



EDA-AppNHp (EDA-AMPPNP) - ATTO-700

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO 700, Triethylammonium salt

Cat. No.	Amount
NU-810-700	40 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 6 months after date of delivery

Molecular Formula: C₁₃H₂₃N₈O₁₃P₃ - ATTO 700 (free acid)

Molecular Weight: 1113.29 g/mol (free acid)

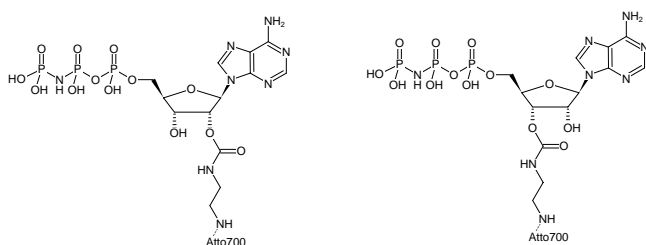
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

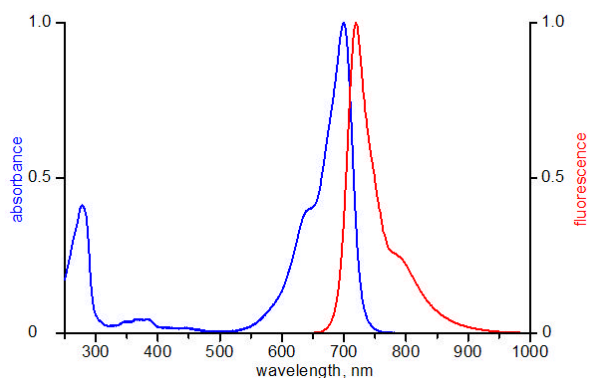
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 700 nm, λ_{em} 716 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-AppNHp (EDA-AMPPNP) - ATTO-700



excitation and emission spectrum of ATTO 700