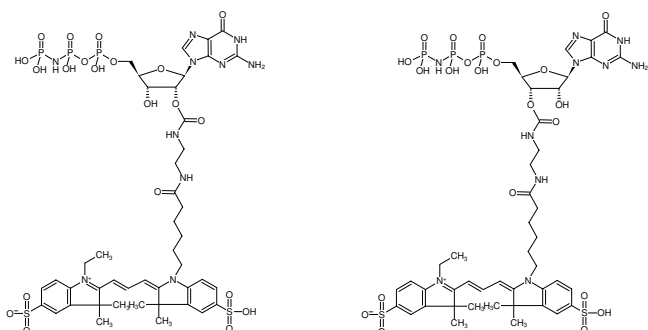




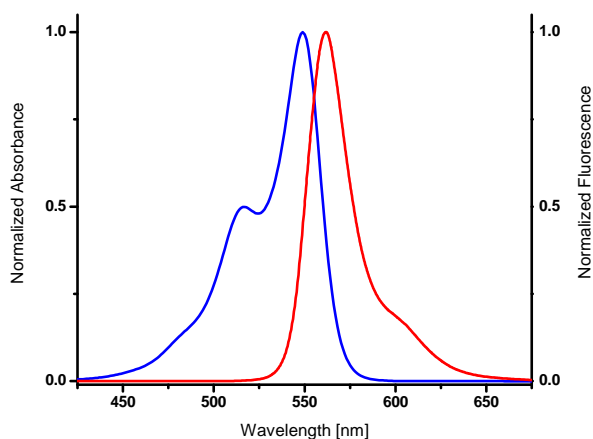
EDA-GppNHp (EDA-GMPPNP)-Cy3

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-[(β,γ)-imido] triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-860-CY3	50 µl (1 mM)



Structural formula of EDA-GppNHp (EDA-GMPPNP)-Cy3



excitation and emission spectrum of Cy3

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₄₄H₅₉N₁₀O₂₁P₃S₂ (free acid)

Molecular Weight: 1221.05 g/mol (free acid)

Exact Mass: 1220.25 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 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.