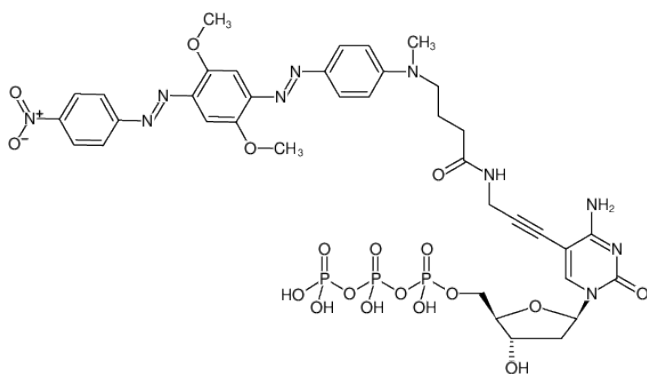


**5-Propargylamino-dCTP-BHQ-2**

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

Cat. No.	Amount
NU-809-BHQ2	30 µl (1 mM)



Structural formula of 5-Propargylamino-dCTP-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:** C₃₇H₄₃N₁₀O₁₈P₃ (free acid)**Molecular Weight:** 1008.72 g/mol (free acid)**Exact Mass:** 1008.20 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)