

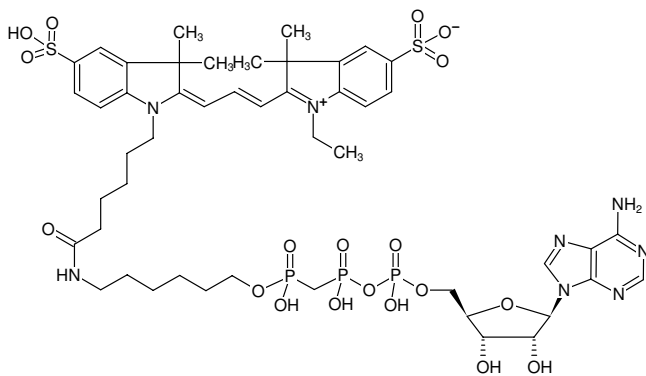


γ -(6-Aminoethyl)-AppCp-Cy3

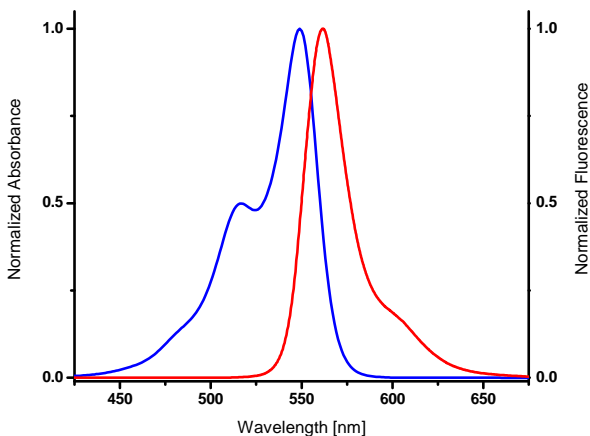
gamma-(6-Aminoethyl)-AMPPCP-Cy3

γ -(6-Aminoethyl)-adenosine-5'-[(beta,gamma)-methylene]triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-279-CY3	60 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-AppCp-Cy3



excitation and emission spectrum of Cy3

For general laboratory use.

Shipping: shipped on gel packs

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: C₄₈H₆₇N₈O₁₉P₃S₂ (free acid)

Molecular Weight: 1217.14 g/mol (free acid)

Exact Mass: 1216.32 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ϵ 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)