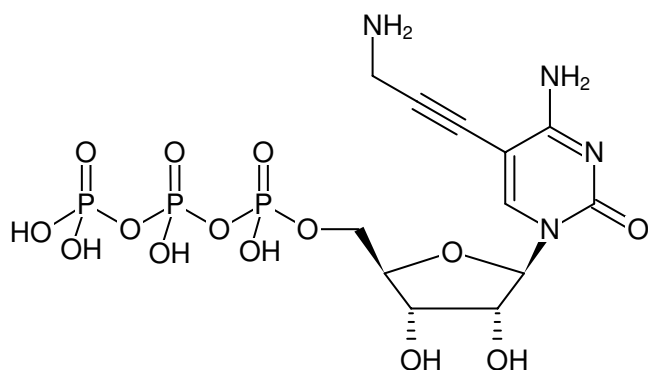


**5-Propargylamino-CTP**

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

Cat. No.	Amount
NU-831S	50 µl (50 mM)
NU-831L	5 x 50 µl (50 mM)



Structural formula of 5-Propargylamino-CTP

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:** C₁₂H₁₉N₄O₁₄P₃ (free acid)**Molecular Weight:** 536.01 g/mol (free acid)**Exact Mass:** 536.22 g/mol (free acid)**CAS#:** 150718-26-6**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 50 mM - 55 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 298 nm, ε 8.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)