



5-Propargylamino-dUTP

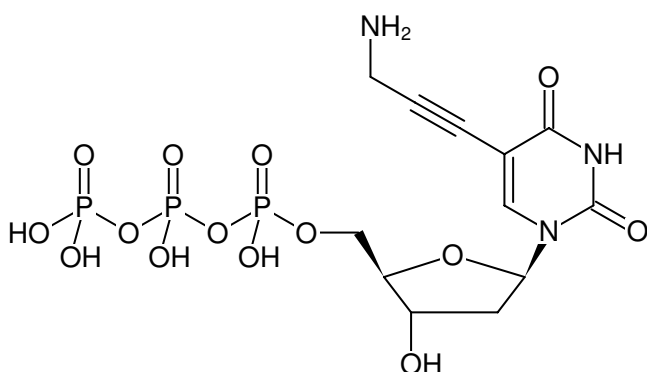
5-Propargylamino-2'-deoxyuridine-5'-triphosphate, Sodium salt
Aminopropargyl-dUTP

Cat. No.	Amount
NU-998S	10 µl (100 mM)
NU-998L	5 x 10 µl (100 mM)

Selected References:

Arango *et al.* (2018) Acetylation of Cytidine in mRNA Promotes Translation Efficiency. *Cell* **175** (7):1872.

Kuwahara *et al.* (2005) Comparison study on PCR amplification of modified DNA by using various kinds of polymerase and modified nucleoside triphosphates. *Nucleic Acids Symp. Ser. (Oxf)*. **49**:275.



Structural formula of 5-Propargylamino-dUTP

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: C₁₂H₁₈N₃O₁₄P₃ (free acid)

Molecular Weight: 521.20 g/mol (free acid)

Exact Mass: 521.00 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 289 nm, ε 10.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)