

Biotin-14-dCTP, S pack

Biotin-14-dCTP is enzymatically incorporated into DNA/cDNA as substitute for its natural counterpart dCTP. The resulting Biotin-labeled DNA/cDNA probes are subsequently detected using streptavidin conjugated with horseradish peroxidase (HRP), alkaline phosphatase (AP), a fluorescent dye or agarose/magnetic beads. Optimal substrate properties for Nick Translation are ensured by a 14-atom linker attached to the N4 position of cytosine. For PCR incorporation experiments e.g. with Taq polymerase Biotin-11-dCTP (#NU-809-BIOX) or Biotin-16-dCTP (#NU-809-BIO16) are recommended whose Biotin moiety is attached to C5 position of cytidine via a 11-atom or 16-atom linker, respectively. Recommended Biotin-14-dCTP/dCTP ratio for Nick Translation: 50% Biotin-14-dCTP/ 50% dCTP Please note: The optimal final concentration of Biotin-14-dCTP may vary depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Biotin-14-dCTP/dCTP ratio is recommended.

Ordering Information

[Order Online »](#)

JBS-NU-956-BIO14-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 Bio-14-dCTP, Triethylammonium salt

Appearance colorless to slightly yellow solution

Application Notes Incorporation into DNA/cDNA by – Nick Translation with DNase I/ DNA Polymerase I
Primer Extension with Klenow fragment

Concentration 1 mM – 1.1 mM

Couple Type Modified nucleotides

Formula $C_{31}H_{54}N_7O_{16}P_3S$ (free acid)

Formulation sterile solution in 10 mM Tris-HCl

MW 905.78 (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