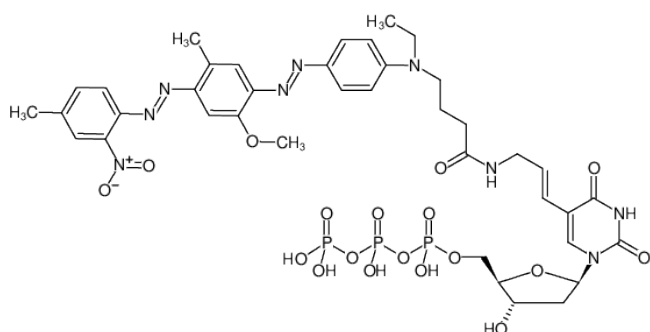




Aminoallyl-dUTP-BHQ-1

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with BHQ-1™, Triethylammonium salt

Cat. No.	Amount
NU-803-BHQ1	30 µl (1 mM)



Structural formula of Aminoallyl-dUTP-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: C₃₉H₄₈N₉O₁₈P₃ (free acid)

Molecular Weight: 1023.78 g/mol (free acid)

Exact Mass: 1023.23 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)