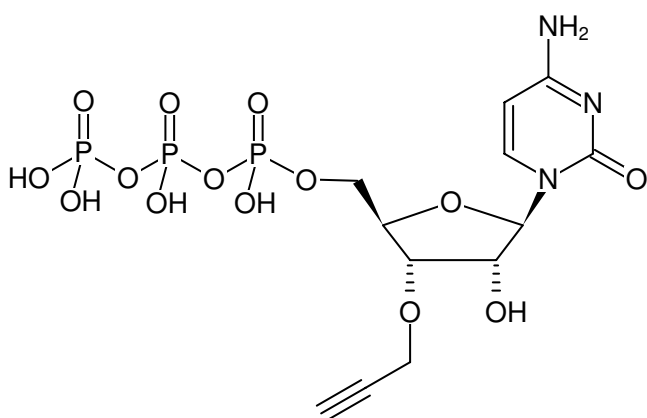


**3'-(O-Propargyl)-CTP**

3'-(O-Propargyl)-cytosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-947	1 μ mol



Structural formula of 3'-(O-Propargyl)-CTP

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** 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**Exact Mass:** 521.00 g/mol**Purity:** \geq 95 % (HPLC)**Form:** clear sticky semisolid**Spectroscopic Properties:** λ_{max} 271 nm, ϵ 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)