



α -[(6-Aminoethyl)imido]-AMP-Cy3

α -[(6-Aminoethyl)imido]-AMP, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-894-CY3	50 μ l (5 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume \leq 2 ml).

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: C₄₇H₆₄N₉O₁₃PS₂ (free acid)

Molecular Weight: 1058.17 g/mol (free acid)

Exact Mass: 1057.38 g/mol (free acid)

Purity: \geq 95 % (HPLC)

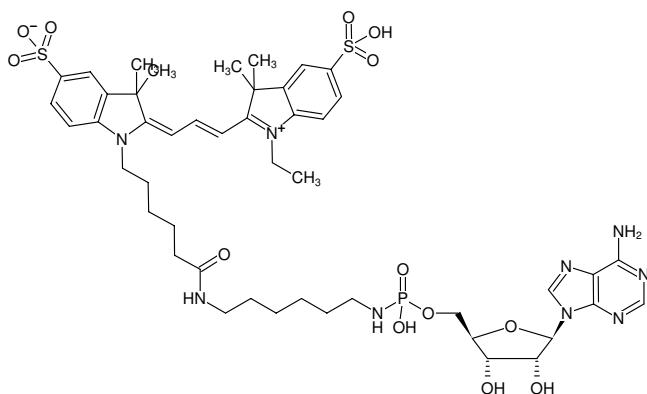
Form: solution in water

Color: pink

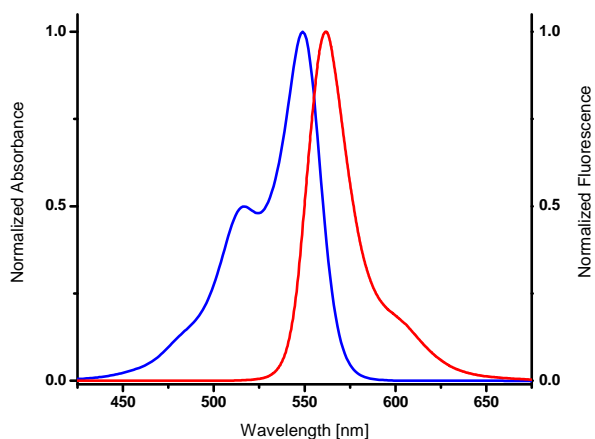
Concentration: 5.0 mM - 5.5 mM

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)



Structural formula of α -[(6-Aminoethyl)imido]-AMP-Cy3



excitation and emission spectrum of Cy3