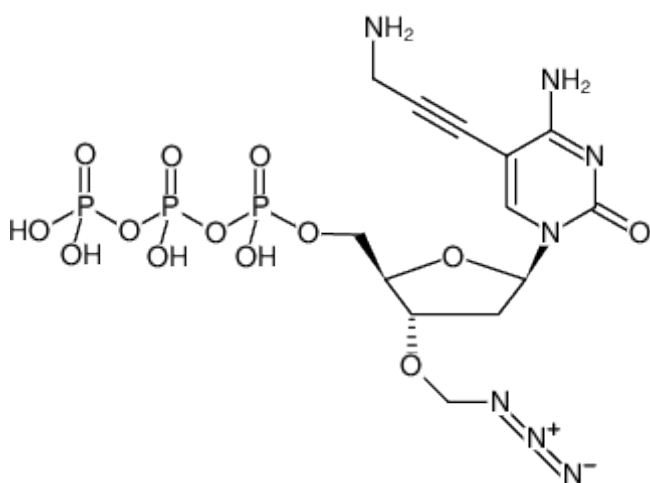




3'-O-Azidomethyl-5-Propargylamino-dCTP

3'-O-Azidomethyl-5-Propargylamino-2'-deoxycytidine-5'-triphosphate, Sodium salt
5-Propargylamino-3'-azidomethyl-dCTP

Cat. No.	Amount
NU-288-1	1 mg



Structural formula of 3'-O-Azidomethyl-5-Propargylamino-dCTP

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: C₁₃H₂₀N₇O₁₃P₃ (free acid)

Molecular Weight: 575.26 g/mol (free acid)

Exact Mass: 575.03 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 294 nm, ε 9.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)