

Thermo Scientific Nalgene Polycarbonate Carboys and Large Bottles with Gasketed Closure

Cat. No. DS2205-0020;

Cat. No. 2251-all sizes; Cat. No. 2317-all sizes;

Cat. Nos. DS2127-0030, DS2127-0250,
and DS2127-2000

Autoclavable Nalgene® polycarbonate carboys and large bottles are supplied with a white polypropylene closure. Available closure sizes are 38-430 (38 mm), 53B (53 mm) or 83B (83 mm). The 53B and 83B closures include a gasket to assure leakproof service, especially during shipment. Firm hand-tightening is all that is needed for a positive seal.

The gasket has chemical resistance comparable to polycarbonate. This allows you to safely expose the gasket to chemicals that are compatible with the polycarbonate container. For complete chemical resistance information, consult the current Nalgene Labware Catalog.

Autoclave the sealing gasket while installed in the closure. Completely disengage threads or remove closure before autoclaving.

If the gasket becomes separated from the closure, re-install with the narrow side down (against the closure). Be sure that the gasket is securely seated in the closure before use.

For replacement gaskets, order part number

71-2160-0053 (53-mm closure) or

71-2162-1830 (83-mm closure).

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC

© 2010 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 3816 3355 Other Asian countries: 65 68729717

Europe: Austria: +43 1 801 40 0; Belgium: +32 53 73 42 41;
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 8-0403-80 1210

Thermo
SCIENTIFIC