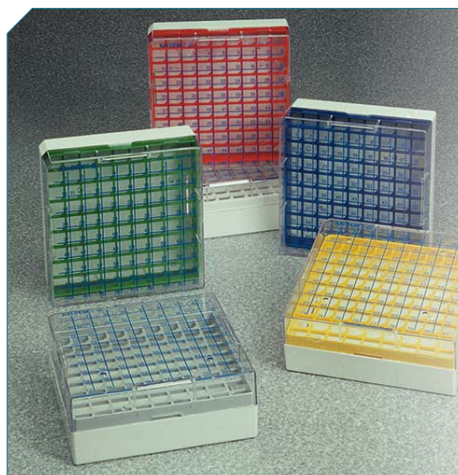


Nalgene Cryobox and Nunc Cryogenic Storage Tubes

Buy 3 cases of Nalgene Polycarbonate 2 in. CryBoxes and get 1 case of Nunc 1.8mL Polypropylene Cryo Tubes at no additional cost!

Break-resistant, durable polycarbonate Thermo Scientific™ Nalgene™ Cryoboxes.

- Durable plastic cryoboxes for short- or long-term storage
- Keyed lid on plastic cryoboxes prevent misalignment
- Printed grid for easy inventory management
- Plastic boxes autoclavable for decontamination and usable at temperature range from -196°C to +121°C



Thermo Scientific™ Nunc™ Cryotubes are a high quality, reliable solution for highly-regulated applications, irreplaceable high-value content storage and tissue, cell, blood or other materials for biobanking.

- Sterility to SAL 10⁻⁶
- Made from noncytotoxic and nonmutagenic
- Large writing area and graduations for easy sample identification

Limit 5 redemptions per lab

Available through December 31, 2017

Redeem now at thermofisher.com/labessentials

Cat. No.	Description	No./Per Case
5026-0909	Nalgene Cryobox, 1.2/2ML PC;9X9 ARRAY	24
377267	Nunc 1.8 mL Cryotube, Round, Internal Thread, Starfoot	50/1800

Terms and Conditions:

This promotion is open to customers in the U.S. and Canada only. Valid for orders placed between July 1, 2017 and December 31, 2017. Redemptions for the promotional case and proof of purchase must be submitted by December 31, 2017. Promotion offers for Nalgene Economy Bottles are limited to 4 redemptions per laboratory. Promotions offers for Nalgene RTU Wash Bottles, Nalgene Sterilizing Pans, Nalgene Rapid-Flow Filter Units, Nalgene Syringe Filters, Nalgene Analytical Filters, Nunc Delta-Treated Products, Nunc Serological Pipettes and Conical Tubes, Nalgene PETG Erlenmeyer Flasks and Nalgene Cryoboxes and Nunc CryoTubes are limited to 5 redemptions per laboratory. Cannot be combined with other discounts or promotion. Offer void where prohibited, licensed or restricted by federal, state, provincial, or local laws or regulation or agency/institutional policy. Other restrictions may apply. U.S. customers acknowledge that this offer may include a discount or other price reduction that must be properly and accurately accounted for and reported by customer in accordance with all federal and state laws, including without limitation the federal antikickback law (42 C.F.R. § 1320a-7b(b)(3)(A)) and regulations thereunder (42 C.F.R. §1001.952(h)).

For Research Use Only. Not for use in diagnostic procedures. © 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details. LPEMultiPromoH2 2017