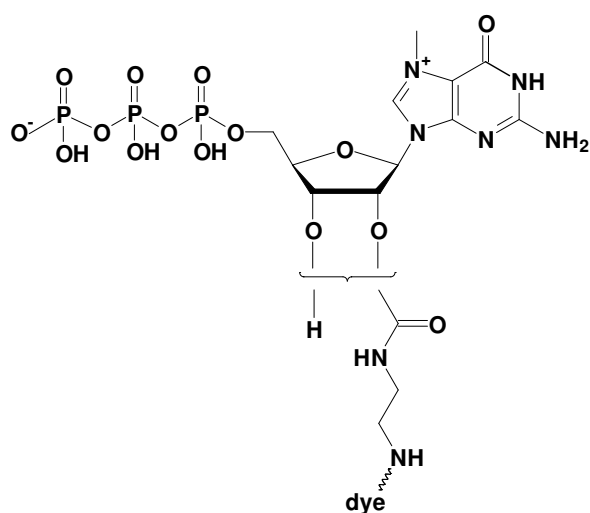


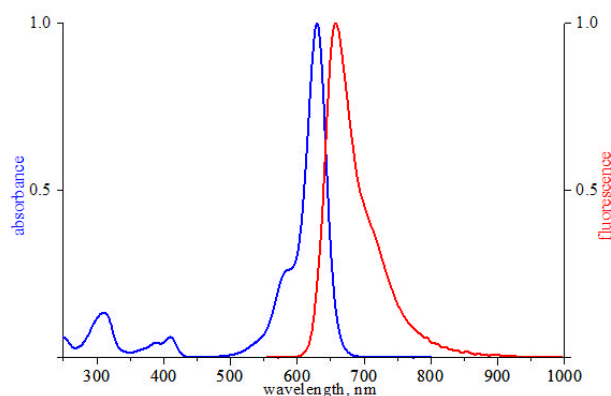
EDA-m⁷GTP - ATTO 633

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with ATTO 633, Triethylammonium salt

Cat. No.	Amount
NU-824-633	10 μ l / 1 mM



structural formula of EDA-m⁷GTP



excitation and emission spectrum of ATTO 633

Cat. No.: NU-824-633

Molecular Formula: C₁₄H₂₄N₇O₁₅P₃ - ATTO 633 (free acid)

Molecular Weight: 1156.30 (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!