

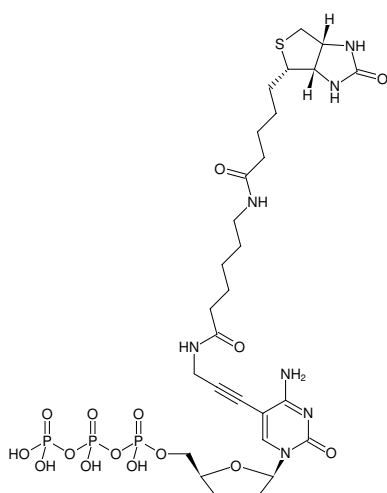
**Biotin-11-ddCTP**

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

Cat. No.	Amount
NU-850-BIOX-S	25 µl (1 mM)
NU-850-BIOX-L	5 x 25 µl (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-ddCTP

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: C₂₈H₄₄N₇O₁₅P₃S (free acid)

Molecular Weight: 843.67 g/mol (free acid)

Exact Mass: 843.18 g/mol (free acid)

Purity: ≥ 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 ±0.5

Spectroscopic Properties: λ_{max} 294 nm, ε 9.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)