

# Bio-16-ddUTP . tetralithium salt

Bio-16-ddUTP (Biotin-16-2',3'-dideoxyuridine-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-16-ddUMP. 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-16-ddUTP is also a substrate for *Taq* DNA Polymerase and other DNA polymerases.

Citations: 1

[View Online »](#)

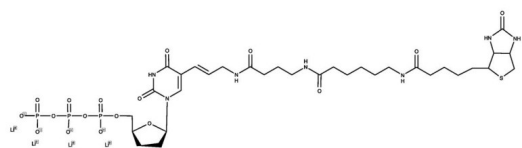
## Ordering Information

[Order Online »](#)

ENZ-42813	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-16-2',3'-dideoxyuridine-5'-triphosphate . 4Li
Appearance	Clear, colorless liquid.
Concentration	1mM
Extinction Coefficient	10,700 M <sup>-1</sup> cm <sup>-1</sup> (240 nm, pH 7.0)
Formula	C <sub>32</sub> H <sub>48</sub> Li <sub>4</sub> N <sub>7</sub> O <sub>17</sub> P <sub>3</sub> S
Formulation	Liquid. Solution in water.
MW	927.7 . 28.0
Purity	≥93% (HPLC)
Purity Detail	Contains <5% Bio-16-ddUDP.

## 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)