

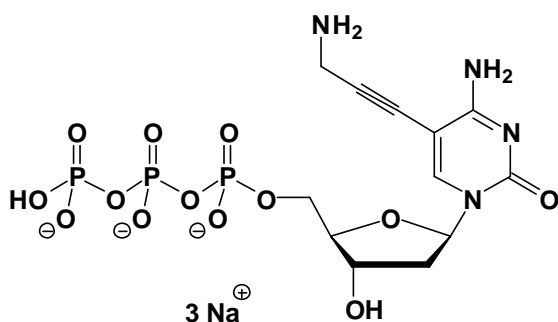
5-Propargylamino-dCTP

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

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

Selected References:

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



Cat. No.: NU-809

Molecular Formula: C₁₂H₁₆N₄O₁₃P₃ (Anion)

Molecular Weight: 517.19 (Anion)

Purity: > 95%, slightly yellowish 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!

1 unit = 1 μl of a 10 mM solution