

Thermo Scientific Nalgene Desiccator

Cat. No. 5311-0250

Features and benefits

- Transparent, unbreakable polycarbonate (PC) construction allows an undistorted view of contents, as well as desiccant levels and colors
- Designed for vacuum or non-vacuum use; this desiccator is designed to hold a vacuum of 28in/Hg (711 mm) for 24 hours
- Silicone O-ring assures an airtight seal; no grease is needed
- Much lighter weight than glass desiccators
- High-domed cover and shape of base provide a generous amount of inside space and stability on the benchtop; dimensions are 33 cm (13 in.) high x 27.9 cm (11 in.) O.D.
- Side-arms on cover and base for maximum convenience

For vacuum use

1. Attach 1/4 in. to 3/8 in. I.D. vacuum tubing (such as Nalgene® Cat. No. 8000-0145) to the barbed end of the enclosed white stopcock. Insert stopcock in the side-arm on the cover or base. Seal the other side-arm with one of the thermoplastic elastomer (TPE) caps provided.
2. Connect the tubing to a vacuum source.
3. Put a dry desiccant in the base. We recommend Drierite® or silica gel.
4. Put a desiccator plate in place. Any standard 230-mm plate can be used. We recommend the unbreakable Nalgene plate (Cat. No. 5312-0230), made of a ceramic-metal composite. Hot crucibles will not harm this plate.
5. Place crucibles or other containers on plate, close cover, and slowly apply a vacuum.
6. If the desiccator is used for a very light ash in crucibles, release the vacuum gently by opening the stopcock slowly.
7. To remove the cover, hold the knob and twist slightly while lifting. If it doesn't come off easily, the vacuum is not completely broken.

For non-vacuum use

1. Close stopcock or remove stopcock assembly, and seal side-arms on cover and base with the TPE caps provided.
2. Place desiccant, plate and samples in desiccator as described on front.

CAUTION:

- This desiccator is designed for room-temperature use only, whether used with or without vacuum.
- DO NOT expose the desiccator directly to the following chemicals or their vapors: strong bases, esters, aromatic or halogenated hydrocarbons or ketones. These chemicals will attack polycarbonate.
- NEVER autoclave the desiccator.

Cleaning

Use a warm soap or neutral pH detergent solution and rinse thoroughly. NEVER use organic solvents, strongly alkaline cleaning agents or abrasives.

For more information contact Nalgene Technical Service at the numbers listed below.

© 2009 Thermo Fisher Scientific Inc. All rights reserved. Drierite is a registered trademark of W. A. Hammond Drierite Company. All other 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

8-0403-96 0909

Thermo
SCIENTIFIC