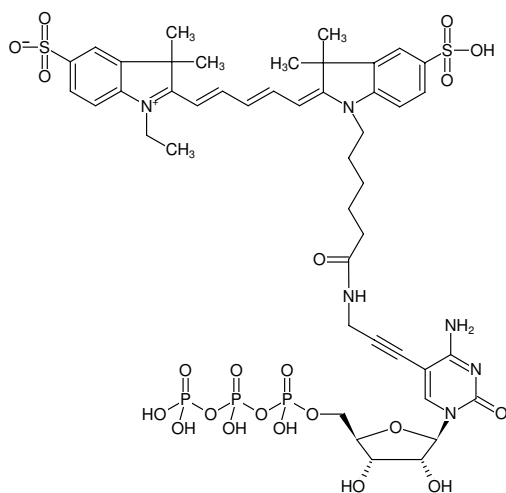




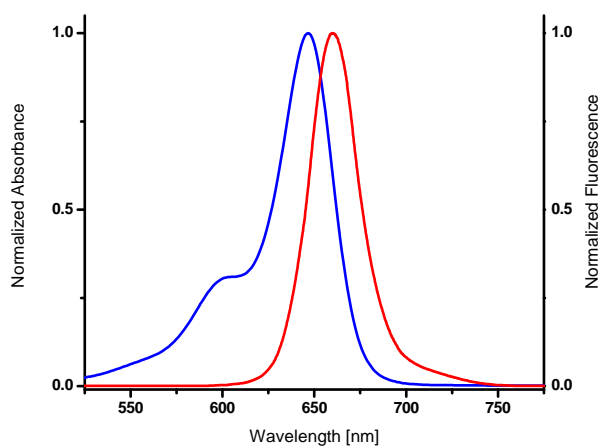
5-Propargylamino-CTP-Cy5 - high concentration

5-Propargylamino-cytidine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-831-CY5-HC	10 µl (10 mM)



Structural formula of 5-Propargylamino-CTP-Cy5 - high concentration



excitation and emission spectrum of Cy5

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: C₄₅H₅₇N₆O₂₁P₃S₂ (free acid)

Molecular Weight: 1175.01 g/mol (free acid)

Exact Mass: 1174.22 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 10 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)