

# 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.

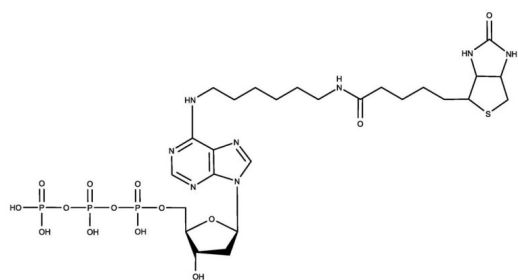
## Ordering Information

[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 <sup>-1</sup> cm <sup>-1</sup> (265 nm, pH 7.0)
Formula	C <sub>26</sub> H <sub>43</sub> N <sub>7</sub> O <sub>14</sub> P <sub>3</sub> S
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.



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)