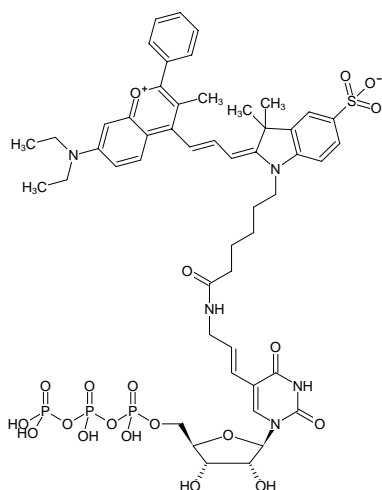




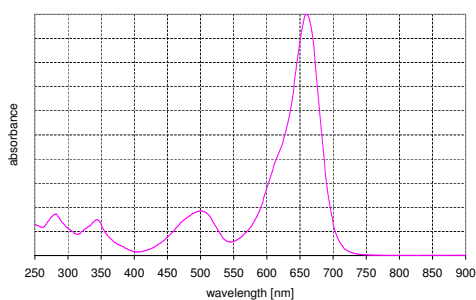
Aminoallyl-UTP - DYQ-660

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with DYQ 660, Triethylammonium salt

Cat. No.	Amount
NU-821-660Q	40 µl (1 mM)



Structural formula of Aminoallyl-UTP - DYQ-660



absorption spectrum of DYQ 660

For research use only!

Shipping: shipped on blue ice

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: C₅₁H₆₂N₅O₂₀P₃S (free acid)

Molecular Weight: 1190.05 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 660 nm; ε 109.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 660 nm; ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)