



**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-3313	-ADCF Silicone
7	12	AP-2170	-ADCF Polypropylene
8	6	AP-0297	-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
- 2 - Updated MYCAP® P1 QS 2" to 3"

**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 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

www.sartorius.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MCMD0250060204A

Drawing Number: MCMD0250060204A

Revision: 2

Released By:

Release Date:

Title:

6x250ml Media; 1/8"x1/4" ASQ