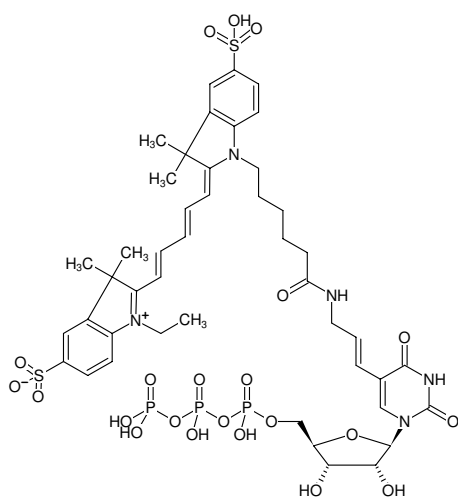




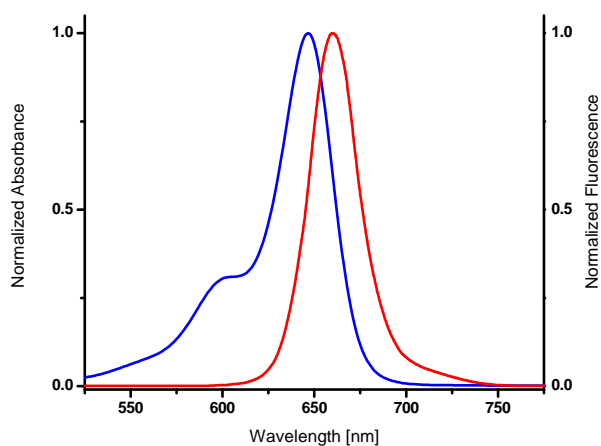
Aminoallyl-UTP-Cy5

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

| Cat. No. | Amount |
|------------|--------------|
| NU-821-CY5 | 30 µl (1 mM) |



Structural formula of Aminoallyl-UTP-Cy5



excitation and emission spectrum of Cy5

For research use only!

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: C₄₅H₅₈N₅O₂₂P₃S₂ (free acid)

Molecular Weight: 1178.01 g/mol (free acid)

Exact Mass: 1177.22 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 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)

Selected References:

Guerra (2006) Analysis of oligonucleotide microarrays by 3' end labeling using fluorescent nucleotides and terminal transferase. *Biotechniques* **41** (1):53.

Hoen et al. (2003) Fluorescent labelling of cRNA for microarray applications. *Nucleic Acids Res.* **31** (5):e20.