

Nalgene® Mr. Frosty® Cryo 1°C Freezing Containers

Cat. No. 5100-0001, -0036, -0050

Polycarbonate container, high-density polyethylene closure, tube holder, foam insert



The Nalgene Mr. Frosty Cryo 1°C Freezing Containers are inexpensive, reusable and convenient devices for freezing biological samples at the recommended controlled rate of cooling, -1°C/minute. Isopropyl alcohol is necessary to achieve the recommended rate of cooling.

DIRECTIONS:

1. Remove high-density polyethylene tube holder and foam insert from polycarbonate unit.
DO NOT discard foam insert.
2. Add 100% isopropyl alcohol to the fill line on the Mr. Frosty container. **DO NOT overfill.**
3. Carefully replace foam insert and tube holder.
4. Place tubes containing sample into holes in tube holder.
5. Place unit in bottom of -70°C to -80°C mechanical freezer. Leave undisturbed for a minimum of 4 hours.
6. Remove frozen tubes from unit and place in a permanent, long-term storage freezer (e.g. -130°C or below).

NOTE: DO NOT store full jar on its side. Store at room temperature when not in use. Replace alcohol every fifth use. Clean with non-abrasive, non-alkaline detergent such as Thermo Scientific Nalgene L-900 Detergent (Cat. No. 900-4000). The tube holder floats with tubes in place and can be used as a floating rack when thawing samples.

© 2010 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

www.thermoscientific.com

8-0404-28 1110