

AMPIGENE[®] Reverse Transcriptase

AMPIGENE[®] Reverse Transcriptase is a highly thermostable reverse transcriptase for the most challenging RNA sample types

AMPIGENE[®] Reverse Transcriptase is a modified MMLV reverse transcriptase engineered for superior cDNA synthesis speed, specificity, and full-length cDNA product.

Ordering Information

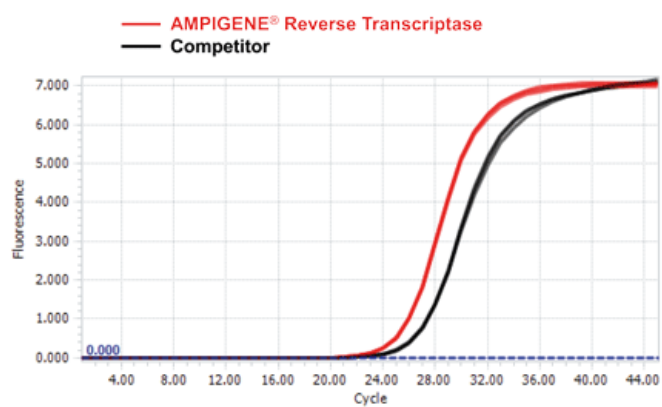
[Order Online »](#)

ENZ-NUC137-10001	10000U
ENZ-NUC137-40001	40000U

Manuals, SDS & CofA

[View Online »](#)

- High cDNA yields from as little as 4 pg total RNA
- Advanced buffer chemistry including Mg and dNTPs
- Thermostable reverse transcriptase 45°C to 55°C
- Reduced RNase H activity



AMPIGENE® Reverse Transcriptase provides higher sensitivity and yield (red amplification plot) compared to a competitor's product (black amplification plot).

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Contents AMPIGENE[®] RT (200 U/μL) with RNase inhibitor, 5x
AMPIGENE[®] RT Buffer

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)