

Biotin-16-ddUTP (L pack)

Ordering Information

[Order Online »](#)

JBS-NU-253-BIO16-L	5x25µl
--------------------	--------

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-(5-aminoallyl)-ddUTP

Appearance colorless to slightly yellow solution

Application Notes Technique for nanoscale patterning of functional molecules on the surface of a DNA origami
3'-end labeling of oligonucleotides with Terminal Transferase, recombinant and labeling DNA
MALDI-TOF MS-based SNP typing assay
Biotinylation of in vitro transcribed tmRNA during surface plasmon resonance (SPR) technique
Use in the standard nick translation reaction for tumor DNA labelling during comparative genomic hybridization

CAS 188755-42-2 (Tetralithium salt), 748750-05-2 (free acid)

Concentration 1.0 mM – 1.1 mM

Couple Type Modified nucleotides

Formula $C_{32}H_{52}N_7O_{17}P_3S$ (free acid)

Formulation sterile solution in 10 mM Tris-HCl

MW 931.78 (free acid)

Purity ≥ 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com