

Bio-7-dATP

Bio-7-dATP (Biotin-7-2'-deoxyadenosine-5'-triphosphate) can replace dATP in reactions in which it serves as a substrate for *E. coli* DNA polymerase (holoenzyme and Klenow fragment), T4 and Taq DNA polymerases, reverse transcriptase (from AMV and M-MuLV) and terminal transferase. Bio-7-dATP can be used to produce biotinylated DNA probes in a variety of labelling reactions including nick translation, random prime labelling, cDNA labelling and 3'-end labelling. The resulting biotin-labeled probe can be used in a variety of hybridization applications including Southern blots, Northern blots, or dot blots. The probes can also be used for *in situ* hybridization procedures on fixed cells and tissues. The biotinylated probes have been shown to hybridize to homologous nucleic acid at the same rate and to the same extent as non-biotinylated probes. The hybridized biotinylated DNA probes can be detected by their interaction with biotin-binding proteins, such as avidin, streptavidin or antibodies coupled to fluorescent dyes or color-producing enzymes.

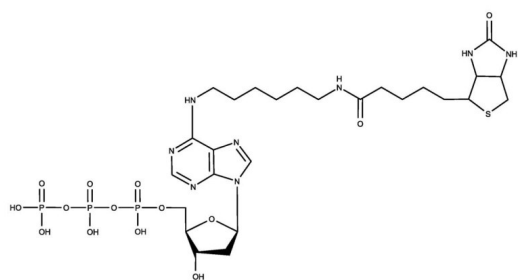
Ordering Information

[Order Online »](#)

ENZ-42819	50nmol
-----------	--------

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-7-2'-deoxyadenosine-5'-triphosphate

Appearance Clear, colorless liquid.

Concentration 1mM

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

Formula C₂₆H₄₃N₇O₁₄P₃S

Formulation Liquid. Solution in water.

MW 816.7 (free acid)

Purity ≥93% (HPLC)

Purity Detail Contains <5% Bio-7-dADP.

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)