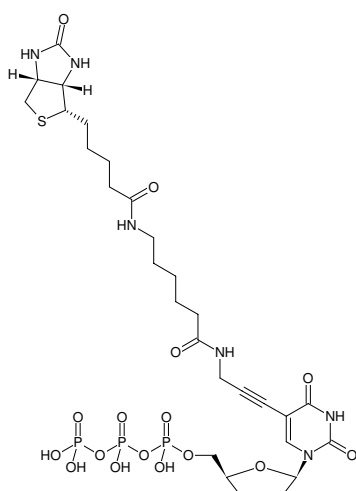


**Biotin-11-ddUTP**

Biotin-X-5-Propargylamino-ddUTP

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

Cat. No.	Amount
NU-1619-BIOX-S	25 $\mu$ l (1 mM)
NU-1619-BIOX-L	5 x 25 $\mu$ 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:** C<sub>28</sub>H<sub>43</sub>N<sub>6</sub>O<sub>16</sub>P<sub>3</sub>S (free acid)**Molecular Weight:** 844.66 g/mol (free acid)**Exact Mass:** 844.17 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** sterile clear aqueous solution in 10 mM Tris-HCl**Concentration:** 1.0 mM - 1.1 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  290 nm,  $\epsilon$  13.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**SNP-analysis<sup>[1]</sup>DNA sequencing<sup>[2]</sup>3'end labeling<sup>[3]</sup>**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.