

Matrix 2D Barcode Storage Solution

Purchase 1 Visionmate Wireless reader, 1 Handheld Decapper and 1 case of Matrix 0.5mL or 1.0mL Screw Top Tubes in barcoded latch racks and get 4 additional cases of tubes at no cost

Accelerate processing while ensuring sample integrity with Thermo Scientific™ 2D barcoded tubes and equipment.

- **VisionMate Wireless Reader** enables sample tracking on the go; in the field, cold room, or around the lab. Store over 600,000 codes in the internal memory!
- **Matrix Handheld Decapper** is a practical addition to benchtop processing of screw top storage tubes. Process a full rack of Matrix tubes in less than one minute.
- **Matrix 0.5 mL or 1.0mL 2D Barcoded Screw top tubes** have a permanent 2D barcode on the rack bottom to enable orientation detection and ID when using automated storage equipment. Ideal for biobanking. Recommended for use at temperatures from -80°C to -196°C.



Cat. No.	Description	No./Per Case
3122-V2	VisionMate Wireless Tube Reader	1
4105MAT	Matrix Handheld Decapper, 8-Ch, Matrix	1
3741-BR	Matrix 1.0 mL Barcoded Screw top tubes w/caps installed, sterile	5 racks
3744-BR	Matrix 0.5 mL, 2D Barcoded Screw top tubes w/caps installed, sterile	5 racks

Available through December 31, 2017

Redeem now at thermofisher.com/labessentials

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. **LPMultiPromoH2 2017**