



Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-1181	-ADCF Polypropylene
2	1	AP-3211	-ADCF C-FLEX®(374) -ADCF Silicone
3	1	AP-1048	-HDPE
4	1	AP-2170	-ADCF Polypropylene
5	1	AP-0639	-Polystyrene
6	1	AP-2767	-Polypropylene, PES
7	2	AP-0133	-Tefzel®

Revision History:

0 - Initial release
 1 - P1: CS 13.25" to 13.5"; BS 124mm to 117mm

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:

MYCAP™: 15ml PS Centrifuge; 2-Ports;
 1/8"x1/4" C-FLEX® w/Dip Tube Tip



AllPure Technologies, LLC
 80 Progress Ave
 New Oxford, PA 17350

biotech
 www.sartorius.com
 www.allpureinc.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MYCAP CPS G0015B

Drawing Number: MYCAP CPS G0015B

Revision: 1

Released By:

Release Date: