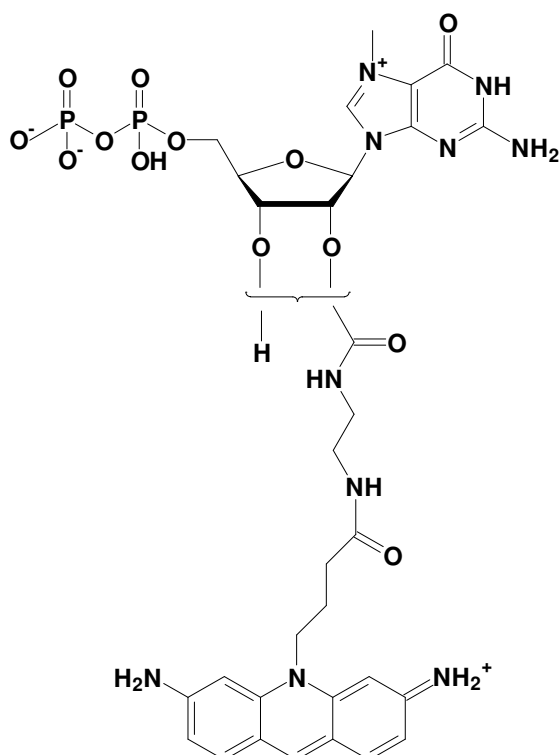


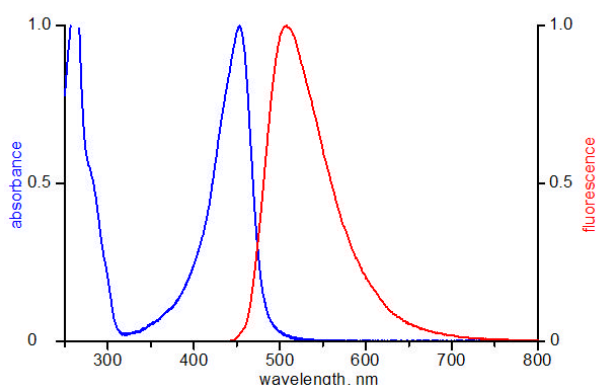
EDA-m⁷GDP - ATTO 465

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 465

Cat. No.	Amount
NU-827-465	120 μ l (1 mM)



structural formula of EDA-m⁷GDP - ATTO 465



excitation and emission spectrum of ATTO 465

Cat. No.: NU-827465

Molecular Formula: C₃₁H₃₈N₁₀O₁₃P₂ (free acid)

Molecular Weight: 820.64 (free acid)

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

Spectroscopic properties:

λ_{exc} 453 nm; λ_{em} 508 nm; ϵ 75,000 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!