

# Biotin-16-dCTP, S pack

## Ordering Information

[Order Online »](#)

JBS-NU-809-BIO16-S	200µl
--------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Biotin-16-Propargylamino-dCTP,  $\gamma$ -[N-(Biotin-6-amino-hexanoyl-6-aminobutanoyl)]-5-(3-propargylamino)-2'-deoxycytidine-5'-triphosphate, Triethylammonium salt

**Appearance** colorless to slightly yellow solution

**Application Notes** Incorporation into DNA/cDNA by – PCR with Taq polymerase  
Nick Translation with DNase I/ DNA Polymerase I  
Primer Extension with Klenow exo-  
3'-End Labeling with Terminal deoxynucleotidyl Transferase (TdT)  
Reverse Transcription with MMLV Reverse Transcriptase

**Couple Type** Modified nucleotides

**Formula**  $C_{32}H_{51}N_8O_{17}P_3S$  (free acid)

**Formulation** sterile solution in 10 mM Tris-HCl

**MW** 944.78 (free acid)

**Purity**  $\geq 95\%$  (HPLC)

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)