



Enhance sample tracking with Matrix Tri-coded Storage Tubes

Avoid taking chances when storing and sharing precious samples

With the rapid growth of biobanking, infectious disease research and development, cancer research and molecular diagnostics, storing valuable samples is more important than it has ever been within the scientific community.

Thermo Scientific™ Matrix™ storage tubes feature permanent high-contrast 2D, 1D linear, and human-readable barcodes, offering three levels of identification to safely track valuable samples. Matrix storage tubes come in sterile and non-sterile options and can withstand many years of freezer storage and repeated freeze/thaw cycles.

Matrix storage tubes enable:

- **Elimination of secondary labeling steps**—2D-barcoded Thermo Scientific™ Matrix™ ScrewTop and Open-Top Storage Tubes are available preprinted with 1D and human-readable codes that match the tubes' 2D barcodes.
- **Enhanced secure tracking**—Permanent 2D barcode is laser-etched onto the base of every tube to securely identify and track samples. The linear and human-readable barcodes match this unique 2D barcode, allowing 1D scanning or visual identification of samples across satellite sites without 2D readers.
- **Stringent quality control**—Every 2D barcoded storage tube is scanned to ensure readability and uniqueness.



Ordering information

Description	Max working volume (µL)	Cap color	Quantity	Sterile/ non-sterile	Cat. No.
500 µL Matrix ScrewTop Tri-coded Tubes					
2D white patch 1D linear barcode, V-bottom	450	Clear	Bulk, 480/case	Sterile	3743-WP1D
2D white patch 1D linear barcode, V-bottom, barcoded rack	450	Clear	5 racks of 96/case	Sterile	3744-WP1D-BR
2D white patch 1D linear barcode, V-bottom	450	Clear	5 racks of 96/case	Sterile	3744-WP1D
2D white patch, V-bottom, barcoded rack	450	Clear	5 racks of 96/case	Sterile	3744-WP-BR
2D white patch, V-bottom	450	Clear	5 racks of 96/case	Sterile	3744-WP
2D white patch 1D linear barcode, V-bottom, barcoded rack	450	Clear	5 racks of 96/case	Non-sterile	3744NS-WP1D-BR
2D white patch 1D linear barcode, V-bottom, barcoded rack	450	No cap	5 racks of 96/case	Sterile	3745-WP1D-BR
2D white patch 1D linear barcode, V-bottom	450	No cap	5 racks of 96/case	Sterile	3745-WP1D
2D white patch 1D linear barcode, V-bottom, barcoded rack	450	No cap	5 racks of 96/case	Non-sterile	3745NS-WP1D-BR
1.0 mL Matrix ScrewTop Storage Tubes					
2D white patch 1D linear barcode, V-bottom	925	Clear	Bulk, 480/case	Sterile	3740-WP1D
2D white patch 1D linear barcode, V-bottom, barcoded rack	925	Clear	5 racks of 96/case	Non-sterile	3741NS-WP1D-BR
2D white patch 1D linear barcode, V-bottom, barcoded rack	925	Clear	5 racks of 96/case	Sterile	3741-WP1D-BR
2D white patch 1D linear barcode, V-bottom	925	Clear	5 racks of 96/case	Sterile	3741-WP1D
2D, white patch, V-bottom, barcoded rack	925	Clear	5 racks of 96/case	Sterile	3741-WP-BR
2D white patch, V-bottom	925	Clear	5 racks of 96/case	Sterile	3741-WP
2D white patch 1D linear barcode, V-bottom	925	No cap	5 racks of 96/case	Non-sterile	3742NS-WP1D
2D white patch 1D linear barcode, V-bottom, barcoded rack	925	No cap	5 racks of 96/case	Sterile	3742-WP1D-BR
2D white patch, V-bottom, barcoded rack	925	No cap	5 racks of 96/case	Sterile	3742-WP-BR
2D white patch 1D linear barcode, V-bottom	925	No cap	5 racks of 96/case	Sterile	3742-WP1D
Description	Total capacity (µL)	Max working volume (µL)	Quantity	Sterile/ non-sterile	Cat. No.
Matrix 2D Barcoded Open-Top Storage Tubes					
0.75 mL, 2D white patch 1D linear barcode, V-bottom	750	670	10 latch racks of 96 tubes/case	Sterile	3732-WP1D
0.75 mL, 2D white patch 1D linear barcode, V-bottom	750	670	10 latch racks of 96 tubes/case	Non-sterile	3732NS-WP1D
1.4 mL, 2D white patch 1D linear barcode, V-bottom	1,400	1,260	Bulk, case of 1,000 tubes	Non-sterile	3790-WP1D
1.4 mL, 2D white patch 1D linear barcode, V-bottom, barcoded rack	1,400	1,260	10 latch racks of 96 tubes/case	Non-sterile	3791-WP1D-BR
1.4 mL, 2D white patch 1D linear barcode, V-bottom, barcoded rack	1,400	1,260	10 latch racks of 96 tubes/case	Sterile	3792-WP1D-BR
Accessories					
Matrix SeptraSeal caps	NA	NA	10 mats of 96 caps/case	Non-sterile	4463
SeptraSeal caps	NA	NA	10 mats of 96 caps/case	Sterile	4464
SeptraSeal, pre-split caps	NA	NA	10 mats of 6 caps/case	Sterile	4466

Find out more at thermofisher.com/tricodedtubes