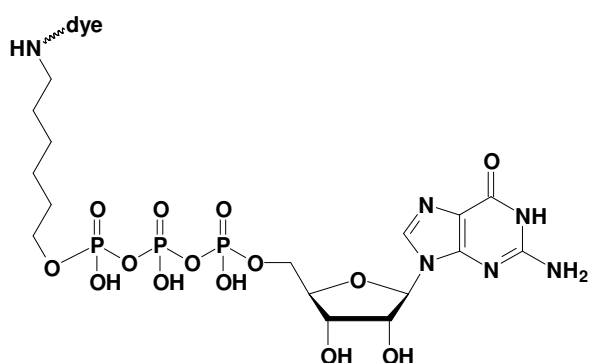


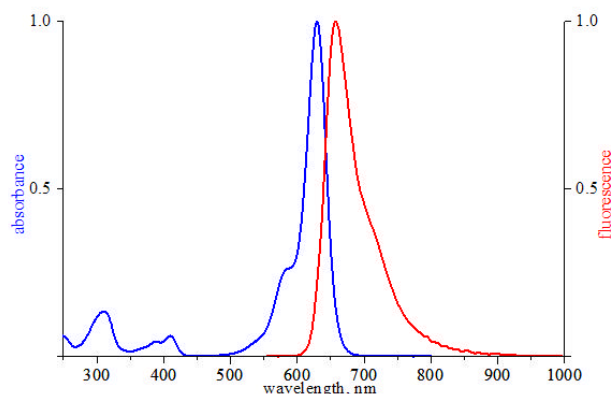
γ -[6-Aminoethyl]-GTP - ATTO 633

γ -[6-Aminoethyl]-guanosine-5'-triphosphate, labeled with ATTO 633, Triethylammonium salt

Cat. No.	Amount
NU-834-633	80 μ l (1 mM)



structural formula of γ -[6-Aminoethyl]-GTP



excitation and emission spectrum of ATTO 633

Cat. No.: NU-834-633

Molecular Formula: C₁₆H₂₉N₆O₁₄P₃ - ATTO 633 (free acid)

Molecular Weight: 1155.35 (free acid)

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

Spectroscopic properties:

λ_{exc} 629 nm; λ_{em} 657 nm; ϵ 130,000 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!