



EDA-m⁷GDP-Cy5

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-827-CY5	160 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: C₄₇H₆₁N₉O₁₉P₂S₂ (free acid)

Molecular Weight: 1182.11 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

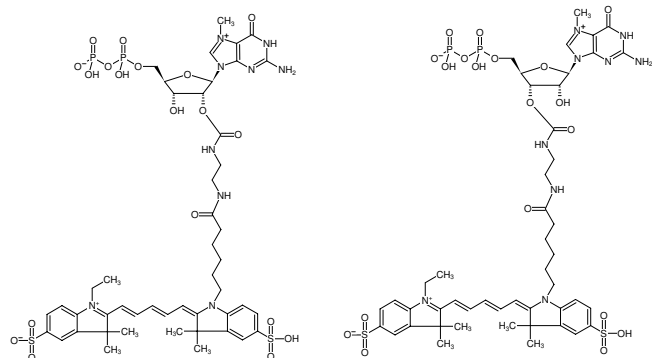
Form: solution in water

Color: blue

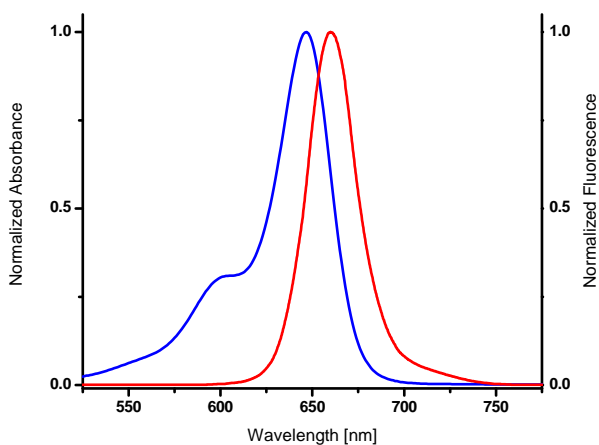
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GDP-Cy5



excitation and emission spectrum of Cy5