



Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-1181	-ADCF Polypropylene
2	1	AP-3219	-ADCF C-FLEX®(374) -ADCF Silicone
3	1	AP-2170	-ADCF Polypropylene
4	1	AP-0202	-Polypropylene
5	1	AP-2767	-Polypropylene, PES
6	2	AP-0133	-Tefzel®

Revision History:

0 - Initial release

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.

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: MYCAP CPP G0050D
 Drawing Number: MYCAP CPP G0050D
 Revision: 0
 Released By:

Title:
 MYCAP™; 50ml PP Centrifuge; 2-Ports;
 1/8"x1/4" C-FLEX® w/Dip Tube Tip; QS on inlet
 Release Date: