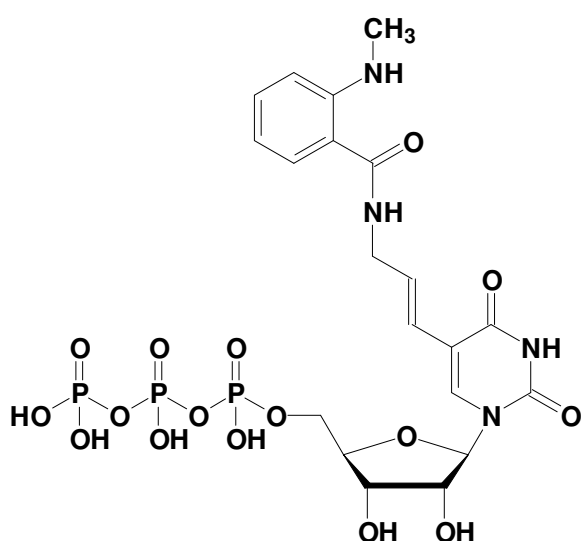


Aminoallyl-UTP - MANT

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

Cat. No.	Amount
NU-821-MNT	100 μ l (5 mM)



structural formula of Aminoallyl-UTP - MANT

Cat. No.: NU-821-MNT

Molecular Formula: C₂₀H₂₇N₄O₁₆P₃ (free acid)

Molecular Weight: 672.37 (free acid)

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

Spectroscopic properties:

λ_{exc} 335 nm; λ_{em} 440 nm; ϵ 2,700 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!