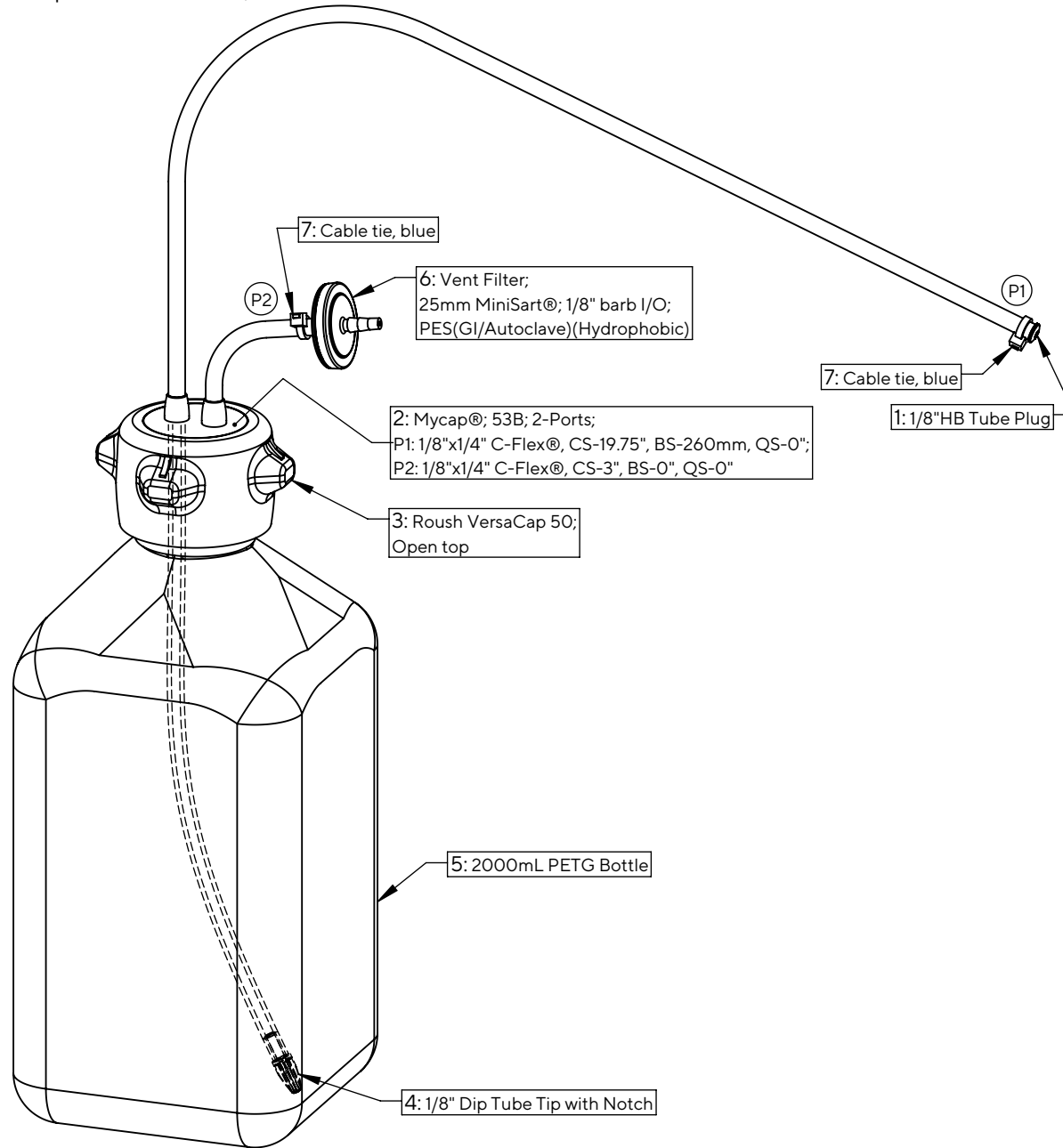


2000mL Nalgene Square Media Bottle, PETG



Bill of Materials:

Item	Qty	AP Num.	Material
1	1	AP-1181	-Polypropylene
2	1	AP-4396	-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"; AP-4544 to AP-4396
- 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.

SARTORIUS

Sartorius Stedim North America, Inc
80 Progress Ave
New Oxford, PA 17350

www.sartorius.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MYCAP NGPG2000B

Drawing Number: MYCAP NGPG2000B

Revision: 1

Released By:

Release Date:

Title:

Mycap®; 2000mL PETG; 2-Ports;
1/8"x1/4" C-Flex® w/Dip Tube