



Thermo Scientific Nalgene Graduated Cylinders Polypropylene (PP)

Cat. No. 3662

This Nalgene® graduated cylinder meets the accuracy requirements of ASTM Class B, E1272 "Cylinder, Graduated, Laboratory, Glass"; and all requirements of ISO Standard 6706 "Plastic Laboratory Ware – Graduated Measuring Cylinders."

Features

- Large, easy-to-read graduations
- Strong blue plastic base provides long life; large size for stability
- Generous pour spout
- Non-wetting so there is no meniscus; reading is easier; calibrated to contain/to deliver at 20°C
- Excellent chemical resistance, especially for short-term contact
- Durable; never slippery, even when wet

Use Guidelines

To obtain accurate results you MUST clean your PP cylinders thoroughly before each use. Nalgene L900 Liquid Detergent is recommended. The cylinders can be cleaned using automatic dishwashers, chemically sterilized or autoclaved, however, autoclaving may affect accuracy.



Visit www.thermoscientific.com product resources pages for complete product use guidelines.

For more information, see the current Nalgene Labware Catalog, or contact Technical Support at:
technical.nalgene@thermofisher.com

© 2011 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

www.thermoscientific.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: +81-3-5826-1616; 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 95059 552 or 434-254-375; Netherlands: +31 76 571 4440;
Nordic/Baltic countries: +358 9 329 10200; 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 2000

8-0402-93 0411

Thermo
SCIENTIFIC