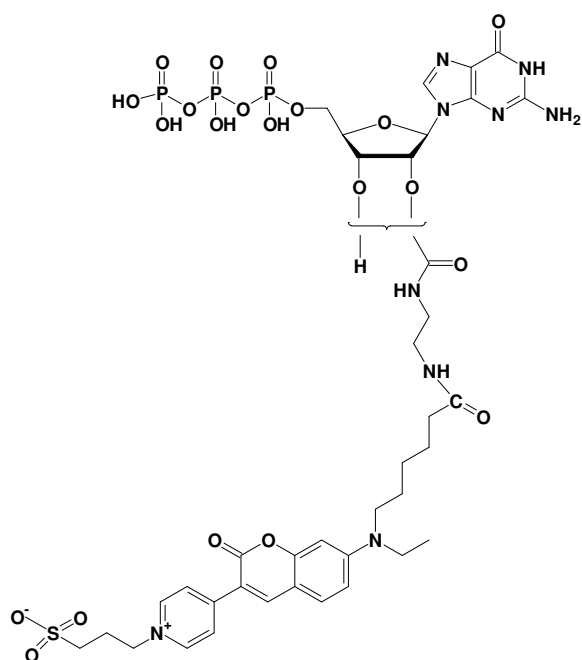


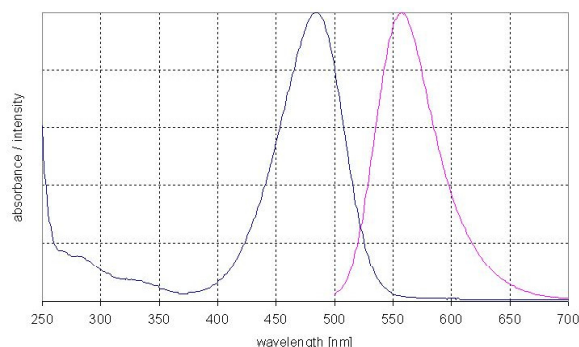
EDA-GTP - DY 485XL

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, labeled with DY 485XL,
 Triethylammonium salt

Cat. No.	Amount
NU-820-485XL	40 μ l (1 mM)



structural formula of EDA-GTP - DY 485XL



excitation and emission spectrum of DY 485XL

Cat. No.: NU-820-485XL

Molecular Formula: C₃₈H₅₀N₉O₂₁P₃S (free acid)

Molecular Weight: 1093,84 (free acid)

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

Spectroscopic properties:

λ_{exc} 485 nm; λ_{em} 560 nm; ϵ 50,000 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!