

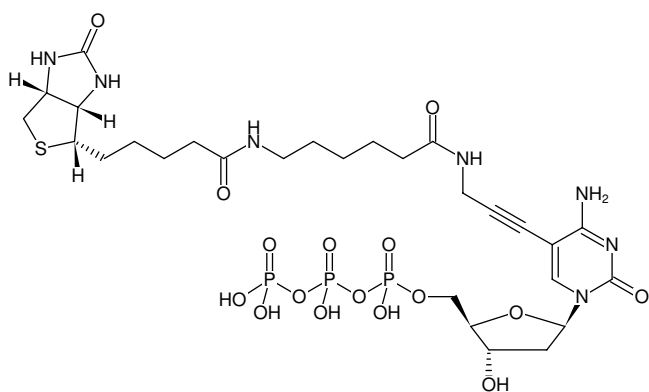


### Biotin-11-dCTP - high concentration

Biotin-X-5-Propargylamino-dCTP

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

Cat. No.	Amount
NU-809-BIOX-HC	500 $\mu$ l (5 mM)



Structural formula of Biotin-11-dCTP - high concentration

#### 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>28</sub>H<sub>44</sub>N<sub>7</sub>O<sub>16</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 859.67 g/mol (free acid)

**Exact Mass:** 859.18 g/mol (free acid)

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

**Form:** filtered solution (30 kDa) in 10 mM Tris-HCl

**Color:** colorless to slightly yellow

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{\max}$  294 nm,  $\epsilon$  9.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)