



EDA-AppNHp (EDA-AMPPNP)-ATTO-465

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-810-465	80 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤ 2 ml).

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₃₀H₃₈N₁₁O₁₄P₃ (free acid)

Molecular Weight: 869.61 g/mol (free acid)

Exact Mass: 869.18 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

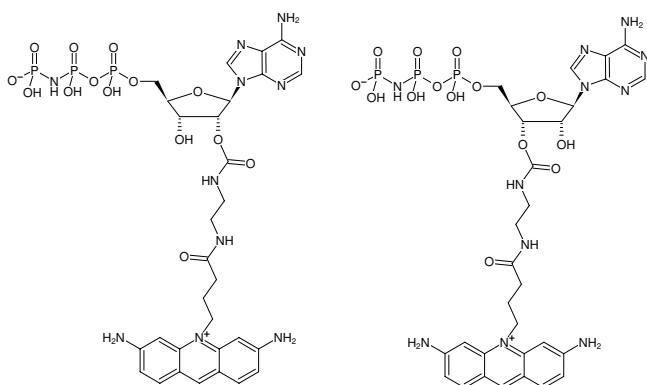
Form: solution in water

Color: yellow

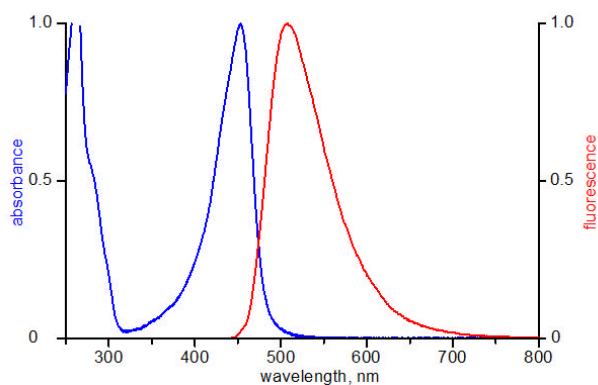
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 453 nm, λ_{em} 506 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-465



excitation and emission spectrum of ATTO 465