

NALGENE



Thermo Scientific Nalgene Transparent Burets

(Acrylic body; PMP stopcock and tip; Teflon $^{\ensuremath{\circledast}}$ TFE plug) Cat. No. 3650

Nalgene[®] burets are individually calibrated and guaranteed to meet accuracy requirements of ASTM E287 Class B. They are compatible with most dilute acids and bases commonly used as titrants (except perchloric acid). They are unaffected by dilute mineral acids and bases, except hydrofluoric acid and ammonium hydroxide. The pH range of 1 to 14 is ideal, and up to 5% wt/wt of acids or 30% bases can be used.

These burets are not intended for use with alcohols or other organic solvents.

The body of the buret is molded of clear acrylic. While resilient, it may crack or shatter from harsh impact or other abuse.

The stopcock body and tip are made of polymethylpentene (PMP). The self-lubricating Teflon TPE stopcock plug will not seize and will never need greasing.

Cleaning

For accuracy in reading, always clean the buret before use, including the first time. Use a mild soap or detergent solution.

Do not use abrasive cleansers, wire brushes or acid cleaning solutions.

A rinse of 1% acetic acid solution will clean burets used with NaOH solution. Thoroughly rinse with distilled water afterwards. The stopcock plug can be removed for easy cleaning by squeezing and slipping off the retaining O-ring.

Do not autoclave the buret.

For further information, see the current Nalgene Labware Catalog or contact Technical Service at the numbers listed below.

© 2010 Thermo Fisher Scientific Inc. All rights reserved. Teflon is a registered trademark of DuPont. All other 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 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
 8-0400-80 0210