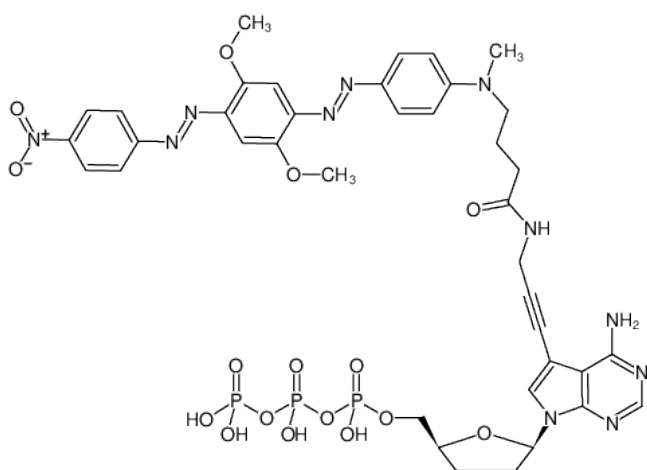




7-Propargylamino-7-deaza-ddATP-BHQ-2

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

| Cat. No. | Amount |
|--------------|--------------|
| NU-1612-BHQ2 | 30 µl (1 mM) |



Structural formula of 7-Propargylamino-7-deaza-ddATP-BHQ-2

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: C₃₉H₄₄N₁₁O₁₆P₃ (free acid)

Molecular Weight: 1015.76 g/mol (free acid)

Exact Mass: 1015.22 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: dark blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 579 nm, ε 38.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)