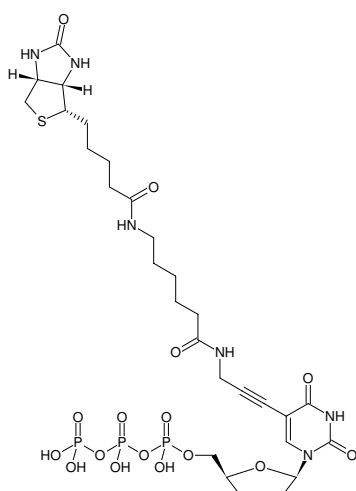


**Biotin-11-ddUTP**

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 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:** C₂₈H₄₃N₆O₁₆P₃S (free acid)**Molecular Weight:** 844.66 g/mol (free acid)**Exact Mass:** 844.17 g/mol (free acid)**Purity:** \geq 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 \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.