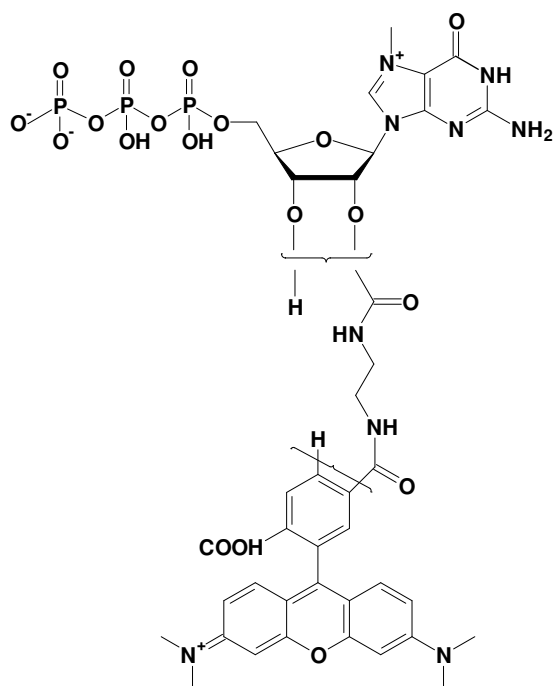


EDA-m⁷GTP - 5/6-TAMRA

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with 5/6-TAMRA, Triethylammonium salt

Cat. No.	Amount
NU-824-TAM	160 μ l (1 mM)



structural formula of EDA-m⁷GTP - 5/6-TAMRA

Cat. No.: NU-824-TAM

Molecular Formula: C₃₉H₄₄N₉O₁₉P₃ (free acid)

Molecular Weight: 1035.74 (free acid)

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

Spectroscopic properties:

λ_{exc} 545 nm; λ_{em} 575 nm; ϵ 90,000 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!