

Thermo Scientific

Nalgene Beakers Made of Teflon PFA

Cat. Nos. 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.
2. 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.
3. 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.

© 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 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

8-0405-66 1210

Thermo
SCIENTIFIC