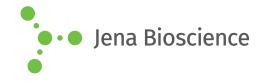
DATA SHEET





Biotin-16-dCTP

Biotin-16-Propargylamino-dCTP

γ-[N-(Biotin-6-amino-hexanoyl-6-aminobutanoyl)]-5-(3-propargylamino)-2'-deoxycytidine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-809-BIO16-S	200 μl (1 mM)
NU-809-BIO16-L	5 x 200 μl (1 mM)

Structural formula of Biotin-16-dCTP

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₂H₅₁N₈O₁₇P₃S (free acid)

Molecular Weight: 944.78 g/mol (free acid)

Exact Mass: 944.23 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: colorless to slightly yellow **Concentration:** 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 294 nm, ϵ 9.3 L mmol $^{-1}$ cm $^{-1}$ (Tris-HCl

pH 7.5

Applications:

Incorporation into DNA/cDNA by

- PCR with Taq polymerase in-house data
- Nick Translation with DNAse I/ DNA Polymerase I $^{\left[1,2\right] }$
- Primer Extension with Klenow exo^{- [3]}
- 3'-End Labeling with Terminal deoxynucleotidyl Transferase (TdT) [4]
- Reverse Transcription with MMLV Reverse Transcriptase [5]

Description:

Biotin-16-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 and thus labeling efficiency as well as an efficient detection of the Biotin moiety is ensured by a 16-atom linker attached to the C5 position of cytidine.

Recommended Biotin-16-dCTP/dCTP ratio for PCR: 50% Biotin-16-dCTP/ 50% dCTP

Please note: The optimal final concentration of Biotin-16-dCTP may very depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Biotin-16-dCTP/dTTP ratio is recommended.

Selected References:

[1] Sugimoto et al. (1999) Isolation and mapping of a putative b subunit of human ATP synthase (ATP-BL) from human leukocytes. DNA Res. 6 (1):29.

[2] Sreedharan *et al.* (1995) Structure, expression, and chromosomal localization of the type I human vasoactive intestinal peptide receptor gene. *Proc. Natl. Acad. Sci. U S A.* **92 (7)**:2939.

[3] Yue *et al.* (1997) Apoptosis and necrosis in the newborn piglet brain following transient cerebral hypoxia-ischaemia. *Neuropathol. Appl. Neurobiol.* **23 (1)**:16.

[4] Zammetto et al. (2005) Unambiguous identification of expressed MAGE-A genes on a DNA microarry. Clin Chem **51**:2420.

[5] Grimmond *et al.* (2001) Expression Profiling with DNA Microarray's: A User's Perspective and Guide. In: DNA Microarrays: Gene Expression Applications (Jordan). Springer Verlag Berlin Heidelberg.