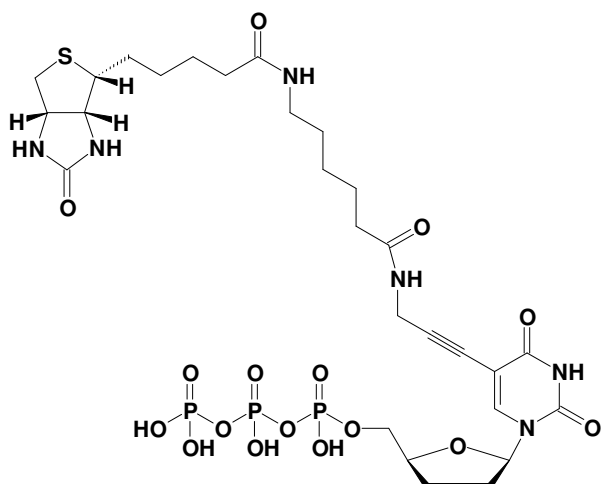


## Biotin-11-ddUTP

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

Cat. No.	Amount
NU-1619-BIOX	25 $\mu$ l / 5 mM



**Cat. No.:** NU-1619-BIOX

**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 (free acid)

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

**Spectroscopic properties:**  $\lambda_{\max}$  290 nm;  
 $\epsilon$  13,000 M<sup>-1</sup>cm<sup>-1</sup> 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:

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.