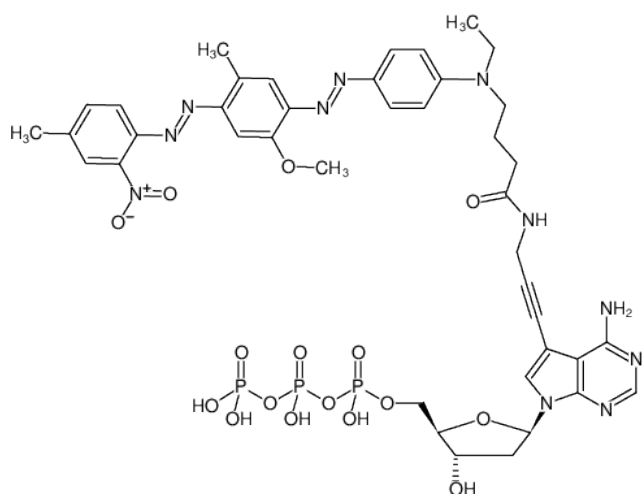




7-Propargylamino-7-deaza-dATP-BHQ-1

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with BHQ-1™, Triethylammonium salt

Cat. No.	Amount
NU-1611-BHQ1	30 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dATP-BHQ-1

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: C₄₁H₄₈N₁₁O₁₆P₃ (free acid)

Molecular Weight: 1043.82 g/mol (free acid)

Exact Mass: 1043.25 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: dark purple

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 534 nm, ε 34.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)