



EDA-m⁷GDP-ATTO-488

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 488, Triethylammonium salt

Cat. No.	Amount
NU-827-488	80 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.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₉H₄₄N₁₀O₂₁P₂S₂ (free acid)

Molecular Weight: 1114.16 g/mol (free acid)

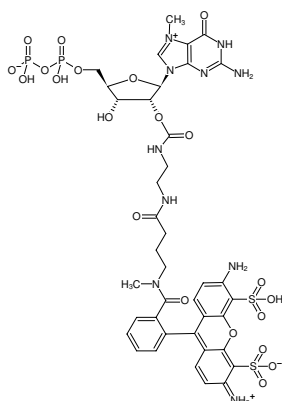
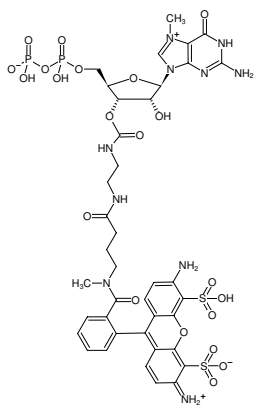
Purity: ≥ 95 % (HPLC)

Form: yellow solution in water

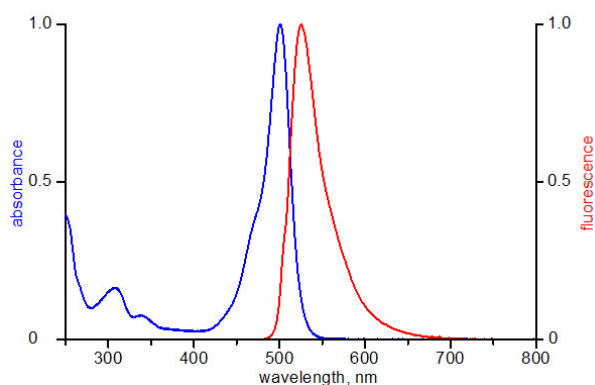
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 500 nm, λ_{em} 520 nm, ε 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GDP-ATTO-488



excitation and emission spectrum of ATTO 488