



1: Sartopore® 2 capsule filter; 0.2µm; 300cm²; 1/4"-1/2" HB I/O (GI)

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

9: Ratchet clamp; 1/4"OD tubing

10: 1/8"HB Straight Connector

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

6: MYCAP™; 38-430 Cap; 3-Ports; P1: 1/8"x1/4" Dow Pharma 50 Sil, CS-9.5", BS-215mm, QS-1.5"; P2: 1/8"x1/4" Dow Pharma 50 Sil, CS-9.5", BS-215mm, QS-0"; P3: 1/8"x1/4" Dow Pharma 50 Sil, CS-3", BS-0", QS-0"

11: 1/8"IDx1/4"OD C-FLEX® 374 tube; 18"OAL

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

8: 1000ml PETG Bottle

7: 1/8" Dip Tube Tip w/Notch

14: 1/8"IDx1/4"OD Dow Pharma50 Silicone tube; 3"OAL

15: 1/8"HB Tee

4: 1/8"IDx1/4"OD Dow Pharma 50 Silicone tube; 6"OAL w/Qs collar at 3"

5: 1/8"HB Cross

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

16: Cable tie, blue; TYP

Bill of Materials:

| Item | Qty | AP Number | Material |
|------|-----|-----------|-------------------------------|
| 1 | 1 | AP-3311 | -PP, PES, Polyester -Silicone |
| 2 | 1 | AP-2119 | -ADCF Silicone |
| 3 | 1 | AP-0029 | -ADCF Polypropylene |
| 4 | 1 | AP-0827 | -ADCF Silicone |
| 5 | 2 | AP-0638 | -ADCF Polypropylene |
| 6 | 6 | AP-3110 | -ADCF Silicone |
| 7 | 12 | AP-2170 | -ADCF Polypropylene |
| 8 | 6 | AP-1330 | -PETG |
| 9 | 6 | AP-0118 | -Polypropylene |
| 10 | 6 | AP-0089 | -ADCF Polypropylene |
| 11 | 6 | AP-0378 | -ADCF C-FLEX®(374) |
| 12 | 6 | AP-0917 | -PC, PE, Silicone |
| 13 | 6 | AP-2767 | -Polypropylene, PES |
| 14 | 2 | AP-2291 | -ADCF Silicone |
| 15 | 1 | AP-1488 | -ADCF Polypropylene |
| 16 | 38 | AP-0133 | -Tefzel® |

Revision History:

- 0 - Initial release
- 1 - Adjusted Sartopore® filter (AP-3311) for improved image scale

Dimensional Tolerances:

- Tubing 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 AllPure Technologies, LLC. Reproduction or distribution is prohibited without written consent from AllPure Technologies, LLC.

Title:

6x1000ml Media; 1/8"x1/4" C-FLEX® w/ASQ



AllPure Technologies, LLC
80 Progress Ave
New Oxford, PA 17350

www.sartorius.com
www.allpureinc.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MCMD1000060204B

Drawing Number: MCMD1000060204B

Revision: 1

Released By:

Release Date: