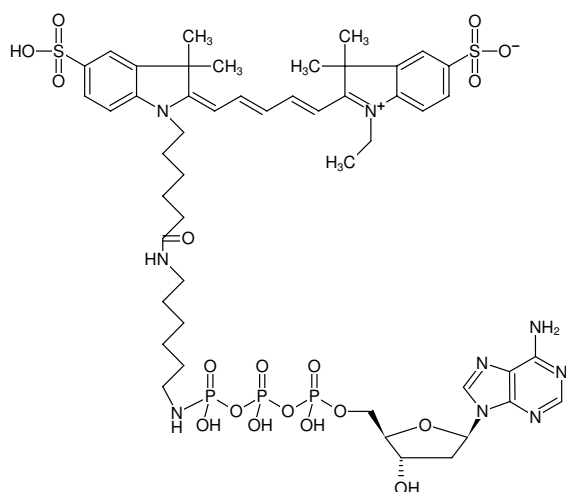




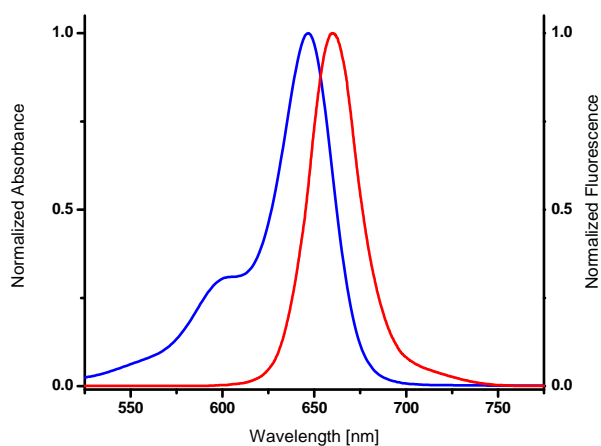
γ -[[(6-Aminoethyl)-imido]-dATP-Cy5

γ -[[(6-Aminoethyl)-imido]-2'-deoxyadenosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-861-CY5	120 μ l (1 mM)



Structural formula of γ -[[6-Aminoethyl)-imido]-dATP-Cy5



excitation and emission spectrum of Cy5

For research use only!

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: C₄₉H₆₈N₉O₁₈P₃S₂ (free acid)

Molecular Weight: 1228.17 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ϵ 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.