Care of Nalgene labware

General Cleaning

For most general applications, plastics may be washed in a mild (non-alkaline is recommended) detergent, followed by a rinse with tap water, and then distilled water. Do not use abrasive cleaners or scouring pads on any plastic labware. Do not use strong alkaline cleaning agents with polycarbonate.

Autoclaving

To prevent baking contaminants onto the surface of the plastic, all items should be carefully cleaned and rinsed in distilled water before autoclaving. Do not autoclave containers (except those made of fluorocarbons) containing detergent or wetting solutions.



A Always completely disengage threads or remove closure before autoclaving.

Polystyrene, polyvinyl chloride, nylon, styrene, acrylonitrile, acrylic, and low- and high-density polyethylene are not autoclavable under any conditions.

Society of Plastics Industry (SPI) Recycling Codes are molded into the bottom of all injection blow-molded or extrusion blow-molded bottles, 500 ml and larger.

For further information on how to safely and effectively use and care for your Nalgene labware, consult the current Nalgene Labware Catalog.

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

www.nalgerelabware.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; Notherlands: +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

8-0401-47 0609 +33 2 2803 2180

Thermo Scientific Nalgene **Bottles**



Thermo

NALGENE'



The anatomy of a Nalgene bottle



Nalgene bottles



Narrow-mouth Boston round

12 sizes from 4 ml to 2 L in: FEP, black FEP, FLPE, HDPE, amber HDPE, LDPE, PFA, PP, PVC



Wide-mouth round

7 sizes from 30 ml to 1.5 L in: LDPE, FEP, FLPE, HDPE, amber HDPE, PMP, PP



Straight-side wide-mouth jars 7 sizes from 15 ml to 1 L in:

clear, gray, blue and green PC, PMP, PP



Mason jars

4 sizes from 500 ml to 3 L in: PP



Rectangular

5 sizes from 125 ml to 2 L in: HDPE and amber HDPE



Wide-mouth square

5 sizes from 60 ml to 4.3 L in: HDPE and PP



Narrow-mouth square

6 sizes from 30 ml to 2 L in: HDPE, PC, PETG, PP