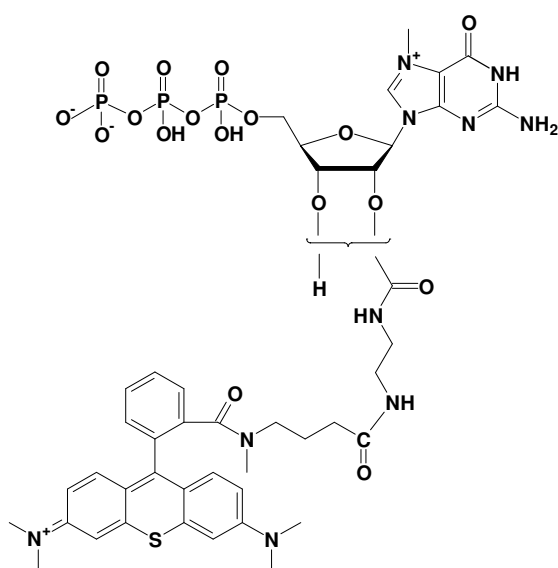


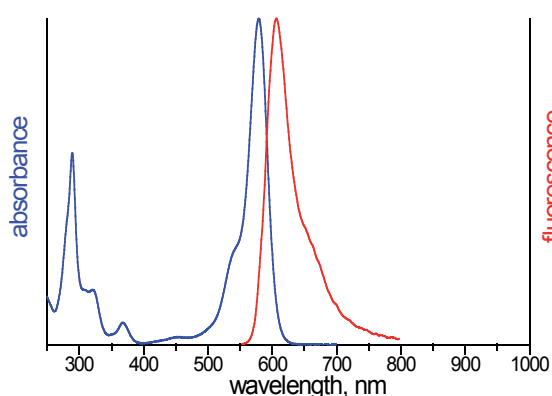
EDA-m⁷GTP - ATTO Thio12

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-824-THIO12	40 μ l (1 mM)



structural formula of EDA-m⁷GTP - ATTO Thio12



excitation and emission spectrum of ATTO Thio12

Cat. No.: NU-824-THIO12

Molecular Formula: C₄₃H₅₃N₁₀O₁₇P₃S (free acid)

Molecular Weight: 1106,93 (free acid)

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

Spectroscopic properties:

λ_{exc} 579 nm; λ_{em} 609 nm; ϵ 1,1 x 10⁵ 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!