BSE Risk

The components and materials used in construction of Mycap® bottles closures 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 and BSE transmitters.

# Mycap<sup>®</sup> GL45 Bottle Closures

## Applications

Mycap<sup>®</sup> is ideal for critical fluid-handling applications. Attach to glass bottles with GL45 threaded closures, autoclave and use for:

- Buffer and media storage
- Buffer and media addition
- Bioreactor feeds
- Sample collection
- Bulk harvest
- Product pooling
- Bulk intermediate hold

## Technical Data

### Specifications

Elastomer seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex <sup>®</sup> formulation 374 or Tuflux <sup>®</sup> TPE
Cap material	Polycarbonate (not fluid contact)
Closure style	GL45
Tube ports	Inlet   Outlet, Air vent - 25 mm PES membrane Minisart <sup>®</sup>
Fittings	Luer <sup>®</sup> lock or tube plug
Sterilization	Autoclavable

C-Flex<sup>®</sup> is a registered trademark of Saint-Gobain Performance Plastics Corporation

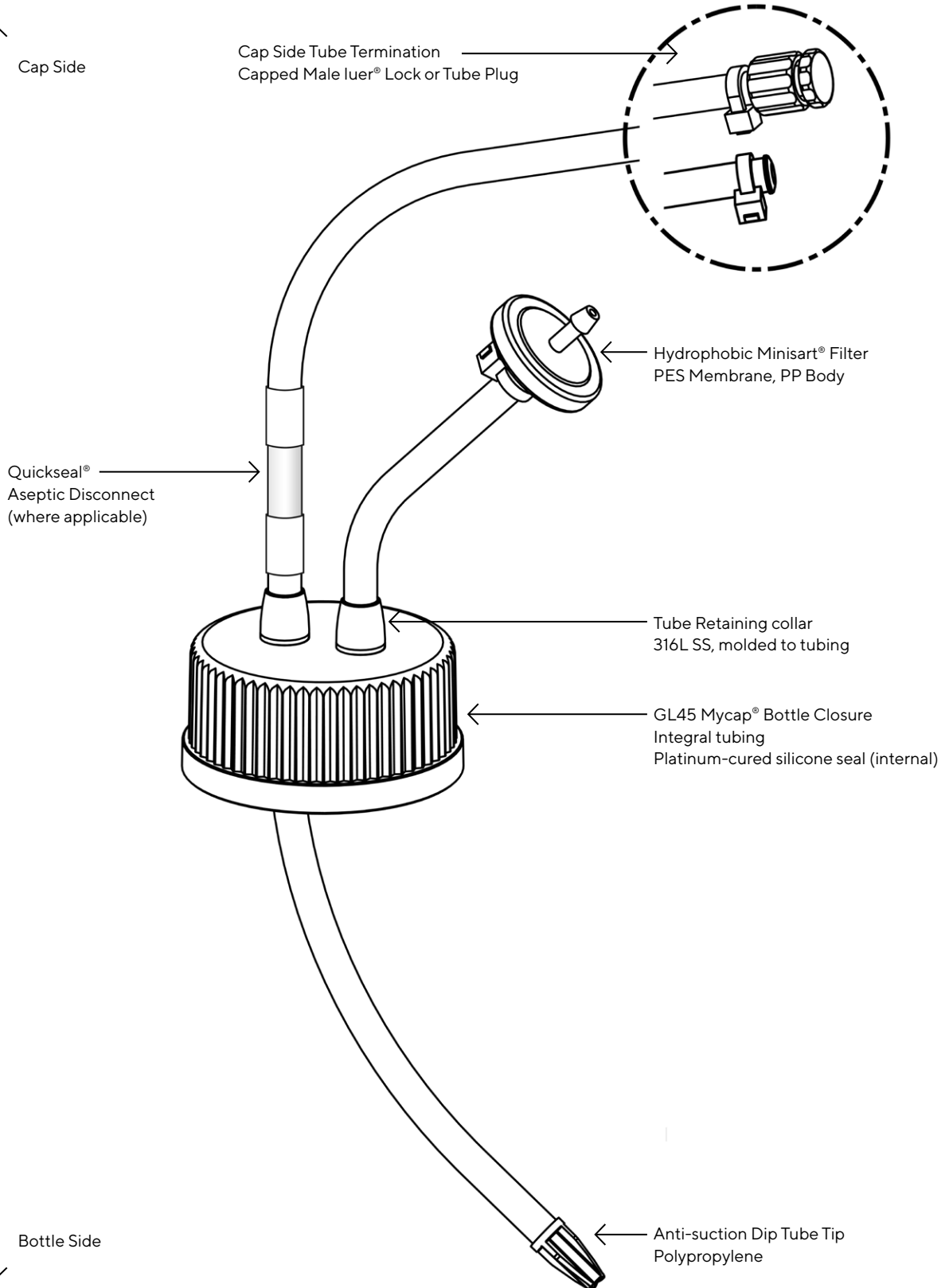
Luer<sup>®</sup> is a registered trademark of Becton Dickinson

Corning<sup>®</sup> and Pyrex<sup>®</sup> are registered trademarks of Corning Incorporated

LpH<sup>®</sup>, Vesphene<sup>®</sup> Spor Klenz<sup>®</sup> are registered trademarks of STERIS Corporation

# Image of Anatomy

↑  
Cap Side



# Ordering Information

## Series A GL45 Mycap® Bottle Closures

### Universal Design Features:

- Optimized tube lengths for Corning® Pyrex® round media storage bottles (PN: 1395-xxx)
- Bottle not included
- Platinum-cured silicone tubing
- Non-sterile
- Autoclavable

Part Number	Description	Bottle Size	Cap Size	Port 1	Port 2	Qty/ Pack
MYCAP13950100A	2 port GL45 Mycap® bottle closure	100 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 8.25"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13950250A	2 port GL45 Mycap® bottle closure	250 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 6.75"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13950500A	2 port GL45 Mycap® bottle closure	500 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 11"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13951000A	2 port GL45 Mycap® bottle closure	1,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 9.25"; capped male, luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13952000A	2 port GL45 Mycap® bottle closure	2,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 14"; capped male luer® fitting <b>Bottle Side:</b> 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	5

## Series B GL45 Mycap® Bottle Closures

### Universal Design Features:

- Optimized tube lengths for Corning® Pyrex® round media storage bottles (PN: 1395-xxx)
- Bottle not included
- C-Flex® tubing (formulation 374)
- Non-sterile
- Autoclavable (tubing may deform or melt during autoclave – exercise caution)

Part Number	Description	Bottle Size	Cap Size	Port 1	Port 2	Qty/ Pack
MYCAP13950100B	2 port GL45 Mycap® bottle closure	100 mL	GL45	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 14.5"; tube plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	5
MYCAP13950250B	2 port GL45 Mycap® bottle closure	250 mL	GL45	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 19"; tube plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	5
MYCAP13950500B	2 port GL45 Mycap® bottle closure	500 mL	GL45	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 17.25"; tube plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	5
MYCAP13951000B	2 port GL45 Mycap® bottle closure	1,000 mL	GL45	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 21.5"; tube plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	5
MYCAP13952000B	2 port GL45 Mycap® bottle closure	2,000 mL	GL45	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 20.25"; tube plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	5

## Series C GL45 Mycap® Bottle Closures

### Universal Design Features:

- Optimized tube lengths for Corning® Pyrex® round media storage bottles (PN: 1395-xxx)
- Bottle not included
- Platinum-cured silicone tubing
- Quickseal® aseptic disconnect
- Non-sterile
- Autoclavable

Part Number	Description	Bottle Size	Cap Size	Port 1	Port 2	Qty/ Pack
MYCAP13950100C	2 port GL45 Mycap® bottle closure	100 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 8.25"; Quickseal® at 2.25"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13950250C	2 port GL45 Mycap® bottle closure	250 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 6.75"; Quickseal® at 2"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13950500C	2 port GL45 Mycap® bottle closure	500 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 11"; Quickseal® at 2"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13951000C	2 port GL45 Mycap® bottle closure	1,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 9.25"; Quickseal® at 2.25"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13952000C	2 port GL45 Mycap® bottle closure	2,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 14"; Quickseal® at 2"; capped male luer® fitting <b>Bottle Side:</b> 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	5



## Series D GL45 Mycap® Bottle Closures

### Universal Design Features:

- Optimized tube lengths for Corning® Pyrex® round media storage bottles (PN: 1395-xxx)
- Bottle not included
- C-Flex® tubing (formulation 374)
- Quickseal® aseptic disconnect
- Non-sterile
- Autoclavable (tubing may deform or melt during autoclave – exercise caution)

Part Number	Description	Bottle Size	Cap Size	Port 1	Port 2	Qty/ Pack
MYCAP13950100D	2 port GL45 Mycap® bottle closure	100 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 14.5"; Quickseal® at 2.25"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13950250D	2 port GL45 Mycap® bottle closure	250 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 19"; Quickseal® at 2"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13950500D	2 port GL45 Mycap® bottle closure	500 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 17.25"; Quickseal® at 2.25"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13951000D	2 port GL45 Mycap® bottle closure	1,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 21.5"; Quickseal® at 2.25"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13952000D	2 port GL45 Mycap® bottle closure	2,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 20.25"; Quickseal® at 2"; tube plug at end <b>Bottle Side:</b> 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	5

## Series E GL45 Mycap® Bottle Closures

### Universal Design Features:

- Optimized tube lengths for Corning® Pyrex® round media storage bottles (PN: 1395-xxx)
- Bottle not included
- Tuflux® TPE
- Non-sterile
- Autoclavable (tubing may deform or melt during autoclave – exercise caution)

Part Number	Description	Bottle Size	Cap Size	Port 1	Port 2	Qty/ Pack
MYCAP13950100E	2 port GL45 Mycap® bottle closure	100 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 8.25"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13950250E	2 port GL45 Mycap® bottle closure	250 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 6.75"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13950500E	2 port GL45 Mycap® bottle closure	500 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 11"; capped male luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13951000E	2 port GL45 Mycap® bottle closure	1,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 9.25"; capped male, luer® fitting <b>Bottle Side:</b> 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	5
MYCAP13952000E	2 port GL45 Mycap® bottle closure	2,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 14"; capped male luer® fitting <b>Bottle Side:</b> 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	5

## Series F GL45 Mycap® Bottle Closures

### Universal Design Features:

- Optimized tube lengths for Corning® Pyrex® round media storage bottles (PN: 1395-xxx)
- Bottle not included
- Tuflux® TPE tubing
- Quickseal® aseptic disconnect
- Non-sterile
- Autoclavable (tubing may deform or melt during autoclave – exercise caution)

Part Number	Description	Bottle Size	Cap Size	Port 1	Port 2	Qty/ Pack
MYCAP13950100F	2 port GL45 Mycap® bottle closure	100 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 14.5"; Quickseal® at 2.25"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13950250F	2 port GL45 Mycap® bottle closure	250 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 19"; Quickseal® at 2"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13950500F	2 port GL45 Mycap® bottle closure	500 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 17.25"; Quickseal® at 2.25"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13951000F	2 port GL45 Mycap® bottle closure	1,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 21.5"; Quickseal® at 2.25"; tube plug at end <b>Bottle Side:</b> 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	5
MYCAP13952000F	2 port GL45 Mycap® bottle closure	2,000 mL	GL45	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 20.25"; Quickseal® at 2"; tube plug at end <b>Bottle Side:</b> 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	5

# Mycap<sup>®</sup> Bottle Assemblies

## Applications

Mycap<sup>®</sup> is ideal for critical fluid-handling applications.

Mycap<sup>®</sup> 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

## Technical Data

### Specifications

Elastomer Seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex <sup>®</sup> formulation 374
Cap Material	Polycarbonate (not fluid contact)
Bottle Style	Nalgene <sup>®</sup> PETG bottle
Tube Ports	Inlet   Outlet, Air vent - 25 mm PES membrane Minisart <sup>®</sup>
Fittings	Luer <sup>®</sup> Lock or tube plug
Sterilization	Gamma Irradiated

C-Flex<sup>®</sup> is a registered trademark of Saint-Gobain Performance Plastics Corporation

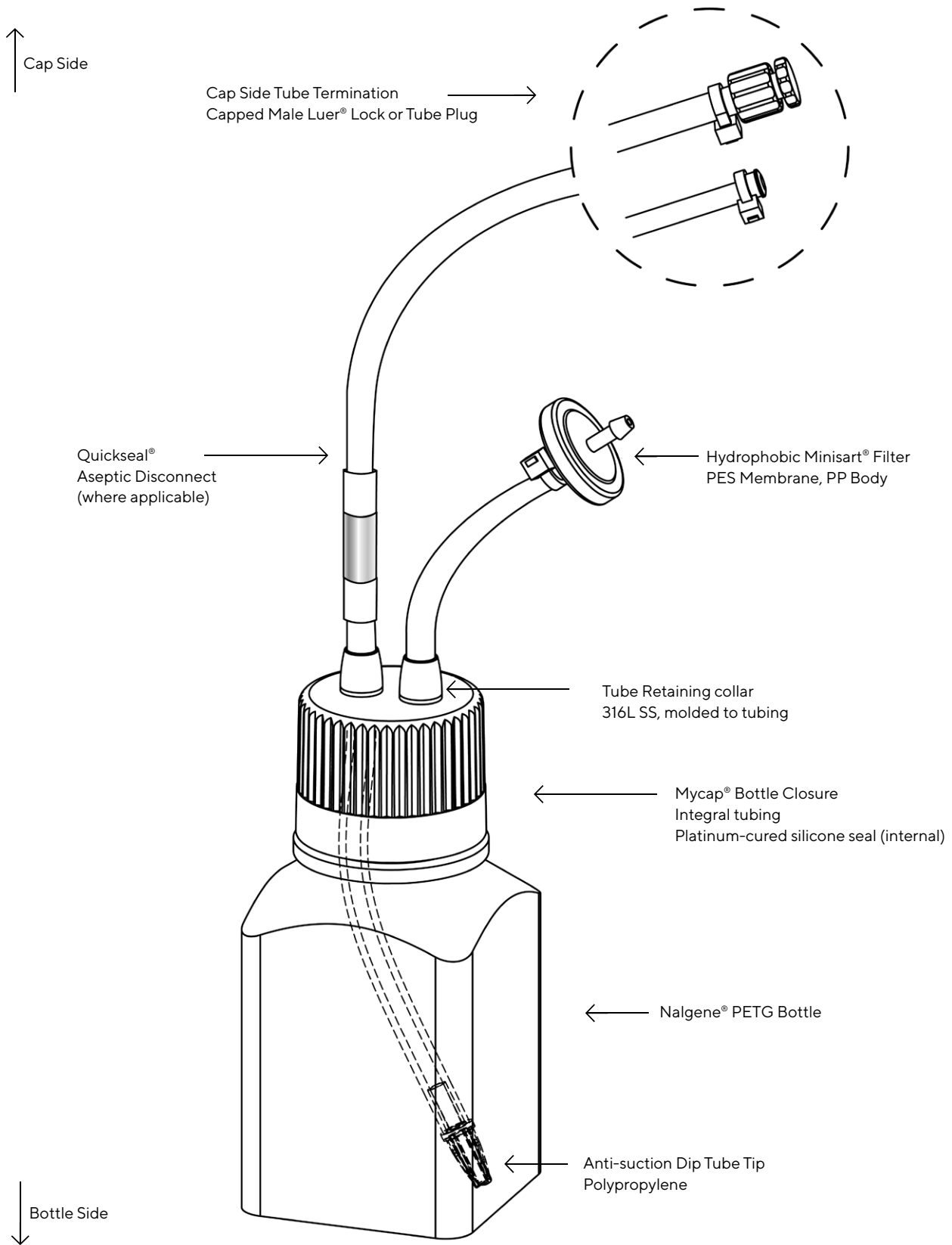
Luer<sup>®</sup> is a registered trademark of Becton Dickinson

Dow Corning<sup>®</sup> is a registered trademark of Dow Corning Corporation

LpH<sup>®</sup>, Vesphene<sup>®</sup> Spor Klensz<sup>®</sup> are registered trademarks of STERIS Corporation

Nalgene<sup>®</sup> is a registered trademark of ThermoFisher Scientific

# Image of Anatomy



# Ordering Information

## Series A Mycap® Bottle Assemblies

### Universal Design Features:

- Optimized Tube Lengths including dip tube for fluid transfer
- Includes Nalgene® PETG Media Bottle
- Platinum-Cured Silicone Tubing (Dow-Corning® PHARMA-50)
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPNGPG0030A	2 port Mycap® bottle closure with 30 mL bottle	30 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 3.75"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0060A	2 port Mycap® bottle closure with 60 mL bottle	60 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 3.25"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0125A	2 port Mycap® bottle closure with 125 mL bottle	125 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 8"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0250A	2 port Mycap® bottle closure with 250 mL bottle	250 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 6.5"; capped male, Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0500A	2 port Mycap® bottle closure with 500 mL bottle	500 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 11.25; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG1000A	2 port Mycap® bottle closure with 1,000 mL bottle	1,000 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 9.5"; capped male, Luer® fitting <b>Bottle Side:</b> 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

## Series B Mycap® Bottle Closures

### Universal Design Features:

- Optimized Tube Lengths including dip tube for fluid transfer
- Includes Nalgene® PETG Media Bottle
- C-Flex® tubing (formulation 374)
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPNGPG0030B	2 port Mycap® bottle closure with 30 mL bottle	30 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 15.75"; capped male Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0060B	2 port Mycap® bottle closure with 60 mL bottle	60 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 15.25"; capped male Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0125B	2 port Mycap® bottle closure with 125 mL bottle	125 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 14"; capped male Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0250B	2 port Mycap® bottle closure with 250 mL bottle	250 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 18.5"; capped male, Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0500B	2 port Mycap® bottle closure with 500 mL bottle	500 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 17.25; capped male Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG1000B	2 port Mycap® bottle closure with 1,000 mL bottle	1,000 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 21.75"; capped male, Luer® fitting <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

## Series C Mycap® Bottle Assemblies

### Universal Design Features:

- Optimized Tube Lengths for including dip tube for fluid transfer
- Includes Nalgene® PETG Media Bottle
- Platinum-Cured Silicone Tubing (Dow-Corning® PHARMA-50)
- Quickseal® aseptic disconnect
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPNGPG0030C	2 port Mycap® bottle closure with 30 mL bottle	30 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 3.75"; Quickseal® at 1.75"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0060C	2 port Mycap® bottle closure with 60 mL bottle	60 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 3.25"; Quickseal® at 2.25"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0125C	2 port Mycap® bottle closure with 125 mL bottle	125 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 8"; Quickseal® at 2"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0250C	2 port Mycap® bottle closure with 250 mL bottle	250 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 6.5"; Quickseal® at 3"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG0500C	2 port Mycap® bottle closure with 500 mL bottle	500 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 11.25"; Quickseal® at 2"; capped male Luer® fitting <b>Bottle Side:</b> 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
MYCAPNGPG1000C	2 port Mycap® bottle closure with 1,000 mL bottle	1,000 mL	<b>Tube Size:</b> 1/8" ID × 1/4" OD <b>Cap Side:</b> 9.5"; Quickseal® at 2"; capped male, Luer® fitting <b>Bottle Side:</b> 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



## Series D Mycap® Bottle Assemblies

### Universal Design Features:

- Optimized Tube Lengths for including dip tube for fluid transfer
- Includes Nalgene® PETG Media Bottle
- C-Flex tubing (formulation 374)
- Quickseal® aseptic disconnect
- Gamma Irradiated
- Individually Double Bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPNGPG0030D	2 port Mycap® bottle closure with 30 mL bottle	30 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 15.75"; Quickseal® at 1.75"; plug at end <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0060D	2 port Mycap® bottle closure with 60 mL bottle	60 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 15.25"; Quickseal® at 2.25"; plug at end <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0125D	2 port Mycap® bottle closure with 125 mL bottle	125 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 14"; Quickseal® at 2"; plug at end <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0250D	2 port Mycap® bottle closure with 250 mL bottle	250 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 18.5"; Quickseal® at 2.5"; plug at end <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG0500D	2 port Mycap® bottle closure with 500 mL bottle	500 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 17.25"; Quickseal® at 2"; plug at end <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPNGPG1000D	2 port Mycap® bottle closure with 1,000 mL bottle	1,000 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 21.75"; Quickseal® at 1.5"; plug at end <b>Bottle Side:</b> To Bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

# Mycap<sup>®</sup> Centrifuge Tubes

## Applications

Mycap<sup>®</sup> is ideal for critical fluid-handling applications.

Mycap<sup>®</sup> 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

## Technical Data

### Specifications

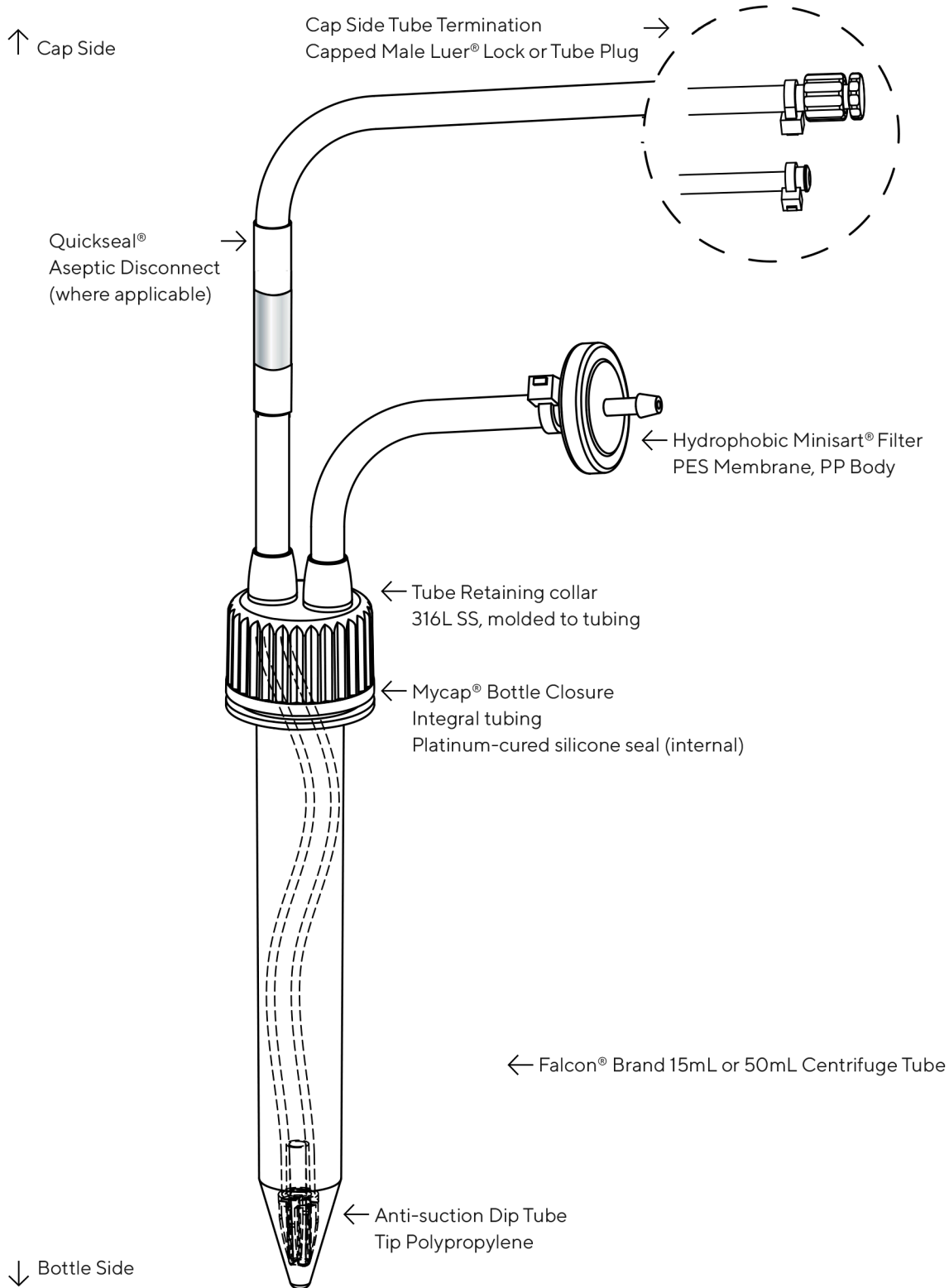
Elastomer seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex <sup>®</sup> formulation 374 or Tuflux <sup>®</sup> TPE
Cap material	Polycarbonate (not fluid contact)
Bottle style	polystyrene and polypropylene centrifuge tube
Tube ports	Inlet   outlet, Air vent - 25 mm PES membrane Minisart <sup>®</sup>
Fittings	luer <sup>®</sup> lock or tube plug
Sterilization	Gamma irradiated

C-Flex<sup>®</sup> is a registered trademark of Saint-Gobain Performance Plastics Corporation

luer<sup>®</sup> is a registered trademark of Becton Dickinson

LpH<sup>®</sup>, Vesphene<sup>®</sup> Spor Klenz<sup>®</sup> are registered trademarks of STERIS Corporation

# Image of Anatomy



# Ordering Information

## Series A Mycap® Bottle Assemblies

Universal design features:

- Optimized tube lengths including dip tube for fluid transfer
- Includes 15 mL polystyrene or 50 mL polypropylene centrifuge tube
- Platinum-cured silicone tubing
- Gamma irradiated
- Individually double bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPCPSPG0015A	2 port Mycap® bottle closure with 15 mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 7.25"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPCPPG0050A	2 port Mycap® bottle closure with 50 mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 7.5"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

## Series B Mycap® Bottle Closures

Universal design features:

- Optimized tube lengths including dip tube for fluid transfer
- Includes 15 mL polystyrene or 50 mL polypropylene centrifuge tube
- C-Flex® tubing (formulation 374)
- Gamma irradiated
- Individually double bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPCPSPG0015B	2 port Mycap® bottle closure with 15 mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.25"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPCPPG0050B	2 port Mycap® bottle closure with 50 mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.5"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

## Series C Mycap® Bottle Assemblies

Universal design features:

- Optimized tube lengths for including dip tube for fluid transfer
- Includes 15 mL polystyrene or 50 mL polypropylene centrifuge tube
- Platinum-cured silicone tubing
- Quickseal® aseptic disconnect
- Gamma irradiated
- Individually double bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015C	2 port Mycap® bottle closure with 15 mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 7.25"; Quickseal® at 2.25"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 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 50 mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 7.5"; Quickseal® at 2.5"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

## Series D Mycap® Bottle Assemblies

Universal design features:

- Optimized tube lengths for including dip tube for fluid transfer
- Includes 15 mL polystyrene or 50 mL polypropylene centrifuge tube
- C-Flex® tubing (formulation 374)
- Quickseal® aseptic disconnect
- Gamma irradiated
- Individually double bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015D	2 port Mycap® bottle closure with 15 mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.25"; Quickseal® at 2.25"; plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPCPPG0050D	2 port Mycap® bottle closure with 50 mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.5"; Quickseal® at 2.5"; plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

## Series E Mycap® Bottle Closures

Universal design features:

- Optimized tube lengths including dip tube for fluid transfer
- Includes 15 mL polystyrene or 50 mL polypropylene centrifuge tube
- Tuflux® TPE tubing (formulation 374)
- Gamma irradiated
- Individually double bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015E	2 port Mycap® bottle closure with 15 mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.25"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPCPPG0050E	2 port Mycap® bottle closure with 50 mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.5"; capped male luer® fitting <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

## Series F Mycap® Bottle Assemblies

Universal design features:

- Optimized tube lengths for including dip tube for fluid transfer
- Includes 15 mL polystyrene or 50 mL polypropylene centrifuge tube
- Tuflux® TPE tubing (formulation 374)
- Quickseal® aseptic disconnect
- Gamma irradiated
- Individually double bagged

Part Number	Description	Bottle Size	Port 1	Port 2	Qty/Pack
MYCAPCPSG0015F	2 port Mycap® bottle closure with 15 mL centrifuge tube	15 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.25"; Quickseal® at 2.25"; plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1
MYCAPCPPG0050F	2 port Mycap® bottle closure with 50 mL centrifuge tube	50 mL	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 13.5"; Quickseal® at 2.5"; plug at end <b>Bottle Side:</b> To bottom; dip tube tip	<b>Tube Size:</b> 1/8" ID x 1/4" OD <b>Cap Side:</b> 3"; Minisart® hydrophobic PES filter <b>Bottle Side:</b> None	1

# Mycap<sup>®</sup> Single-Bottle Assemblies for Media

## Applications

Mycap<sup>®</sup> for Media make media distribution and storage quicker and easier than ever before.

Bottle assemblies include a Sartopore<sup>®</sup> 2 filter for sterile filtration directly into easy-to-handle bottles. Simply connect the filter inlet to your bulk media source to sterilize media as the bottle fills.

Cut the Quickseal<sup>®</sup> aseptic disconnect after filling and your aseptic media distribution is complete, without ever having to go into a biosafety cabinet.

Choose tube welding or aseptic connector for convenient aseptic connection and dispense sterile media at point-of-use, including bioreactors.

## Technical Data

### Specifications

Elastomer Seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex <sup>®</sup> formulation 374
Cap Material	Polycarbonate (not fluid contact)
Bottle Style	PETG Square bottle
Tube Ports	Inlet: Sartopore <sup>®</sup> 2 Media Filter, Air vent - 25 mm PES membrane Minisart <sup>®</sup>
Fittings	Outlet: Aseptiquik <sup>®</sup> S or tube plug
Sterilization	Gamma Irradiated

C-Flex<sup>®</sup> is a registered trademark of Saint-Gobain Performance Plastics Corporation

Luer<sup>®</sup> is a registered trademark of Becton Dickinson

LpH<sup>®</sup>, Vesphene<sup>®</sup> Spor Klenz<sup>®</sup> are registered trademarks of STERIS Corporation

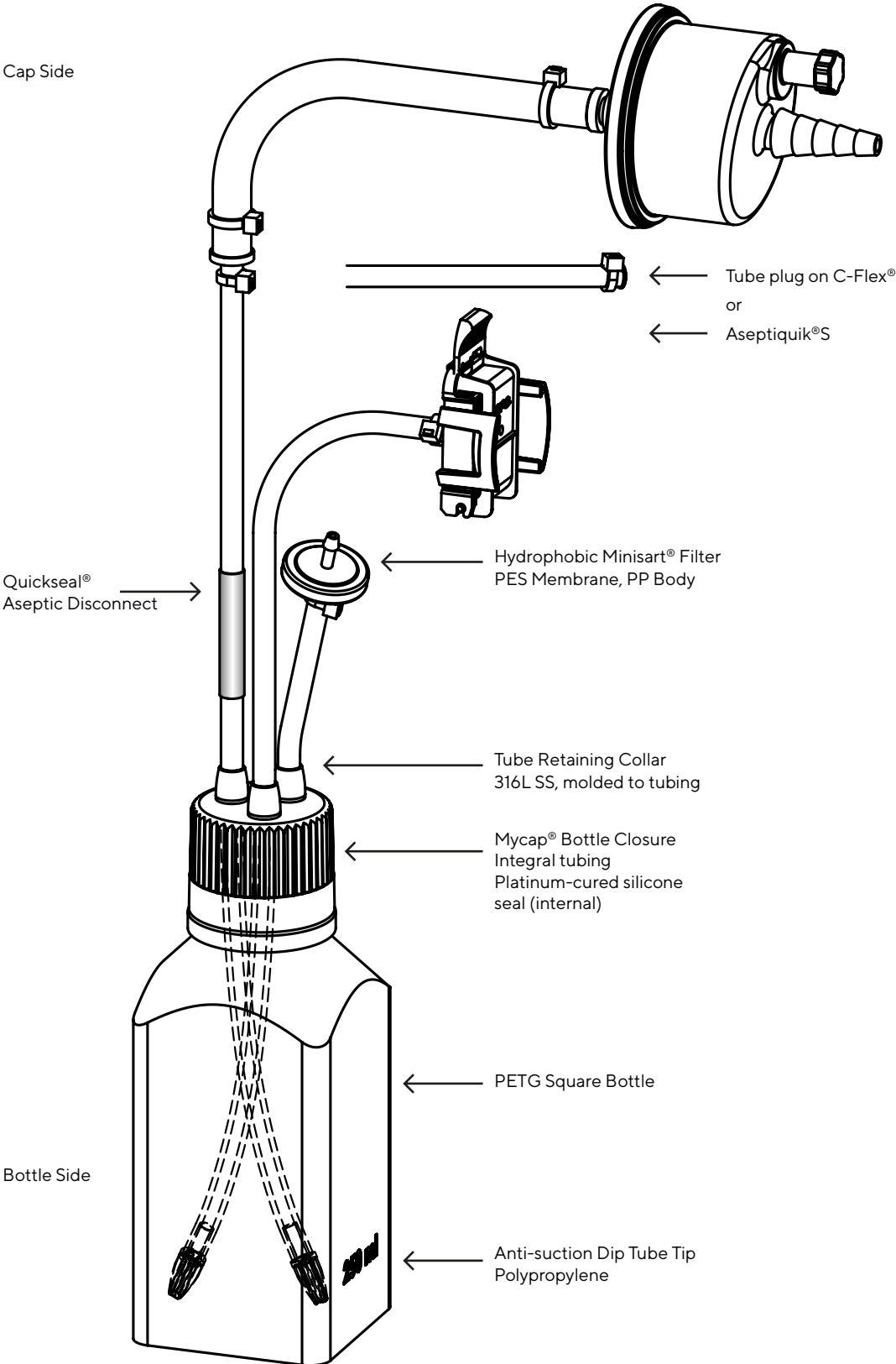
Sartopore<sup>®</sup>, Minisart<sup>®</sup>, Tuflux<sup>®</sup> TPE, Quickseal<sup>®</sup> and Mycap<sup>®</sup> is a registered trademark of Sartorius-Stedim Biotech

AseptiQuik<sup>®</sup>, is a registered trademark of Colder Products Company

# Image of Anatomy

## Single Bottle with Aseptiquik® S

↑  
Cap Side



← Tube plug on C-Flex®  
or  
← Aseptiquik® S

Quickseal®  
Aseptic Disconnect →

← Hydrophobic Minisart® Filter  
PES Membrane, PP Body

← Tube Retaining Collar  
316L SS, molded to tubing

← Mycap® Bottle Closure  
Integral tubing  
Platinum-cured silicone  
seal (internal)

← PETG Square Bottle

↓  
Bottle Side

← Anti-suction Dip Tube Tip  
Polypropylene



# Ordering Information

## Single Bottle; Aseptic Connection by Tube Welding

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000010204T	1,000 mL Square Media Bottle; 1/8 x 1/4 Tuflux TPE® Tubing	1,000 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500010204T	500 mL Square Media Bottle; 1/8 x 1/4 Tuflux TPE®	500 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250010204T	250 mL Square Media Bottle; 1/8 x 1/4 Tuflux TPE® Tubing	250 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1

## Single Bottle; Aseptic Connection by Tube Welding

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000010204W	1,000 mL Square Media Bottle; 1/8" x 1/4" C-Flex® Tubing	1,000 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4" - 1/2" barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500010204W	500 mL Square Media Bottle; 1/8" x 1/4" C-Flex® Tubing	500 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4" - 1/2" barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250010204W	250 mL Square Media Bottle; 1/8" x 1/4" C-Flex® Tubing	250 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4" - 1/2" barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1

## Single Bottle; Aseptic Connection by Aseptic Connector

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000010204A	1,000 mL Square Media Bottle; ⅝ × ¼ Silicone Tubing; Aseptiquik® S	1,000 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; ¼–½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500010204A	500 mL Square Media Bottle; ⅝ × ¼ Silicone Tubing; Aseptiquik® S	500 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; ¼–½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250010204A	250 mL Square Media Bottle; ⅝ × ¼ Silicone Tubing; Aseptiquik® S	250 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; ¼–½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1

## Single Bottle; Aseptic Connection by Tube Welding or Aseptic Connector

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000010204D	1,000 mL Square Media Bottle; 1/8 x 1/4 Tuflux TPE® Tubing w/Aseptiquik® S	1,000 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500010204D	500 mL Square Media Bottle; 1/8 x 1/4 Tuflux TPE® Tubing w/Aseptiquik® S	500 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250010204D	250 mL Square Media Bottle; 1/8 x 1/4 Tuflux TPE® Tubing w/Aseptiquik® S	250 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD1000010204B	1,000 mL Square Media Bottle; 1/8 x 1/4 C-Flex® Tubing w/Aseptiquik® S	1,000 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500010204B	500 mL Square Media Bottle; 1/8 x 1/4 C-Flex® Tubing w/Aseptiquik® S	500 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250010204B	250 mL Square Media Bottle; 1/8 x 1/4 C-Flex® Tubing w/Aseptiquik® S	250 mL	Size 4 Sartopore® 2 capsule filter; 0.2 µm; 1/4 - 1/2 barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1

# Mycap<sup>®</sup> Manifold Bottle Assemblies for Media

## Applications

Mycap<sup>®</sup> for Media make media distribution and storage quicker and easier than ever before. Mycap<sup>®</sup> for Media includes six bottles.

Manifolds include a Sartopore<sup>®</sup> 2 filter for sterile filtration directly into easy-to-handle bottles. Simply connect the filter inlet to your bulk media source to sterilize media as the bottle fills.

Cut the Quickseal<sup>®</sup> aseptic disconnect after filling and your aseptic media distribution is complete, without ever having to go into a biosafety cabinet.

Choose tube welding or aseptic connector for convenient aseptic connection and dispense sterile media at point-of-use, including bioreactors.

## Technical Data

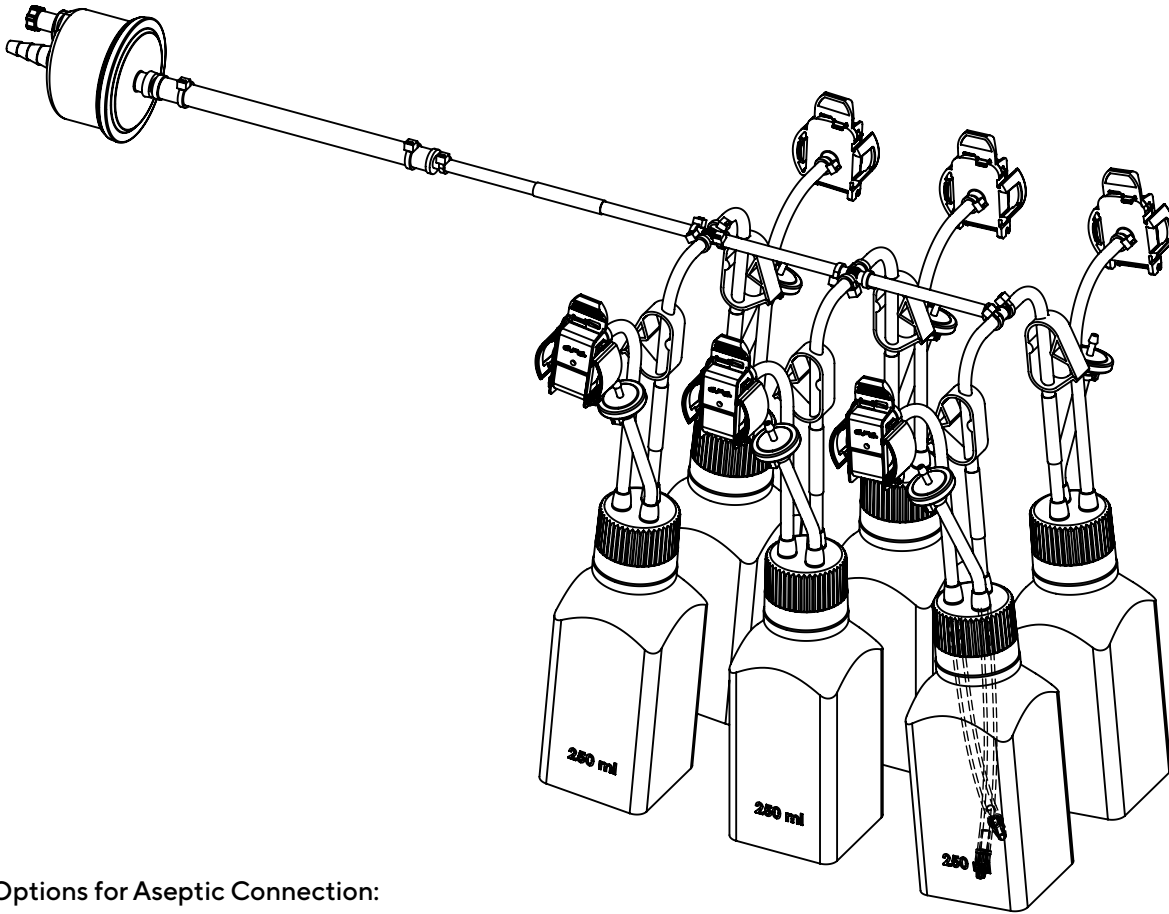
### Specifications

Elastomer Seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex <sup>®</sup> formulation 374
Cap Material	Polycarbonate (not fluid contact)
Bottle Style	PETG Square bottle
Tube Ports	Inlet: Sartopore <sup>®</sup> 2 Media Filter, Air vent - 25 mm PES membrane Minisart <sup>®</sup>
Fittings	Outlet: Aseptiquik <sup>®</sup> S or tube plug
Sterilization	Gamma Irradiated

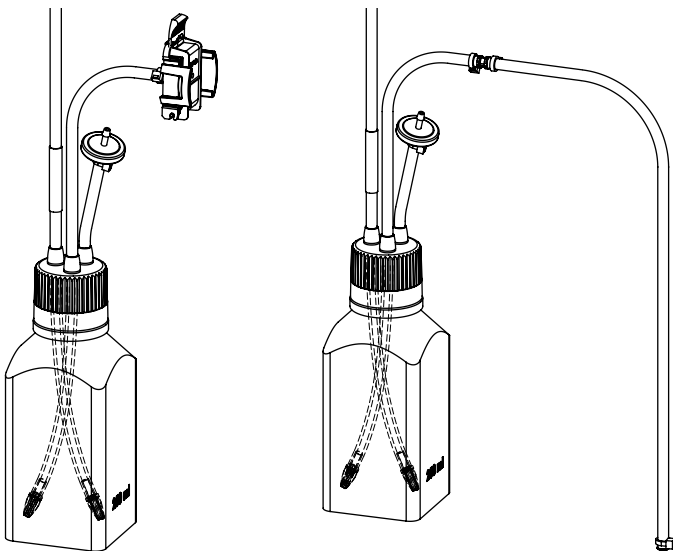
C-Flex<sup>®</sup> is a registered trademark of Saint-Gobain Performance Plastics Corporation  
LpH<sup>®</sup>, Vesphene<sup>®</sup> Spor Klens<sup>®</sup> are registered trademarks of STERIS Corporation  
Sartopore<sup>®</sup>, Minisart<sup>®</sup>, Quickseal<sup>®</sup> and Mycap<sup>®</sup> is a registered trademark of Sartorius-Stedim Biotech  
Aseptiquik<sup>®</sup>, is a registered trademark of Colder Products Company

# Image of Anatomy

## Six-Bottle Manifold With Aseptiquik® S



## Two Options for Aseptic Connection:



# Ordering Information

## Six-Bottle Manifold; Aseptic Connection by Tube Welding or Aseptic Connectors

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Manifold Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000060204B	6 × 1,000 mL Square Media Bottle; ½ × ¼ C-Flex® Tubing w/Aseptiquik® S	1,000 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼ - ½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500060204B	6 × 500 mL Square Media Bottle; ½ × ¼ C-Flex® Tubing w/Aseptiquik® S	500 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼ - ½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250060204B	6 × 250 mL Square Media Bottle; ½ × ¼ C-Flex® Tubing w/Aseptiquik® S	250 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼ - ½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1

## Six-Bottle Manifold; Aseptic Connection With Aseptic Connector

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Manifold Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000060204A	6 × 1,000 mL Square Media Bottle; ½ × ¼ Silicone Tubing; Aseptiquik® S	1,000 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼–½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500060204A	6 × 500 mL Square Media Bottle; ½ × ¼ Silicone Tubing; Aseptiquik® S	500 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼–½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250060204A	6 × 250 mL Square Media Bottle; ½ × ¼ Silicone Tubing; Aseptiquik® S	250 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼–½ barb inlet; Quickseal®; dip tube to bottom	Aseptiquik® S; dip tube to bottom	Minisart® hydrophobic PES vent filter	1



## Six-Bottle Manifold; Aseptic Connection by Tube Welding

### Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 µm Filter at Manifold Inlet for Sterile Media Fill
- Includes PETG Square Media Bottle
- Gamma Irradiated
- Individually Double Bagged


Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/ Pack
MCMD1000060204W	6 × 1,000 mL Square Media Bottle; ½ × ¼ C-Flex® Tubing	1,000 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼ – ½ barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500060204W	6 × 500 mL Square Media Bottle; ½ × ¼ C-Flex® Tubing	500 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼ – ½ barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250060204W	6 × 250 mL Square Media Bottle; ½ × ¼ C-Flex® Tubing	250 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¼ – ½ barb inlet; Quickseal®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1

**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](http://www.sartorius.com)