

Thermo Scientific Nalgene Polycarbonate Containers

This Nalgene® container is precision-molded of polycarbonate resin (PC). Extremely tough and durable, PC is glass-clear and lightweight. These autoclavable containers are ideal for storing, preparing and dispensing distilled water, compatible reagents or media. They are excellent for preparing or using large-scale cultures, especially where light is required (e.g. algae). All items except the sanitary carboy include linerless, non-contaminating screw closures of polypropylene (PP). Select items include a TPE gasket in the closure.

Chemical Resistance

PC will resist aliphatic hydrocarbons, alcohols, vegetable oils, neutral and acid salts and dilute acids. DO NOT expose PC to strong bases, DMSO, halogenated hydrocarbons, aromatic hydrocarbons and their derivatives, esters, ethers, ketones or amines. These chemicals will attack PC containers.

Cleaning

PC containers can be cleaned with any mild, non-alkaline detergent, even in a laboratory dishwasher. DO NOT use abrasives, such as scouring powders or steel wool, or any alkaline detergent.

Sterilizing

Before autoclaving, be sure to rinse any residual chemicals or detergents thoroughly from PC containers. Before autoclaving, remove closure from container to prevent implosion or collapse of bottle or carboy when cooling. Always use minimum autoclaving temperatures and cycle times. DO NOT exceed 121°C/15 psi. Constant exposure to water at elevated temperatures will eventually damage PC labware. PC labware can be gas or chemically sterilized, but NEVER dry heat sterilized.

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC

Temperature Limits

PC bottles or carboys: -135°C to 135°C

PP screw closures: 0°C to 135°C

TPE gasket -50°C to 121°C

Note: Polycarbonate bottles and carboys are not suitable for vacuum service.

For more information on these and other products, refer to the current Nalgene Labware Catalog, or contact Technical Support at the numbers shown below.

© 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

8-0401-99 1109

Thermo
SCIENTIFIC