

Thermo Scientific Nalgene Graduated Cylinders

Polymethylpentene (PMP)

Cat. No. 3663

Nalgene® graduated cylinders meet the accuracy requirements of ASTM Class B, E 1272, "Cylinder, Graduated, Laboratory, Glass" (except Cat. Nos. 3663-2000 and -4000); and all requirements of ISO Standard 6706 "Plastic Laboratory Ware – Graduated Measuring Cylinders".

Features:

- Wrap-around major graduations for easy reading
- 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 20°C
- Excellent chemical resistance, especially for short-term contact
- Durable
- Never slippery, even when wet
- Transparent

Use Guidelines

To obtain accurate results clean your PMP cylinders thoroughly before each use. The cylinder can be cleaned using automatic dishwashers, chemically sterilized or autoclaved and will tolerate temperatures as high as 135°C, however, autoclaving may affect accuracy.

For more information on these and other products, refer to 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.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 8-0402-92 0609

NALGENE™

Thermo
SCIENTIFIC

Thermo
SCIENTIFIC