



EDA-GppNHp (EDA-GMPPNP)-ATTO-488

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO 488, Triethylammonium salt

Cat. No.	Amount
NU-860-488	50 μ l (1 mM)

For research use only!

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₃₈H₄₄N₁₁O₂₃P₃S₂ (free acid)

Molecular Weight: 1179.87 g/mol (free acid)

Exact Mass: 1179.13 g/mol (free acid)

Purity: \geq 95 % (HPLC)

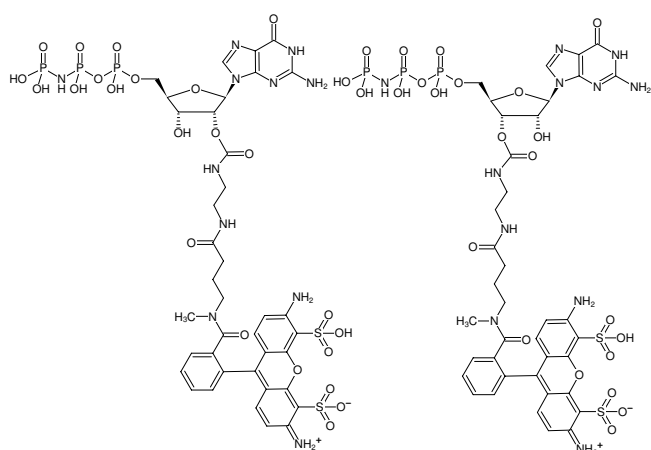
Form: solution in water

Color: yellow

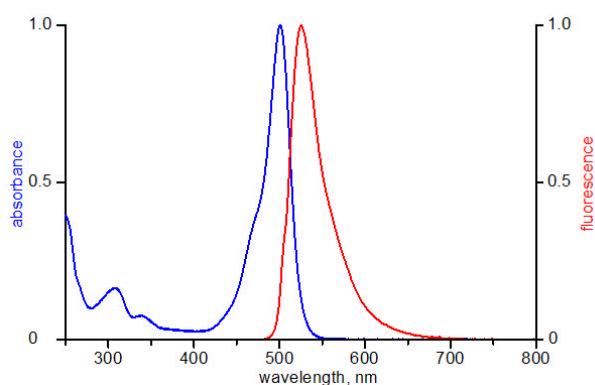
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 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-GppNHp (EDA-GMPPNP)-ATTO-488



excitation and emission spectrum of ATTO 488

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.

Selected References:

Lee *et al.* (2017) Mechanism of SOS PR-domain autoinhibition revealed by single-molecule assays on native protein from lysate. *Nat. Commun.* **8**:15061.

Christensen *et al.* (2016) One-way membrane trafficking of SOS in receptor-triggered Ras activation. *Nat. Struct. Mol. Biol.* **23** (9):838.

Lin *et al.* (2014) H-Ras forms dimers on membrane surfaces via a protein-protein interface. *Proc. Natl. Acad. Sci. USA* **111** (8):2996.