

# Thermo Scientific Nalgene Beakers made of Teflon® PFA

Contact us for Sales and Service thermoscientific.com/contactus \*Contact information contained within this document may be incorrect.

Cat. No. 1510, DS1511

These Nalgene® beakers are made of Teflon PFA (perfluoroalkoxy). They are extremely durable and chemical resistant.

### **Temperature Limits and Chemical Resistance**

Teflon PFA Beakers have the widest temperature range of all Nalgene

fluoropolymer products: -270°C to +250°C. Because of their excellent mechanical properties at low temperatures, they are well-suited to cryogenic work. At room temperature, these beakers are resistant to most chemicals, including strong oxidizing mineral acids such as Aqua Regia, 98% sulfuric acid, 50% chromic acid, 70% nitric acid, 85% phosphoric acid, 48% hydrofluoric acid and 35% hydrochloric acid.

Teflon PFA is non-cytotoxic, based on USP and ASTM biocompatibility testing standards utilizing MEM elution techniques on a WI28 human diploid cell line.

Caution: Do not put these beakers in a flame or on a hot plate.

## Cleaning

Wash in a warm, soapy detergent solution, followed by a rinse with tap water, then distilled water. Do not use brushes, abrasive cleaners or scouring pads.

## They can be washed in a labware washing machine if the following procedures are observed:

- 1. **Do not** wash in a machine equipped with brushes.
- In machines using a high-pressure water spray, place the beakers in a basket and cover them with a screen or cover similar to that used in test tube baskets. Otherwise, the water pressure may cause the beakers to tumble and become scratched.
- Teflon PFA beakers placed on spindles should be weighted or covered. If spindles are made of uncoated metal, a section of vinyl tubing placed over them will cushion and protect them.

#### Sterilizing

Before autoclaving, rinse the beakers thoroughly in distilled water. They may also be chemically sterilized. Beakers can be dry heat sterilized occasionally at 170°C (338°F) for 60 minutes, but this will cause discoloration and possible loss of mechanical strength. Do not radiation sterilize.

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** 











Printed on Rolland Enviro100 Print, which contains 100% post-consumer fiber, manufactured using renewable biogas energy and is certified EcoLogo, Processed Chlorine Free and FSC Recycled.

© 2011 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

