



γ -(6-Aminoethyl)-GTP-DY-480XL

γ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-834-480XL	80 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: 12 months after date of delivery

Molecular Formula: C₄₂H₅₇N₈O₂₀P₃S (free acid)

Molecular Weight: 1118.93 g/mol (free acid)

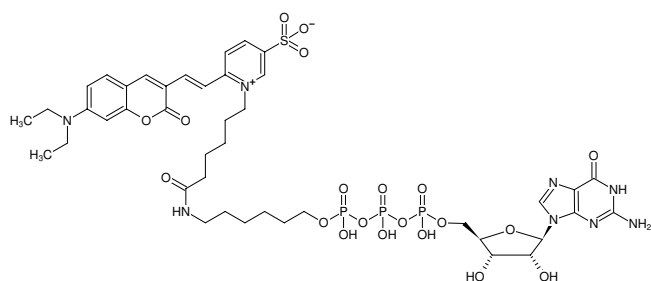
Purity: \geq 95 % (HPLC)

Form: clear aqueous solution

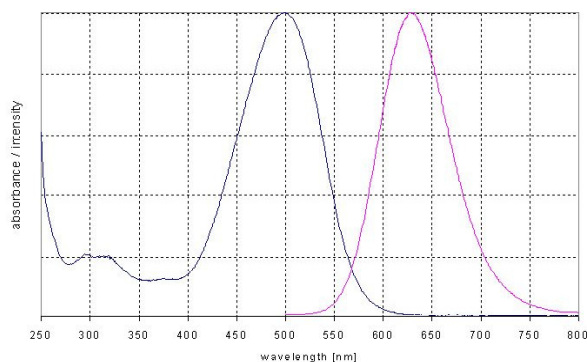
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 500 nm, λ_{em} 630 nm, ϵ 50.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of γ -(6-Aminoethyl)-GTP-DY-480XL



excitation and emission spectrum of DY 480XL