

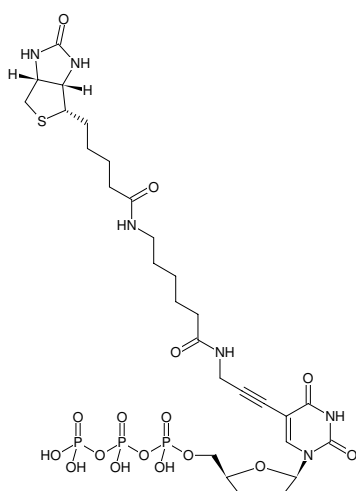


Biotin-11-ddUTP

Biotin-X-5-Propargylamino-ddUTP

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

Cat. No.	Amount
NU-1619-BIOX-S	25 μ l (1 mM)
NU-1619-BIOX-L	5 x 25 μ l (1 mM)



Structural formula of Biotin-11-ddUTP

For research use only!

Shipping: shipped on blue ice

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: C₂₈H₄₃N₆O₁₆P₃S (free acid)

Molecular Weight: 844.66 g/mol (free acid)

Exact Mass: 844.17 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: sterile solution in 10 mM Tris-HCl

Color: colorless to slightly yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{max} 290 nm, ϵ 13.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

SNP-analysis^[1]

DNA sequencing^[2]

3'end labeling^[3]

Selected References:

[1] Tong *et al.* (2002) Single nucleotide polymorphism detection by combinatorial fluorescence energy transfer tags and biotinylated dideoxynucleotides. *Nucleic Acids Res.* **30** (5):e19.

[2] Ju (2002) DNA sequencing with solid-phase-capturable dideoxynucleotides and energy transfer primers. *Anal Biochem.* **309** (1):35.

[3] Wang *et al.* (2009) Role of sgcR3 in positive regulation of enediyne antibiotic C-1027 production of *Streptomyces globisporus* C-1027. *BMC Microbiol.* **9**:14.

Edwards *et al.* (2001) DNA sequencing using biotinylated dideoxynucleotides and mass spectrometry. *Nucleic Acids Res.* **29** (21):E104.