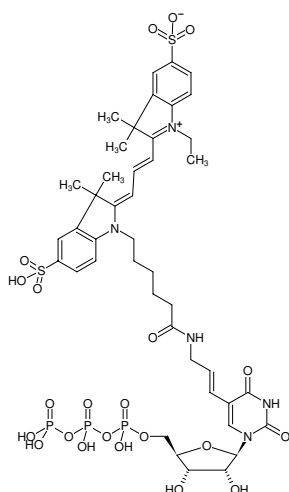




Aminoallyl-UTP-Cy3

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-821-CY3	30 µl (1 mM)



Structural formula of Aminoallyl-UTP-Cy3

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: C₄₃H₅₆N₅O₂₂P₃S₂ (free acid)

Molecular Weight: 1151.97 g/mol (free acid)

Exact Mass: 1151.21 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

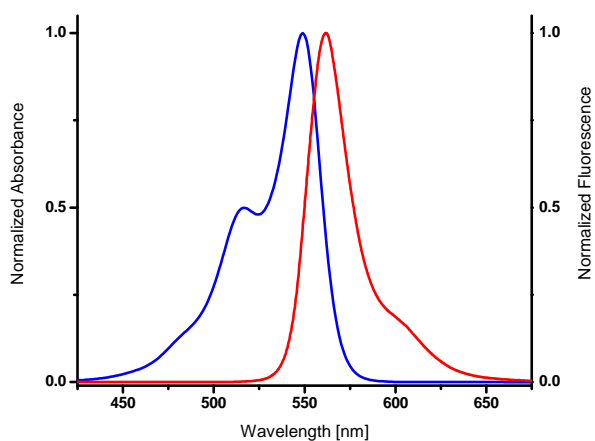
Form: solution in water

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of Cy3