



EDA-GppNHp (EDA-GMPPNP)-ATTO-647N

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-860-647N	50 μ l (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: 6 months after date of delivery

Molecular Formula: C₅₅H₇₂N₁₁O₁₆P₃ (free acid)

Molecular Weight: 1236.15 g/mol (free acid)

Exact Mass: 1235.44 g/mol (free acid)

Purity: \geq 95 % (HPLC)

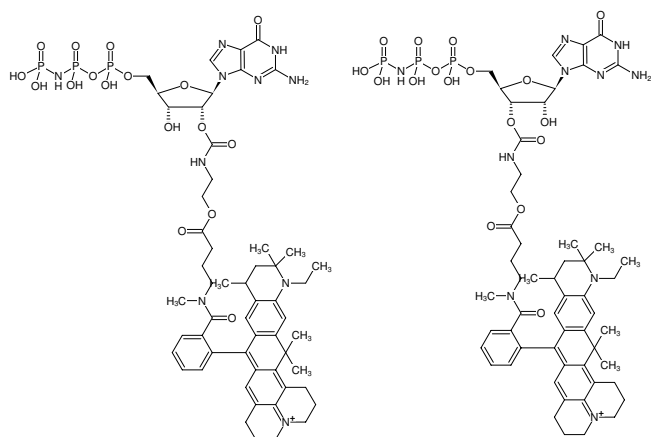
Form: blue solution in water

Concentration: 1.0 mM - 1.1 mM

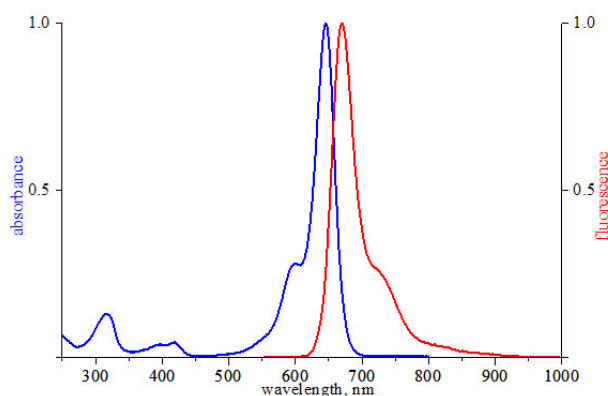
pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 646 nm, λ_{em} 664 nm, ϵ 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

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.



Structural formula of EDA-GppNHp (EDA-GMPPNP)-ATTO-647N



excitation and emission spectrum of ATTO 647N