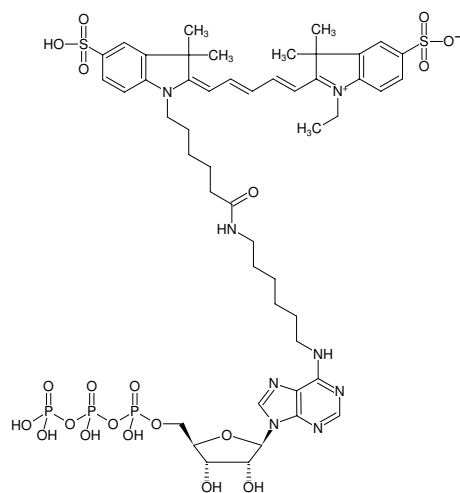




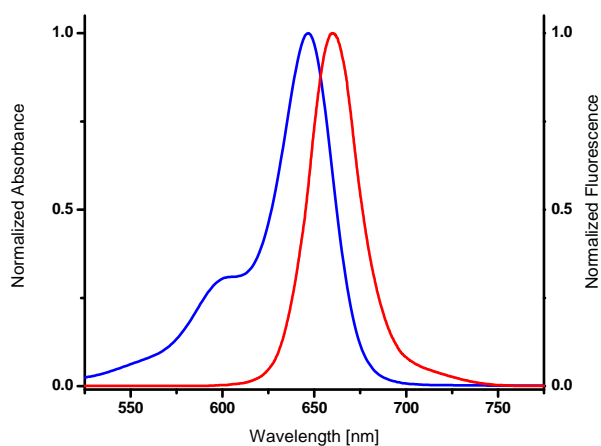
N⁶-(6-Aminoethyl)-ATP-Cy5

N⁶-(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-805-CY5	160 µl (1 mM)



Structural formula of N⁶-(6-Aminoethyl)-ATP-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: C₄₉H₆₇N₈O₂₀P₃S₂ (free acid)

Molecular Weight: 1245.15 g/mol (free acid)

Exact Mass: 1244.31 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 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)

Selected References:

Evans *et al.* (2015) Investigating Apoptozole as a Chemical Probe for HSP70 Inhibition. *PLoS One*. **10** (10):e0140006.