

Thermo Scientific Nalgene Calibrated Ware Guarantee

Nalgene® **Graduated Cylinders** meet the accuracy requirements of ASTM E 1272 Class B “Cylinders, Graduated Laboratory, Glass” (except Cat. Nos. 3663-2000 and -4000) and all requirements of ISO Standard 6706 “Plastic Laboratory Ware - Graduated Measuring Cylinders.”

Nalgene **Volumetric Flasks** meet the accuracy requirements of ASTM E 288, Class B, “Volumetric Flasks” and ISO Standard 384.

Nalgene **Burets** meet the accuracy requirements of ASTM E 287.

Nalgene **Double-Scale Pharmaceutical Graduates** 125mL (8 oz.) and larger conform to the volume and accuracy requirements of ASTM E 1094.



Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.



© 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

8-0404-02 0609

Thermo
S C I E N T I F I C