



EDA-AppNHp (EDA-AMPPNP)-6-ROX

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

| Cat. No. | Amount |
|------------|---------------|
| NU-810-ROX | 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: C₄₆H₅₁N₁₀O₁₇P₃ (free acid)

Molecular Weight: 1108.88 g/mol (free acid)

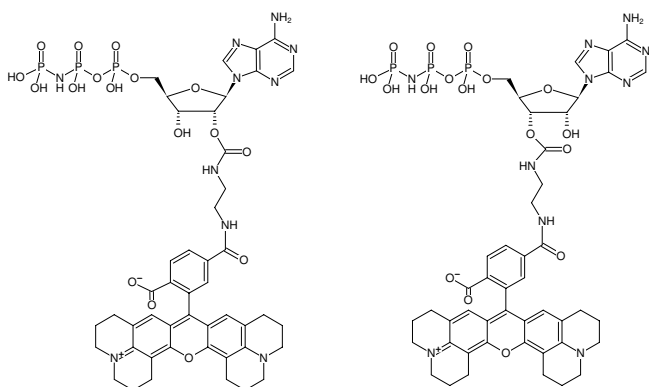
Purity: ≥ 95 % (HPLC)

Form: solution in water

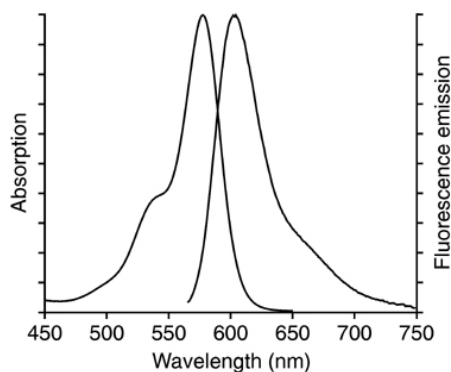
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 575 nm, λ_{em} 600 nm, ε 82.0 L mol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



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



excitation and emission spectrum of 6-ROX