



EDA-GTPyS-6-JOE

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with 6-JOE, Triethylammonium salt

Cat. No.	Amount
NU-1610-JOE	50 µl (1 mM)

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: C₃₆H₃₄Cl₂N₇O₂₂P₃S (free acid)

Molecular Weight: 1111.01 g/mol (free acid)

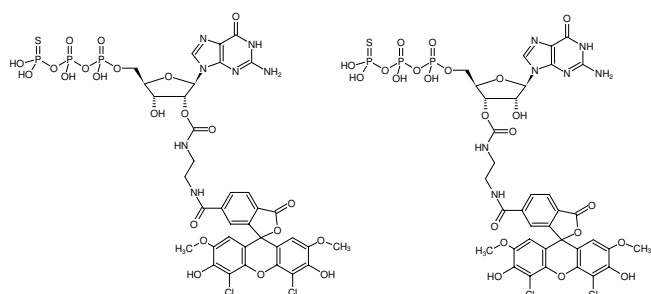
Purity: ≥ 90 % (HPLC)

Form: solution in water

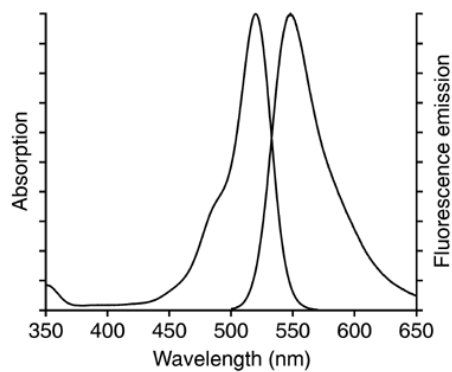
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 520 nm, λ_{em} 548 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-6-JOE



excitation and emission spectrum of 6-JOE