

# AMPIGENE<sup>®</sup> dNTP

## Mix

Ultra-pure, stable, and suitable for a variety of PCR applications

AMPIGENE<sup>®</sup> dNTP Mix is an ultrapure mix of deoxyribonucleotide triphosphates to use as building blocks for a wide variety of PCR applications.

It can withstand multiple freeze-thaw cycles and shown to retain 95% dNTPs in triphosphate form after 5 weeks at room temperature.

AMPIGENE<sup>®</sup> dNTP Mix has been tested on many molecular biology applications including standard PCR, fast PCR, 1-step PCR, high-fidelity PCR, real-time PCR, lon-range PCR, MDA, RDA, LAMP, and DNA sequencing.

### Ordering Information

[Order Online »](#)

ENZ-NUC102-0500	500µl
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)

- Ultra-pure at ≥99%
- Stable after multiple freeze-thaw cycles
- Ideal for standard PCR, real-time PCR, high-fidelity PCR, 1-step PCR, and long range PCR

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Notes      For a wide variety of PCR applications

Concentration      10mM

Purity      ≥99%

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)