



EDA-AppNHp (EDA-AMPPNP)-6-FAM

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with 6 FAM, Triethylammonium salt

Cat. No.	Amount
NU-810-6FM	120 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: C₃₄H₃₃N₈O₁₉P₃ (free acid)

Molecular Weight: 950.59 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

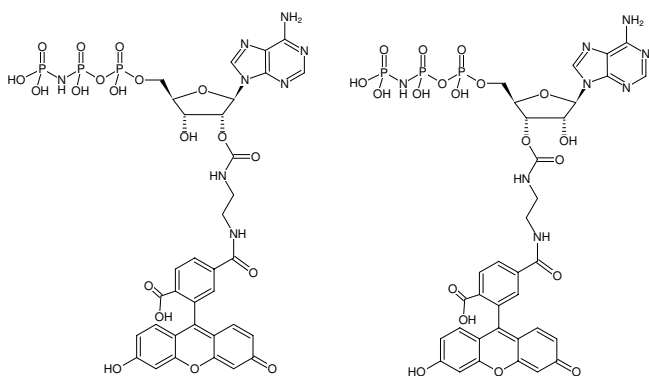
Form: solution in water

Color: yellow

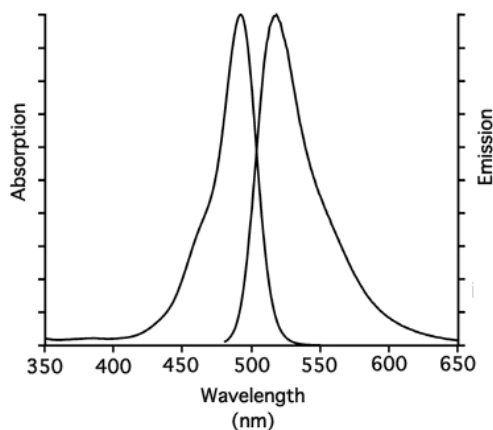
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-6-FAM



excitation and emission spectrum of 6-FAM