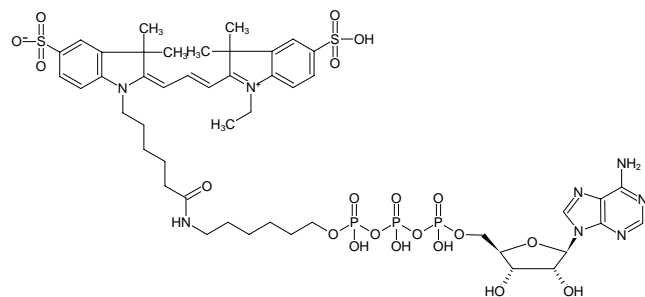




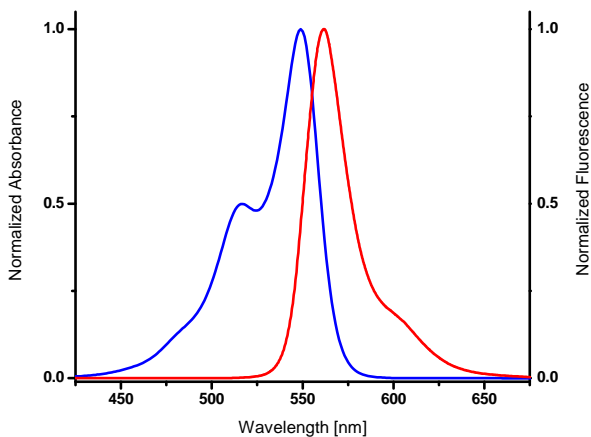
γ-(6-Aminoethyl)-ATP - Cy3

γ-(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-833-CY3	120 µl (1 mM)



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



excitation and emission spectrum of Cy3

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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula: C₄₇H₆₅N₈O₂₀P₃S₂ (free acid)

Molecular Weight: 1219.11 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: pink solution in water, pH 7.5 ±0.5

Concentration: 1 mM - 1.1 mM

pH: 7.5 ±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)

Selected References:

Schax *et al.* (2014) Microarray-based screening of heat shock protein inhibitors. *J. Biotechnol.* **180**:1.