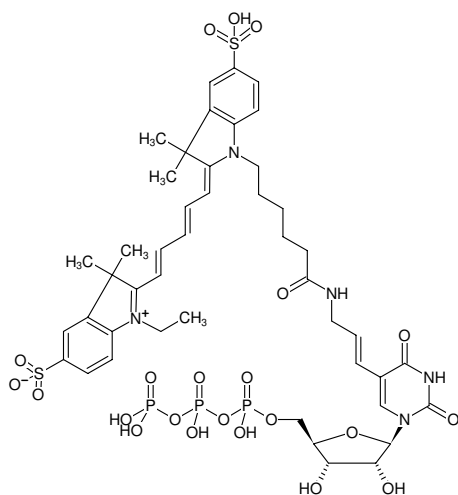




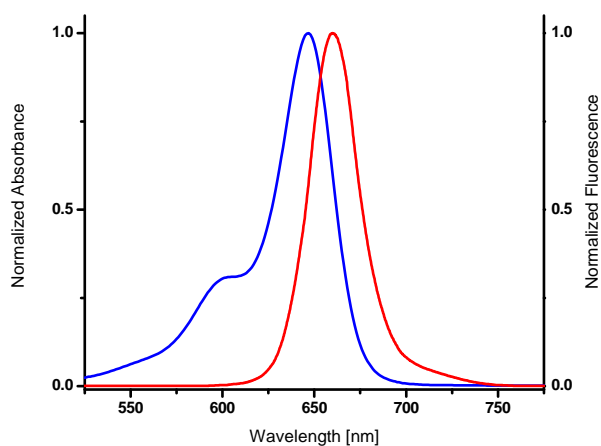
## Aminoallyl-UTP-Cy5

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 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:** C<sub>45</sub>H<sub>58</sub>N<sub>5</sub>O<sub>22</sub>P<sub>3</sub>S<sub>2</sub> (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<sub>exc</sub> 649 nm, λ<sub>em</sub> 670 nm, ε 250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (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.