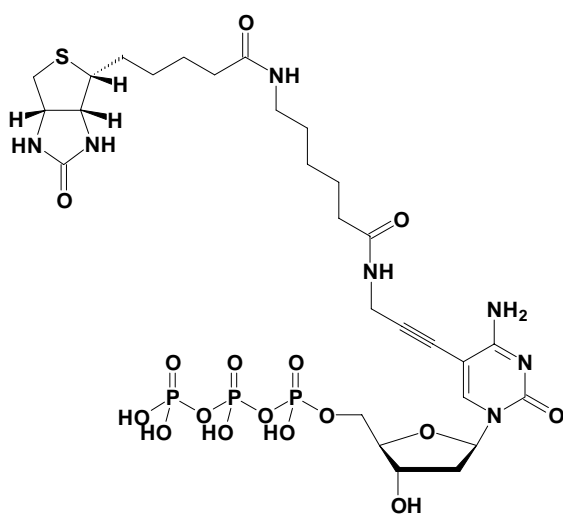


## Biotin-11-dCTP

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

Cat. No.	Amount
NU-809-BIOX	15 $\mu$ l / 5 mM



**Cat. No.:** NU-809-BIOX

**Molecular Formula:**  $C_{28}H_{44}N_7O_{16}P_3S$  (Anion)

**Molecular Weight:** 859.67 (Anion)

**Purity:** > 95%, clear aqueous solution, 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!**