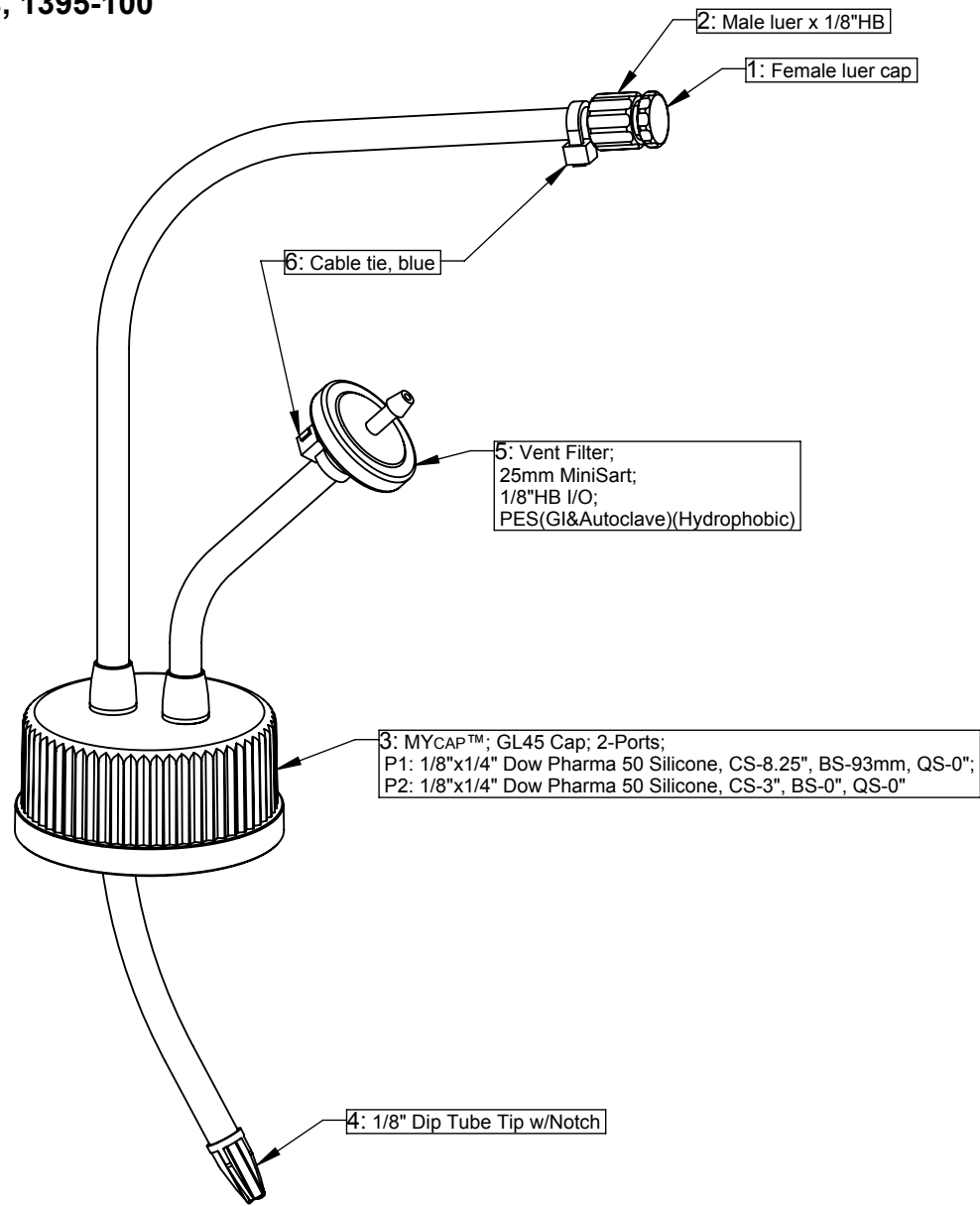


# 100ml Corning Pyrex Glass, 1395-100



### Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-99	-ADCF Polypropylene
2	1	AP-31	-ADCF Polypropylene
3	1	AP-2778	-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 - 100ml 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 13950100A

Drawing Number: MYCAP 13950100A

Revision: 0

Released By:

Release Date: