

Thermo Scientific Nalgene Volumetric Flasks

Cat. No. 4000, polypropylene - translucent

Cat. No. 4001, polymethylpentene - transparent

These flasks are:

- Guaranteed to meet accuracy requirements of volumetric flasks, Class B; ASTM Standard E288 and ISO Standard 384
- Individually calibrated gravimetrically TC/TD at 20°C
- Break-resistant
- Can be autoclaved, but accuracy will be affected
- Supplied with autoclavable, linerless, one-piece PP screw closure that won't trap contaminants; provides leakproof service

Guidelines for Use

Before using for the first time and after each use, wash in a warm detergent solution and rinse thoroughly. This will help ensure the accuracy of your flask. **Do not use abrasive cleaners or wire brushes.**

Before autoclaving, make sure that the flask is cleaned and rinsed thoroughly and just set cap or closure on top of the container without engaging the threads.

Make sure closure is tightly screwed onto flask before mixing its contents.

Chemical Resistance

These flasks resist most chemicals at room temperature, including most mineral acids and alkalies, even in concentrated form. They also resist many organic solvents, including ethyl alcohol, acetone and methyl ethyl ketone.

Do not store aromatic, aliphatic, or halogenated hydrocarbons or their derivatives in these flasks.

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC

© 2009 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

www.nalgenelabware.com

Asia: China Toll-free: 800-810-5118 or 400-650-5118;
India: +91 22 6716 2200, India Toll-free: 1 800 22 8374;
Japan: +81 3 3816 3355 Other Asian countries: 65 68729717
Europe: Austria: +43 1 801 40 0; Belgium: +32 2 482 30 30;
Denmark: +45 4631 2000; France: +33 2 2803 2180;
Germany: +49 6184 90 6940, Germany Toll-free: 08001-536 376;
Italy: +39 02 02 95059 or 434-254-375;
Netherlands: +31 76 571 4440;
Nordic/Baltic countries: +358 9 329 100;
Russia/CIS: +7 (812) 703 42 15;
Spain/Portugal: +34 93 223 09 18;
Switzerland: +41 44 454 12 12;
UK/Ireland: +44 870 609 9203
North America: USA/Canada +1 585 586 8800;
USA Toll-free: 800 625 4327
South America:
USA sales support: +1 585 899 7198
Countries not listed: +49 6184 90 6940 or
+33 2 2803 2180 8-0402-89 1109

Thermo
SCIENTIFIC