



Bill of Materials:			
Item	Qty	AP Num.	Material
1	1	AP-1181	-Polypropylene
2	1	AP-4837	-C-Flex®(374), Silicone
3	1	AP-1184	-Polypropylene
4	1	AP-2170	-Polypropylene
5	1	AP-1393	-PETG
6	1	AP-2767	-Polypropylene, PES
7	2	AP-0133	-Tefzel®

Revision History:
 0 - Initial release
 1 - Update P1 CS 20.25" to 19.75"
 - Remove 'ADCF' from BOM Material list

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.

Title:
 Mycap®; 2000mL PETG; 2-Ports;
 1/8"x1/4" C-Flex® w/Dip Tube;
 QS on inlet port