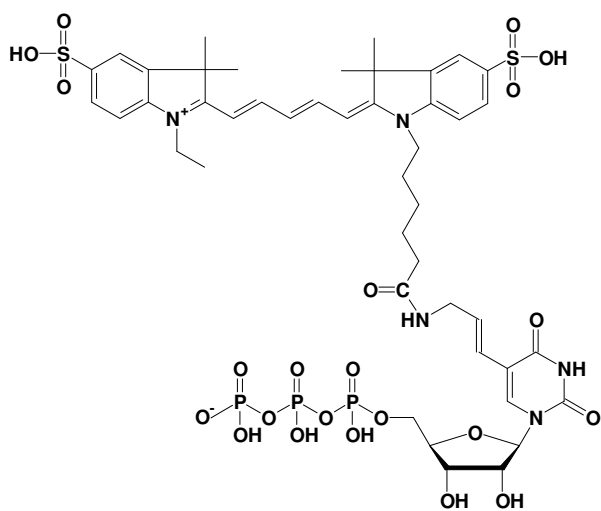


Aminoallyl-UTP - Cy5

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

Cat. No.	Amount
NU-821-CY5	60 μ l (0.5 mM)



structural formula of Aminoallyl-UTP - CY3

Cat. No.: NU-821-CY5

Molecular Formula: C₄₅H₅₈N₅O₂₂P₃S₂ (free acid)

Molecular Weight: 1178,01 (free acid))

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 649 nm; λ_{em} 670 nm; ϵ 2,5 x 10⁵ cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

Selected References:

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

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