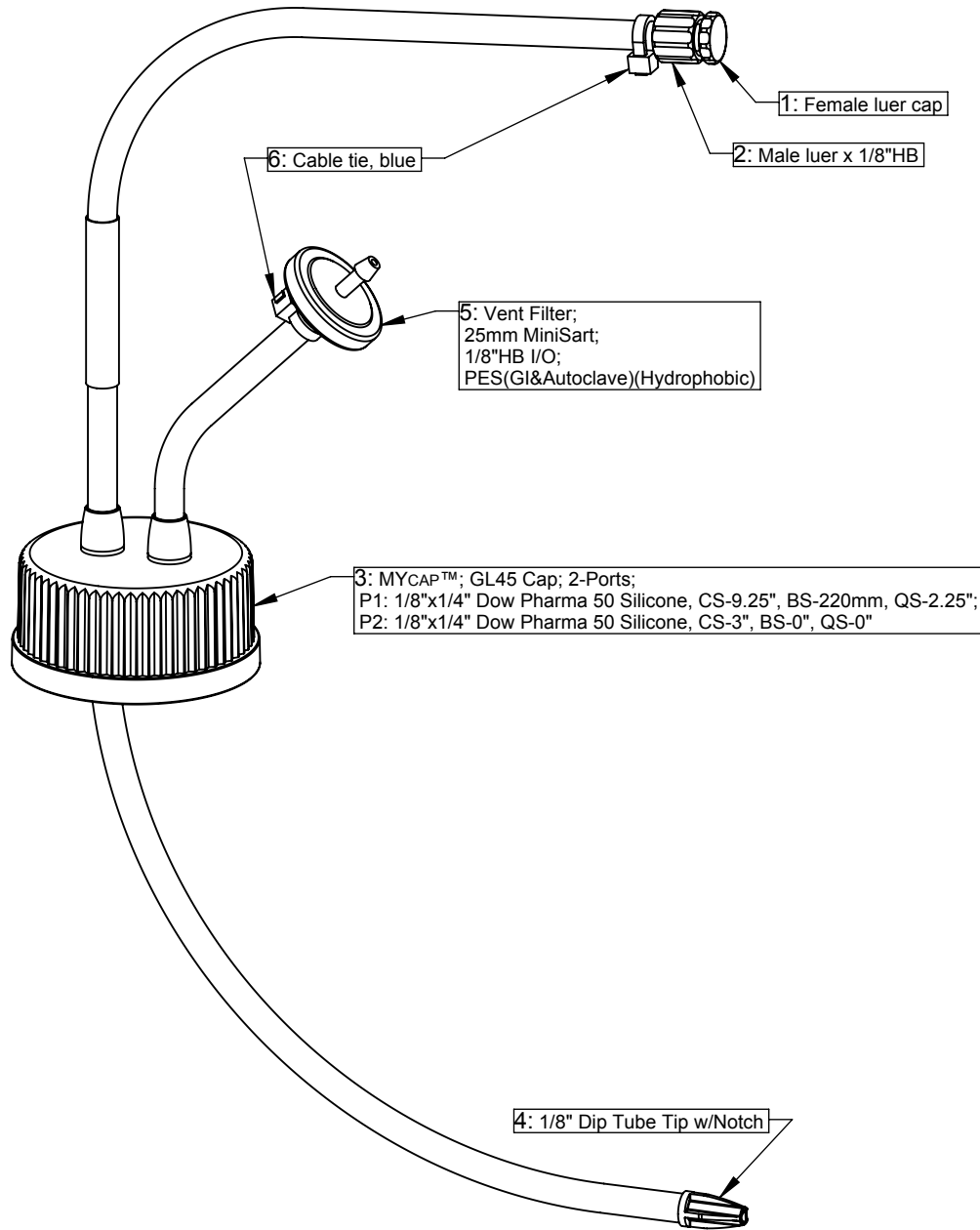


1000ml Corning Pyrex Glass, 1395-1000



Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-0099	-ADCF Polypropylene
2	1	AP-0031	-ADCF Polypropylene
3	1	AP-3021	-ADCF Silicone
4	1	AP-2170	-ADCF Polypropylene
5	1	AP-2767	-Polypropylene, PES
6	2	AP-133	-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. Individually bagged then double bagged, 5 per bag.
3. Fluid contact materials meet criteria for USP Class VI Testing. Consult Validation Guide for additional tests and details.
4. Drawing not to scale.

PROPRIETARY AND CONFIDENTIAL

The information contained herein is the sole property of AllPure Technologies, Inc. Reproduction or distribution is prohibited without written consent from AllPure Technologies, Inc.

Title: MYCAP™; GL45 Cap; 2-Ports;
 1/8"x1/4" Silicone w/Dip Tube;
 QS on inlet port - 1000ml Corning Glass Bottle;
 NOT IRRADIATED



80 Progress Ave
 New Oxford, PA 17350
 (717) 624-3241
 (717) 624-3051 (fax)

Customer Approval: Approval Not Required

Approval Date:

Part Number: MYCAP 13951000C

Drawing Number: MYCAP 13951000C

Revision: 0

Released By:

Release Date: