



EDA-AppNHp (EDA-AMPPNP)-DYQ-661

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-810-661Q	40 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: C₅₂H₆₄N₁₀O₂₁P₃S₂ (Anion)

Molecular Weight: 1322.17 g/mol (Anion)

Purity: ≥ 95 % (HPLC)

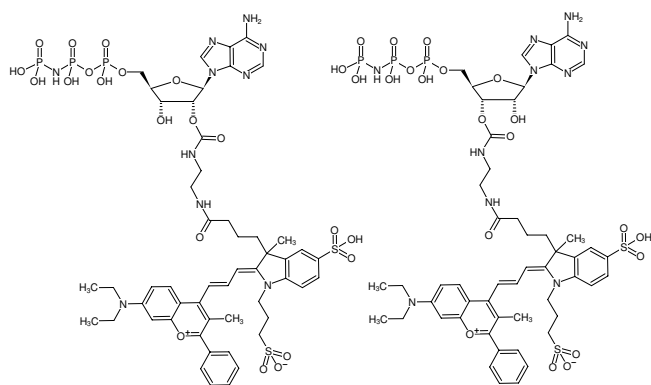
Form: solution in water

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)

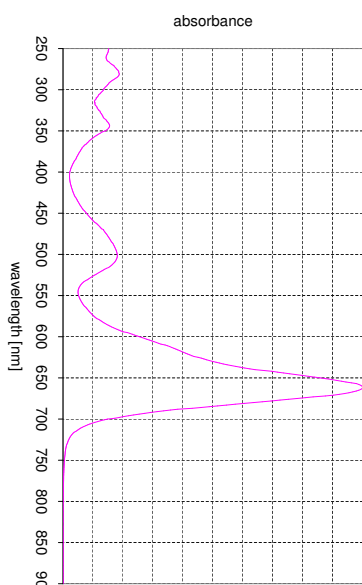


Structural formula of EDA-AppNHp (EDA-AMPPNP)-DYQ-661

Please note: DYQ661 is not a dark quencher, λ_{em} 690 nm^[1]

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.



absorption spectrum of DYQ 661