



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-0917	-PC, PE, Silicone
9	5	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:

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

Customer Approval: Approval Not Required
Approval Date:
Part Number: MCMD0250010204A
Drawing Number: MCMD0250010204A

Title: 250ml Media; 1/8"x1/4" ASQ
Revision: 2
Released By:
Release Date: