



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

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

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

12: 1/8"IDx1/4"OD Dow Pharma50
Silicone tube; 3"OAL

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

5: 1/8"HB Cross

9: Ratchet clamp;
1/4"OD tubing

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

11: 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"

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

8: 1000ml PETG Bottle

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-0917	-PC, PE, Silicone
11	6	AP-2767	-Polypropylene, PES
12	2	AP-2291	-ADCF Silicone
13	1	AP-1488	-ADCF Polypropylene
14	26	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" ASQ

sartorius stedim
biotech
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: MCMD1000060204A
 Drawing Number: MCMD1000060204A

Revision: 1 Released By: Release Date: