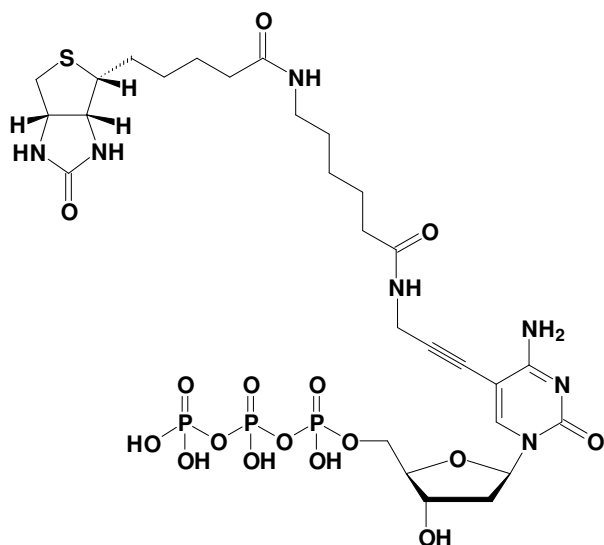


Biotin-11-dCTP

γ -[N-(Biotin-6-amino-hexanoyl)]-5-propargylamino-2'-deoxy-cytidine-5'-triphosphate,
Triethylammonium salt

Cat. No.	Amount
NU-809-BIOX	200 μ l / 5 mM



Cat. No.: NU-809-BIOX

Molecular Formula: C₂₈H₄₄N₇O₁₆P₃S (free acid)

Molecular Weight: 859.67 (free acid)

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

Spectroscopic properties: λ_{\max} 294 nm;
 ϵ 9,300 M⁻¹cm⁻¹ Tris HCl pH 7.5

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:

FISH^[1, 2]

SNP-analysis^[1]

Nick-translation^[2]

TUNEL^[3]

Selected References:

[1] Ye *et al.* (2001) Fluorescent microsphere-based readout technology for multiplexed human single nucleotide polymorphism analysis and bacterial identification. *Human Mutation* **17**:305.

[2] Backx *et al.* (2008) Direct fluorescent labelling of clones by DOP PCR. *Molecular Cytogenetics* **1**:3.

[3] Jones *et al.* (2003) Herpes simplex virus type 2 induces rapid cell death and functional impairment of murine dendritic cells in vitro. *J. Virology* **77**:11139.

Bishop *et al.* (2008) APOBEC3G Inhibits Elongation of HIV-1 Reverse Transcripts. *PLoS Pathogens* **4**:e1000231.