



EDA-m⁷GTP - DY-485XL

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-824-485XL	40 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula: C₃₉H₅₂N₉O₂₁P₃S (free acid)

Molecular Weight: 1107.87 g/mol (free acid)

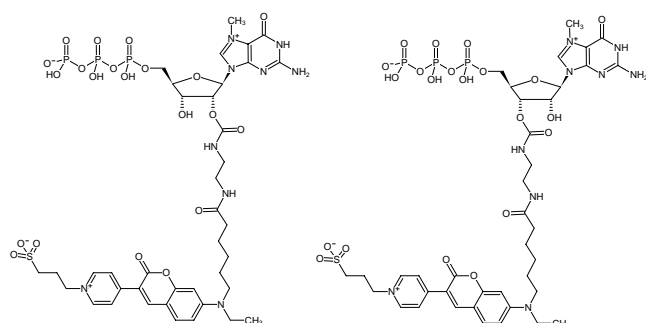
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution, pH 7.5 ± 0.5

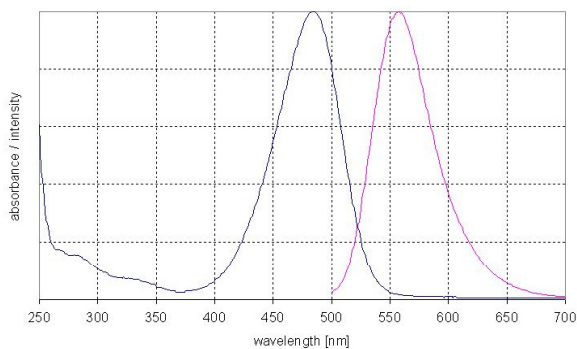
Concentration: 1 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 485 nm; λ_{em} 560 nm; ε 50.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GTP - DY-485XL



excitation and emission spectrum of DY 485XL