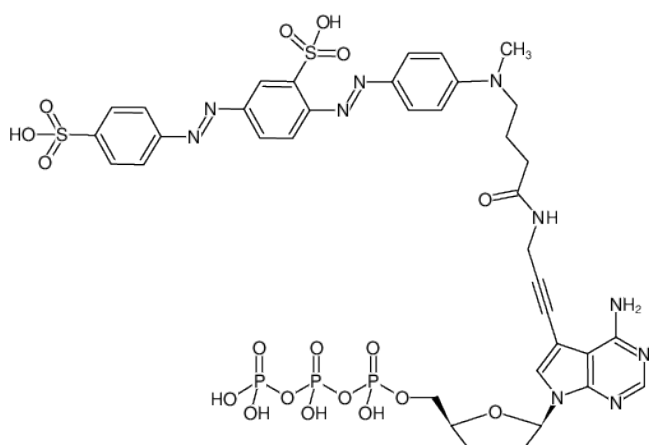




7-Propargylamino-7-deaza-ddATP-BHQ-10

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with BHQ-10™, Triethylammonium salt

Cat. No.	Amount
NU-1612-BHQ10	30 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-ddATP-BHQ-10

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: C₃₇H₄₁N₁₀O₁₈P₃S₂ (free acid)

Molecular Weight: 1070.83 g/mol (free acid)

Exact Mass: 1070.13 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: dark red

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 516 nm, ε 28.7 L mmol⁻¹ cm⁻¹ (PBS pH 7.0)