

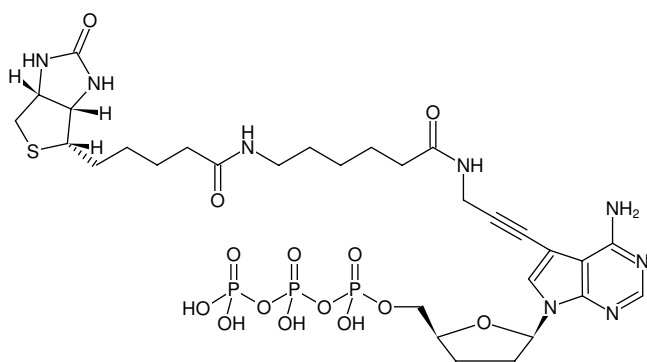
**Biotin-11-ddATP**

$\gamma$ -[N-(Biotin-6-amino-hexanoyl)]-7-propargylamino-2',3'-dideoxy-deaza-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1612-BIOX	25 $\mu$ l (1 mM)

**Selected References:**

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



Structural formula of Biotin-11-ddATP

**For general laboratory use.**

**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:** C<sub>30</sub>H<sub>45</sub>N<sub>8</sub>O<sub>14</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 866.71 g/mol (free acid)

**Exact Mass:** 866.20 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** 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:**  $\lambda_{\text{max}}$  280 nm,  $\epsilon$  12.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)