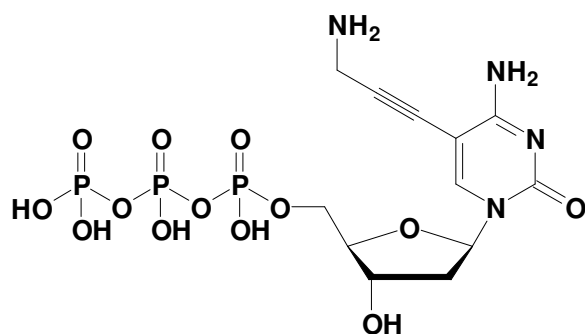


5-Propargylamino-dCTP

5-Propargylamino-2'-deoxy-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-809S	30 Units
NU-809L	150 Units



Cat. No.: NU-809

Molecular Formula: $C_{12}H_{19}N_4O_{13}P_3$ (free acid)

Molecular Weight: 520.22 (free acid)

Purity: > 95%, slightly yellowish aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 294 nm;
 ϵ 9,300 $M^{-1}cm^{-1}$ 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!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Cherkasov *et al.* (2010) New Nucleotide Analogues with Enhanced Signal Properties. *Bioconjugate Chem.* **21** (1):122.

Hobbs (1989) Palladium-catalyzed synthesis of alkynylamino nucleosides - a universal linker for nucleic-acids. *J. Org. Chem.* **54** (14):3420.