

Nalgene Mr. Frosty Freezing Container

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

- **Optimal cooling:** provides critical, repeatable $-1^{\circ}\text{C}/\text{minute}$ cooling rate required for successful cell cryopreservation and recovery.
- **Easy and convenient:** 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 - minimizing 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 Specification

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 information

Description	Quantity	Cat. No
Mr. Frosty, $-1^{\circ}\text{C}/\text{minute}$ Controlled Rate Freezing Vessel, 1.0-2.0 mL tubes	1 unit	5100-0001
Mr. Frosty, $-1^{\circ}\text{C}/\text{minute}$ Controlled Rate Freezing Vessel, 3.6 mL tubes	1 unit	5100-0036
Mr. Frosty, $-1^{\circ}\text{C}/\text{minute}$ Controlled Rate Freezing Vessel, 4.5-5.0 mL tubes	1 unit	5100-0050

Learn more at thermofisher.com/mrfrosty

thermo scientific