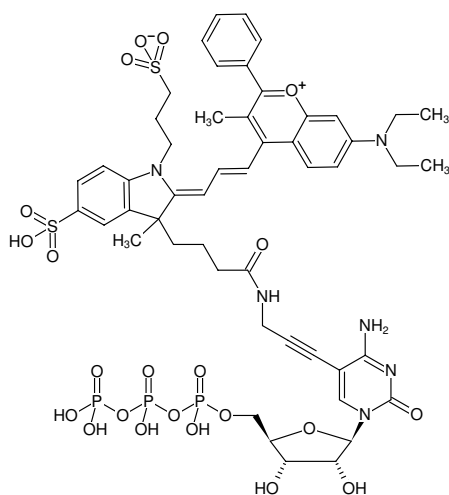




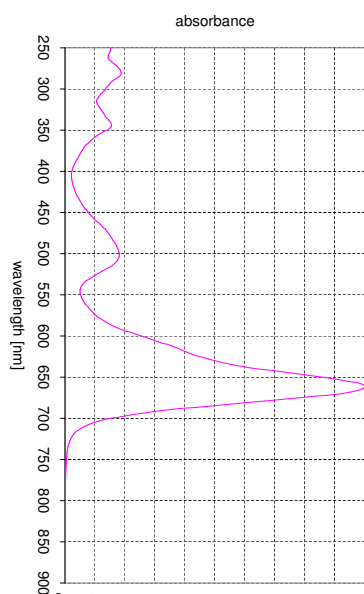
5-Propargylamino-CTP-DYQ-661

5-Propargylamino-cytidine-5'-triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-831-661Q	40 µl (1 mM)



Structural formula of 5-Propargylamino-CTP-DYQ-661



absorption spectrum of DYQ 661

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: C₅₁H₆₀N₆O₂₂P₃S₂ (Anion)

Molecular Weight: 1266.10 g/mol (Anion)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 662 nm, ε 104.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 662 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)

Please note: DYQ661 is not a dark quencher, λ_{em} 690 nm^[1]

Selected References:

[1] Tan *et al.* (2010) Direct detection of Delta9-tetrahydrocannabinol in saliva using a novel homogeneous competitive immunoassay with fluorescence quenching. *Anal. Chim. Acta.* **658 (2)**:187.