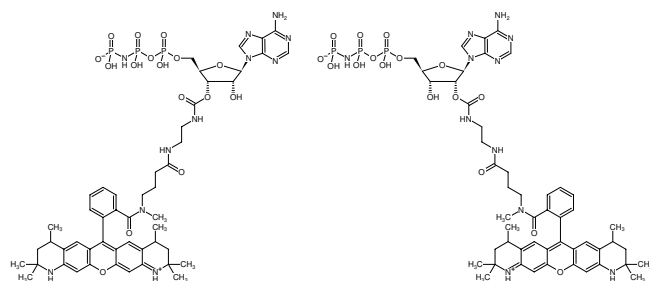


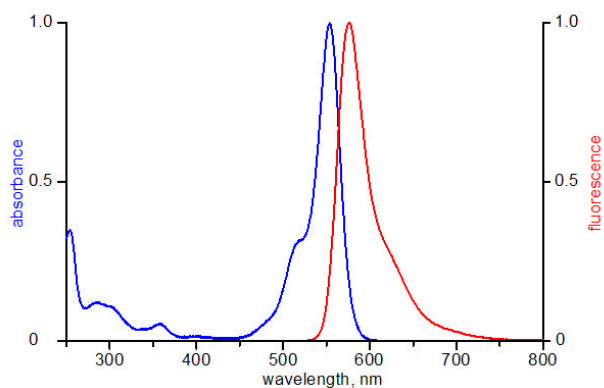

EDA-AppNHp (EDA-AMPPNP)-ATTO-550

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido]triphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-810-550	40 µl (1 mM)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-550



excitation and emission spectrum of ATTO 550

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: C₅₀H₆₄N₁₁O₁₆P₃ (free acid)

Molecular Weight: 1168.04 g/mol (free acid)

Exact Mass: 1167.37 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} 554 nm, λ_{em} 576 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)