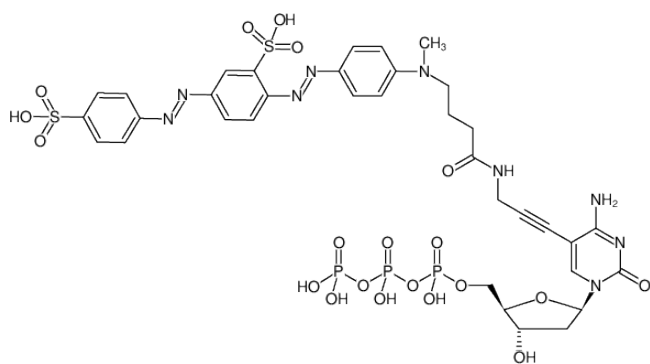




5-Propargylamino-dCTP-BHQ-10

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with BHQ-10™, Triethylammonium salt

Cat. No.	Amount
NU-809-BHQ10	30 µl (1 mM)



Structural formula of 5-Propargylamino-dCTP-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: C₃₅H₄₀N₉O₂₀P₃S₂ (free acid)

Molecular Weight: 1063.79 g/mol (free acid)

Exact Mass: 1063.10 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)