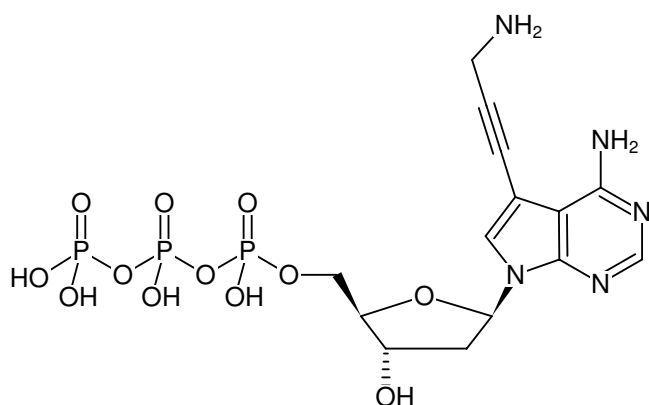




7-Deaza-7-propargylamino-dATP

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1611-1	1 mg (1,84 µmol)



Structural formula of 7-Deaza-7-propargylamino-dATP

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: C₁₄H₂₀N₅O₁₂P₃ (free acid)

Molecular Weight: 543.26 g/mol (free acid)

Exact Mass: 543.03 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: colorless to white

Spectroscopic Properties: λ_{abs} 280 nm, ε 12.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)