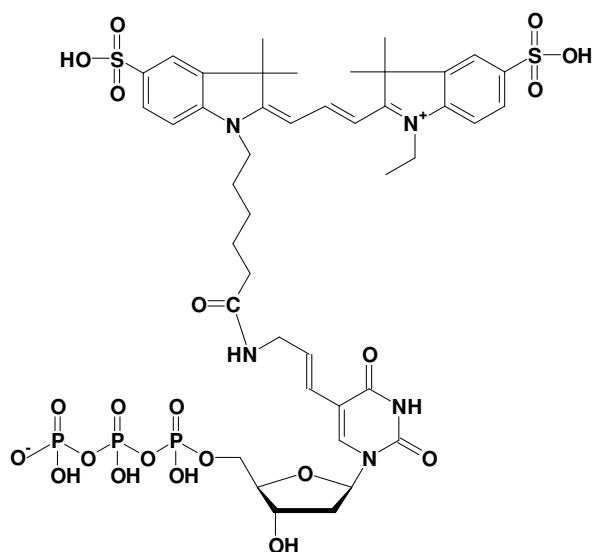


Aminoallyl-dUTP - Cy3

5-(3-Aminoallyl)-2'-deoxy-uridine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-803-CY3	30 μ l (1 mM)



structural formula of Aminoallyl-dUTP - CY3

Cat. No.: NU-803-CY3

Molecular Formula: C₄₃H₅₆N₅O₂₁P₃S₂ (free acid)

Molecular Weight: 1135,97 (free acid)

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

Spectroscopic properties:

λ_{exc} 550 nm; λ_{em} 570 nm; ϵ 1,5 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!