



EDA-GTPyS-DYQ-661

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-1610-661Q	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: C₅₂H₆₃N₉O₂₂P₃S₃ (Anion)

Molecular Weight: 1355.22 g/mol (Anion)

Purity: ≥ 90 % (HPLC)

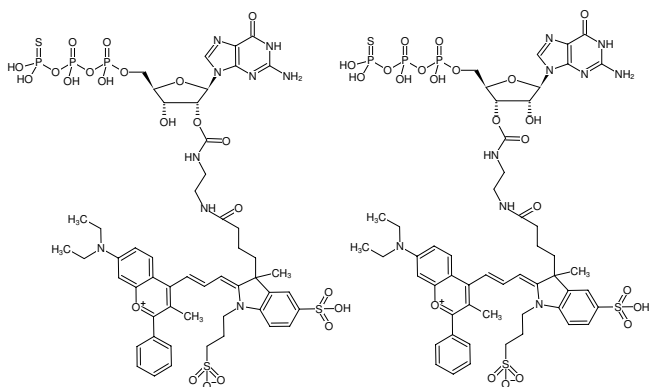
Form: solution in water

Concentration: 1.0 mM - 1.1 mM

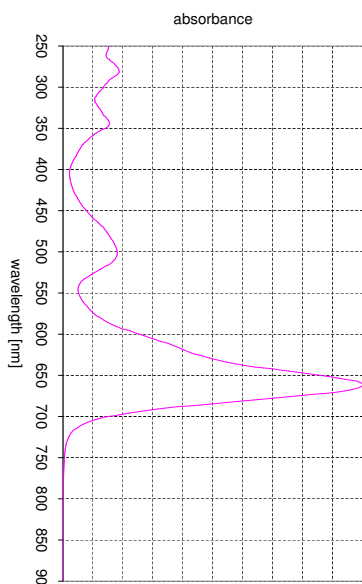
pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 662 nm, ε 104.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 662 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)



Structural formula of EDA-GTPyS-DYQ-661



absorption spectrum of DYQ 661

Selected References:

[1] Tan *et al.* (2010) Direct detection of Delta9-tetrahydrocannabinol in saliva using a novel homogeneous competitive immunoassay with fluorescence quenching. *Anal. Chim. Acta.* **658** (2):187.