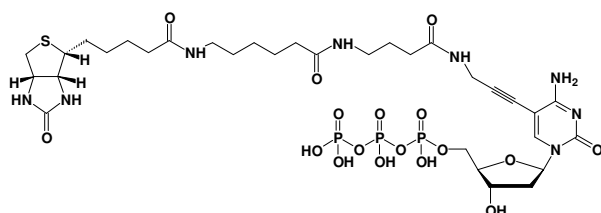


Biotin-16-dCTP

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

Cat. No.	Amount
NU-809-BIO16	200 μ l / (5 mM)



Cat. No.: NU-809-BIO16

Molecular Formula: C₃₂H₅₁N₈O₁₇P₃S (free acid)

Molecular Weight: 944.78 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{\max} 294 nm; ϵ 9300

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

Applications:

Nick translation^[1, 2]

DNA end labeling^[3, 4]

Primer extension^[5, 6]

Substrate:

DNA polymerase I ^[1, 2]

Terminal deoxynucleotidyltransferase^[3]

Klenow fragment^[4]

Taq Pol (inhouse data)

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 1 human vasoactive intestinal peptide receptor gene. *Proc. Natl. Acad. Sci. U S A.* **92 (7)**:2939.

[3] Aune *et al.* (1999) Costimulation reverses the defect in IL-2 but not effector cytokine production by T cells with impaired I κ B α degradation. *J. Immunol.* **162 (10)**:5805.

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

[5] De Rienzo *et al.* (2001) Loss of heterozygosity analysis defines a 3-cM region of 15q commonly deleted in human malignant mesothelioma. *Oncogene* **20 (43)**:6245.

[6] Lee *et al.* (1996) Loss of heterozygosity analysis defines a critical region in chromosome 1p22 commonly deleted in human malignant mesothelioma. *Cancer Res.* **56 (19)**:4297.