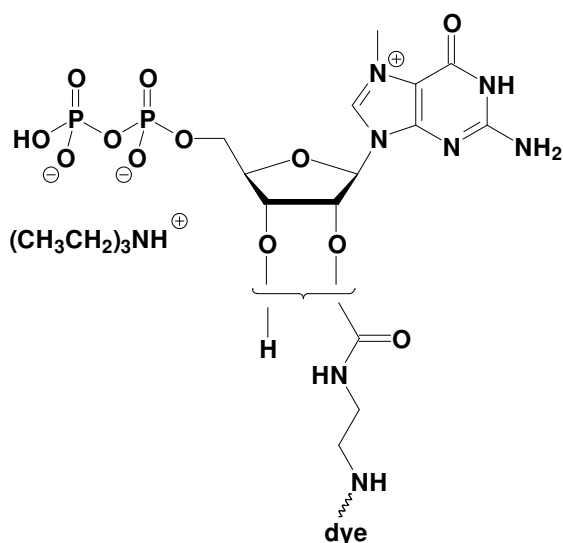


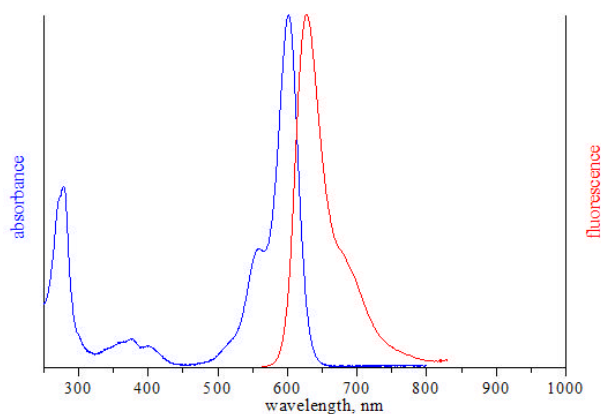
EDA-m⁷GDP - ATTO 594

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 594, Triethylammonium salt

Cat. No.	Amount
NU-827-594	10 μ l / 1 mM



structural formula of EDA-m⁷GDP



excitation and emission spectrum of ATTO 594

Cat. No.: NU-827-594

Molecular Formula: C₁₄H₂₁N₇O₁₂P₂ - ATTO 594 (Anion)

Molecular Weight: 1330.33 (Anion)

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

Spectroscopic properties:

λ_{exc} 601 nm; λ_{em} 627 nm; ϵ 120,000 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.