



4: MYCAP®; 38-430 Cap; 3-Ports;  
 P1: 1/8"x1/4" Dow Pharma 50 Sil, CS-6.5", BS-140mm, QS-3";  
 P2: 1/8"x1/4" Dow Pharma 50 Sil, CS-6.5", BS-140mm, QS-0";  
 P3: 1/8"x1/4" Dow Pharma 50 Sil, CS-3", BS-0", QS-0"

2: 1/4"IDx7/16"OD Dow Pharma 50  
 Silicone tube; 6"OAL

3: 1/4"x1/8"HB Reducer

7: Vent Filter;  
 25mm MiniSart;  
 1/8"HB I/O;  
 PES(GI/Autoclave)  
 (Hydrophobic)

11: Cable tie, blue;  
 TYP

6: 250ml PETG Bottle

5: 1/8" Dip Tube Tip w/Notch

1: Sartopore® 2 capsule filter;  
 0.2µm; 150cm²;  
 1/4"-1/2" HB I/O; (GI)

8: 1/8"HB Straight Connector

9: 1/8"IDx1/4"OD C-FLEX® 374 tube;  
 18"OAL

10: Colder AseptiQuik® S;  
 1/8"HB; Not Autoclaveable

**Bill of Materials:**

Item	Qty	AP Number	Material
1	1	AP-2365	-PP, PES, Polyester -Silicone
2	1	AP-2119	-ADCF Silicone
3	1	AP-0029	-ADCF Polypropylene
4	1	AP-3313	-ADCF Silicone
5	2	AP-2170	-ADCF Polypropylene
6	1	AP-0297	-PETG
7	1	AP-2767	-Polypropylene, PES
8	1	AP-0089	-ADCF Polypropylene
9	1	AP-0378	-ADCF C-FLEX®(374)
10	1	AP-0917	-PC, PE, Silicone
11	7	AP-0133	-Tefzel®

**Revision History:**

- 0 - Initial release
- 1 - Adjusted Sartopore® filter (AP-2365) for improved image scale
- 2 - Updated MYCAP® P1 QS 2" to 3"

**Dimensional Tolerances:**

Tube Lengths:  
 0-12" ±0.25"  
 0-36" ±0.5"  
 36"+ ±1.0"

**Notes:**

1. No assembly required.
2. Certificate of Conformance & Certificate of Irradiation provided for each lot.
3. Individually double bagged.
4. Fluid contact materials meet criteria for USP Class VI Testing. Consult Validation Guide for additional tests and details.
5. Gamma Irradiated.
6. Drawing not to scale.

**PROPRIETARY AND CONFIDENTIAL**

The information contained herein is the sole property of Sartorius Stedim North America. Reproduction or distribution is prohibited without written consent from Sartorius Stedim North America.

**sartorius stedim**  
 biotech  
 Sartorius Stedim North America, Inc  
 80 Progress Ave  
 New Oxford, PA 17350  
 www.sartorius.com

Customer Approval: Approval Not Required	Revision: 2	Released By:	Release Date:
Approval Date:			
Part Number: MCMD0250010204B			
Drawing Number: MCMD0250010204B			

Title:  
 250ml Media; 1/8"x1/4" C-FLEX® w/ASQ