



EDA-AppNHp (EDA-AMPPNP)-ATTO-425

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO 425, Triethylammonium salt

Cat. No.	Amount
NU-810-425	80 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: C₃₅H₄₈N₉O₁₈P₃ (free acid)

Molecular Weight: 975.73 g/mol (free acid)

Exact Mass: 975.23 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

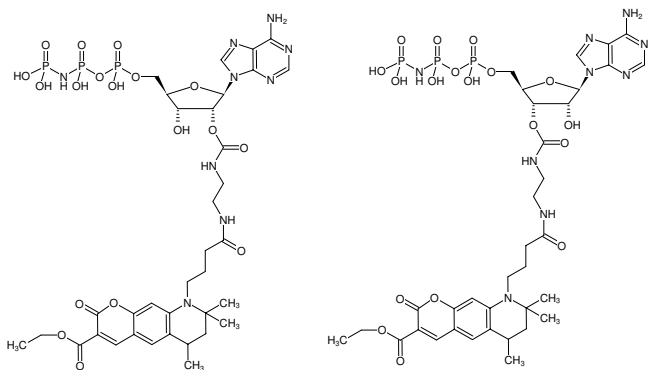
Form: solution in water

Color: green-yellow

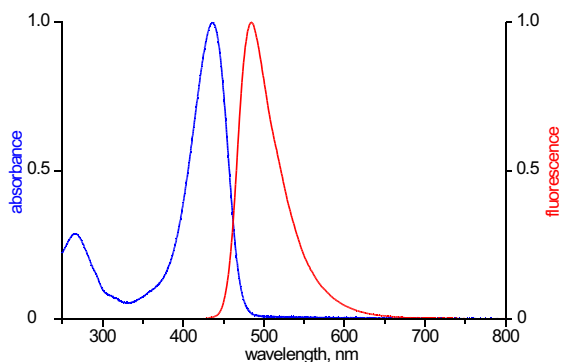
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 439 nm, λ_{em} 485 nm, ε 45.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-425



excitation and emission spectrum of ATTO 425