



EDA-GTPyS-DY-480XL

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-1610-480XL	50 µl (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: C₃₉H₅₀N₉O₂₀P₃S₂ (free acid)

Molecular Weight: 1121.91 g/mol (free acid)

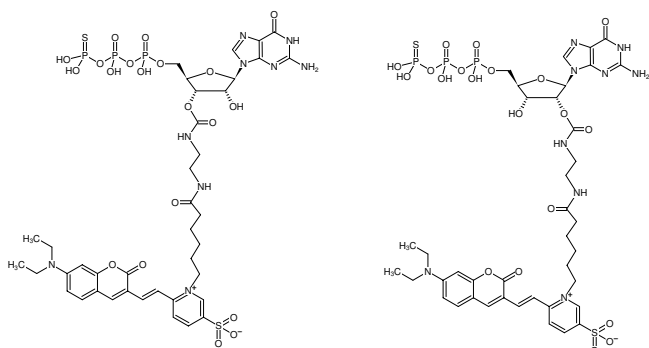
Purity: ≥ 90 % (HPLC), contains approx. 5 % EDA-GDP-dye

Form: solution in water

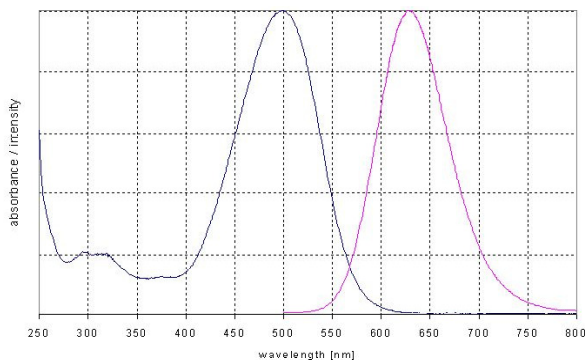
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

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



Structural formula of EDA-GTPyS-DY-480XL



excitation and emission spectrum of DY 480XL