

Biotin-16-UTP

Ordering Information

[Order Online »](#)

JBS-NU-821-BIO16	30µl
------------------	------

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-UTP, γ -[N-(Biotin-6-amino-hexanoyl-6-aminobutanoyl)]-5-(3-aminoallyl)-uridine-5'-triphosphate, Triethylammonium salt

Appearance colorless to slightly yellow solution

Application Notes Microarray based hybridization detection
Northern Blot

Couple Type Modified nucleotides

Formula $C_{32}H_{52}N_7O_{19}P_3S$ (free acid)

Formulation sterile solution in 10 mM Tris-HCl

MW 963.78 (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)