



EDA-AppNHp (EDA-AMPPNP)-DY-751

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-[(β,γ)-imido] triphosphate, labeled with DY 751, Triethylammonium salt

Cat. No.	Amount
NU-810-751	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: C₅₅H₇₃N₁₀O₂₁P₃S₂ (free acid)

Molecular Weight: 1366.36 g/mol (free acid)

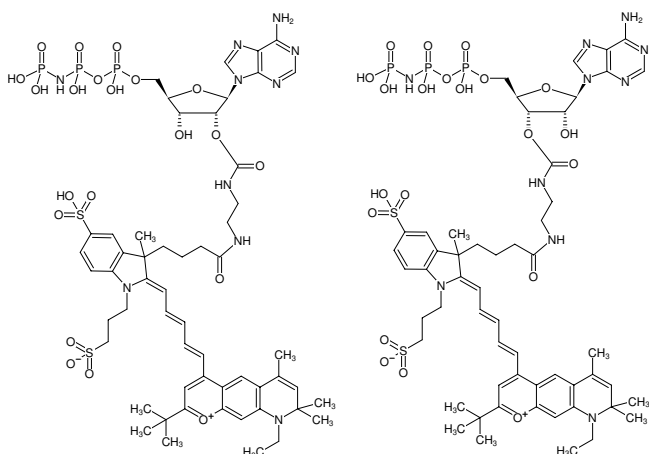
Purity: ≥ 95 % (HPLC)

Form: solution in water

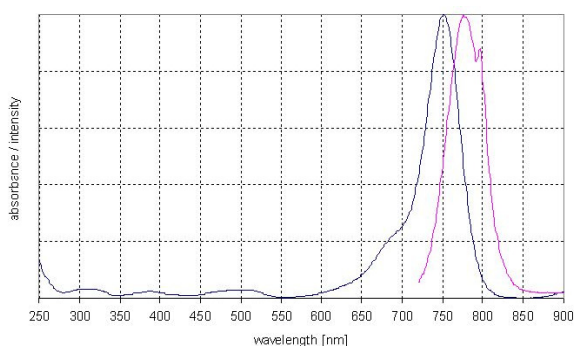
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 751 nm, λ_{em} 779 nm, ε 270.0 L mmol⁻¹ cm⁻¹ (Ethanol)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-DY-751



excitation and emission spectrum of DY 751