

Designed for use with the entire portfolio of Thermo Scientific Nunc and Nalgene cryogenic storage tubes, Nalgene® Mr. Frosty® freezing containers provide simple cooling of samples at the optimal rate to preserve cells.

Thermo Scientific Nalgene Mr. Frosty Freezing Container



- **Optimal cooling:** provides critical, repeatable $-1^{\circ}\text{C}/\text{minute}$ cooling rate required for successful cell cryopreservation and recovery.
- **Easy and safe:** simply fill to the designated line with 100% isopropyl alcohol and place in your mechanical freezer. Each unit is labeled with graphic, step-by-step instructions so there is no fumbling for manuals or printouts. The samples never touch the alcohol - no risk of label or printing removal.
- **Flexibility:** with units and inserts available for 1.0, 1.2, 1.8, 2.0, 3.6, 4.5 and 5.0 ml tubes, the Mr. Frosty freezing container provides the widest tube compatibility available. Holds 12 or 18 tubes.

Product Specifications



Thermo Scientific Nalgene Mr. Frosty Freezing Container

Materials: Polycarbonate jar, high-density polyethylene closure and tube holder, foam insert

Specifications	5100-0001	5100-0036	5100-0050
Closure Size, mm/in:	120 / 4-3/4	120 / 4-3/4	120 / 4-3/4
Height, mm/in:	86 / 3-7/16	151 / 5-5/16	151 / 5-5/16
Diameter, mm/in:	117 / 4-5/8	117 / 4-5/8	117 / 4-5/8
Compatible Nalgene and Nunc Tubes, ml	1.0 - 2.0	3.6	4.5 - 5.0
Number of tubes held	18	12	12
Volume Isopropyl Alcohol, ml	250	500	500

Ordering Guide

Cat. No.	Description	Qty/Case
5100-0001	Mr. Frosty, -1°C/minute Controlled Rate Freezing Vessel, 1.0-2.0 mL tubes	1 unit
5100-0036	Mr. Frosty, -1°C/minute Controlled Rate Freezing Vessel, 3.6 mL tubes	1 unit
5100-0050	Mr. Frosty, -1°C/minute Controlled Rate Freezing Vessel, 4.5-5.0 mL tubes	1 unit

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

Asia: China Toll-free: 800-810-5118, 400-650-5118, India toll-free: 1 800 22 8374,

India: +91 22 6716 2200, Japan: +81 3 3816 3355

Europe: Austria +43 1 801 40 0, Belgium +32 53 73 42 41, Denmark +45 4631 2000; France +33 2 2803 2180,

Germany national toll free 08001-536 376, Germany international +49 6184 90 6940, Italy +39 02 02 95059 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

Other Asian countries 65 68729717

Countries not listed: +49 6184 90 6940 or +33 2 2803 2180

www.thermoscientific.com/mrfrosty

SSLSPMRFROSTY 1110

Thermo
SCIENTIFIC