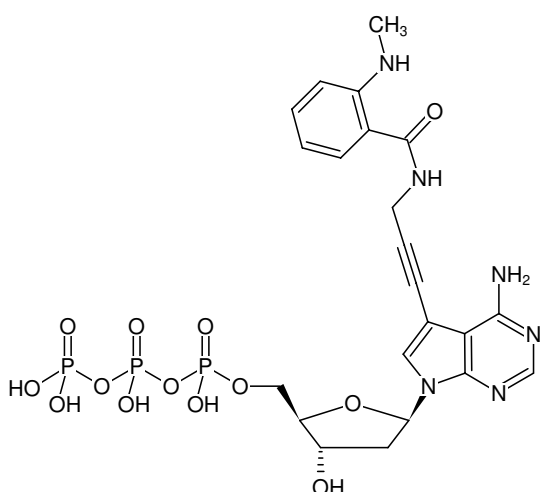




7-Propargylamino-7-deaza-dATP-MANT

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

Cat. No.	Amount
NU-1611-MNT	100 µl (5 mM)



Structural formula of 7-Propargylamino-7-deaza-dATP-MANT

For research use only!

Shipping: shipped on blue ice

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: C₂₂H₂₇N₆O₁₃P₃ (free acid)

Molecular Weight: 676.41 g/mol (free acid)

Exact Mass: 676.08 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 335 nm, λ_{em} 440 nm, ε 2.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)