



EDA-AppNHp (EDA-AMPPNP)-ATTO-594

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO 594, Triethylammonium salt

| Cat. No. | Amount |
|------------|--------------|
| NU-810-594 | 40 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: C₅₄H₆₈N₁₁O₂₂P₃S₂ (free acid)

Molecular Weight: 1380.23 g/mol (free acid)

Exact Mass: 1379.32 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

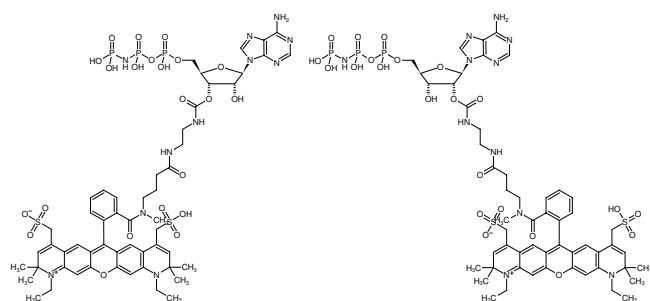
Form: solution in water

Color: red-violet

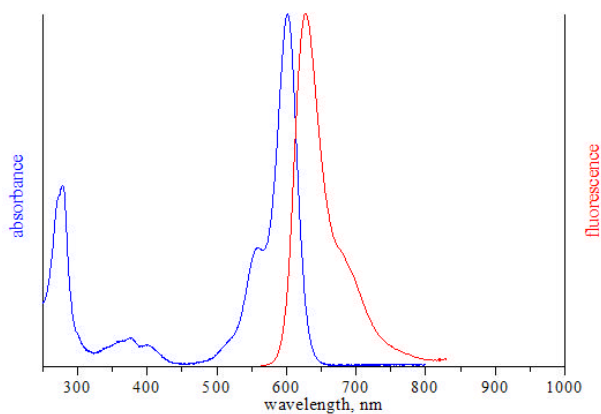
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 626 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-594



excitation and emission spectrum of ATTO 594