



EDA-AppNHp (EDA-AMPPNP)-DY-485XL

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-810-485XL	40 μ l (1 mM)

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at $-20\text{ }^{\circ}\text{C}$

Shelf Life: 6 months after date of delivery

Molecular Formula: $\text{C}_{38}\text{H}_{51}\text{N}_{10}\text{O}_{19}\text{P}_3\text{S}$ (free acid)

Molecular Weight: 1076.85 g/mol (free acid)

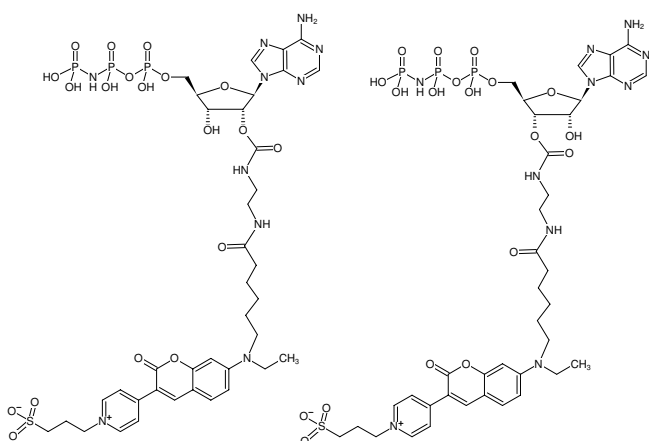
Purity: $\geq 95\%$ (HPLC)

Form: solution in water

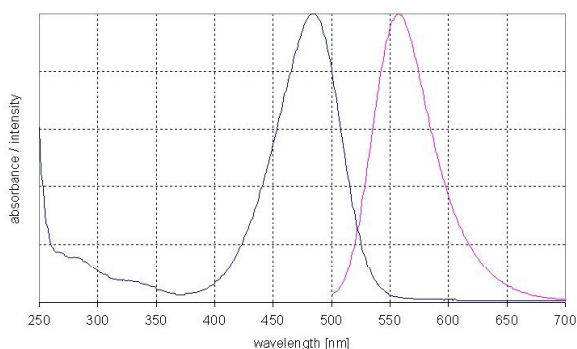
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 485 nm, λ_{em} 560 nm, ϵ 50.0 L $\text{mmol}^{-1}\text{ cm}^{-1}$ (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-DY-485XL



excitation and emission spectrum of DY 485XL