

Bio-N⁶-ddATP

Bio-N⁶-ddATP (Biotin-N⁶-(2',3'-dideoxy)adenosine-5' triphosphate) can be used for 3'-end labeling of single-stranded or double-stranded DNA. It is especially useful as a substrate for terminal transferase for the 3'-terminal addition of a single Bio-N⁶-ddAMP. Biotin-labeled oligonucleotides produced by this reaction are suitable as hybridization probes. Biotin-labeled cDNA produced by this reaction can be used in nucleic acid array assays. Biotin-labeled DNA in hybridization assays can be detected by a reporter molecule linked to streptavidin, avidin, or anti-biotin antibody. Such a complex can be detected directly, e.g. by the excitation of a fluorophore conjugated to streptavidin, or indirectly, e.g. using an enzyme conjugate that can produce an insoluble colored precipitate. Bio-N⁶-ddATP is also a substrate for *Taq* DNA Polymerase and other DNA polymerases.

Citations: 3

[View Online »](#)

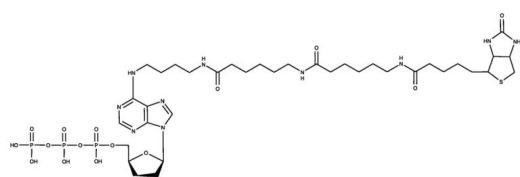
Ordering Information

[Order Online »](#)

ENZ-42809	25nmol
-----------	--------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Biotin-N6-(2',3'-dideoxy)-adenosine-5'-triphosphate

Appearance Clear, colorless liquid.

Concentration 1mM

Extinction Coefficient 17,000 M⁻¹ cm⁻¹ (265 nm, pH 7.0)

Formula C₃₆H₆₁N₁₀O₁₅P₃S

Formulation Liquid. Solution in water.

MW 998.9 (free acid)

Purity ≥93% (HPLC)

Purity Detail Contains <5% Bio-N6-ddADP.

Technical Info / Product Notes Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.

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-
fr@enzolifesciences.com](mailto:info-fr@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)