



γ -(6-Aminoethyl)-AppCp-ATTO-488

gamma-(6-Aminoethyl)-AMPPCP-ATTO-488

γ -(6-Aminoethyl)-adenosine-5'-[(beta,gamma)-methylene]triphosphate, labeled with ATTO-488, Triethylammonium salt

Cat. No.	Amount
NU-279-488	60 μ l (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: 12 months after date of delivery

Molecular Formula: C₄₂H₅₂N₈O₂₁P₃S₂ (free acid)

Molecular Weight: 1175.96 g/mol (free acid)

Exact Mass: 1175.19 g/mol (free acid)

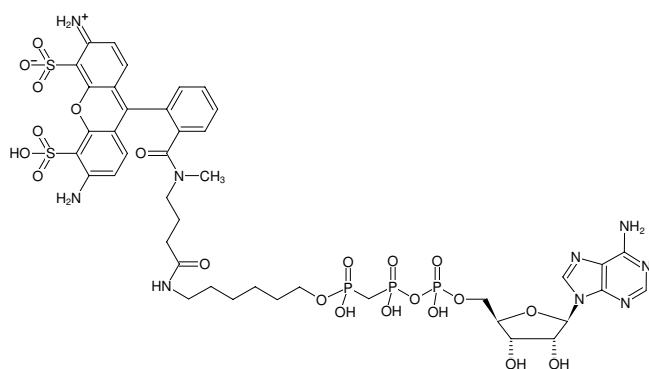
Purity: \geq 95 % (HPLC)

Form: yellow solution in water

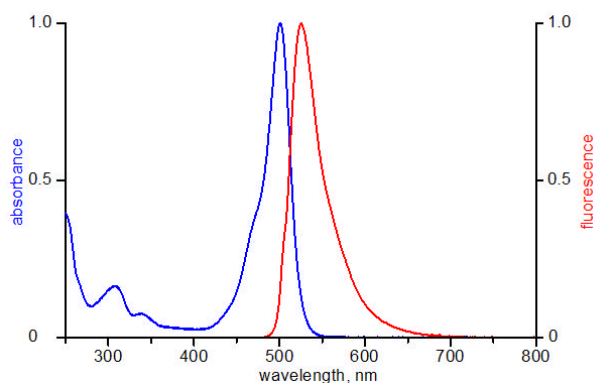
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 500 nm, λ_{em} 520 nm, ϵ 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of γ -(6-Aminoethyl)-AppCp-ATTO-488



excitation and emission spectrum of ATTO 488