



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

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

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"

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: 1/8"HB Tube Plug

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-1181	-ADCF Polypropylene
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 North America, Inc
 80 Progress Ave
 New Oxford, PA 17350

biotech
 www.sartorius.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MCMD0250010204W

Drawing Number: MCMD0250010204W

Revision: 2

Released By:

Release Date:

Title:

250ml Media; 1/8"x1/4" C-FLEX®