



EDA-ATP-Cy5

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-808-CY5	160 µl (1 mM)

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: C₄₆H₆₀N₉O₂₁P₃S₂ (free acid)

Molecular Weight: 1232.07 g/mol (free acid)

Exact Mass: 1231.26 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)

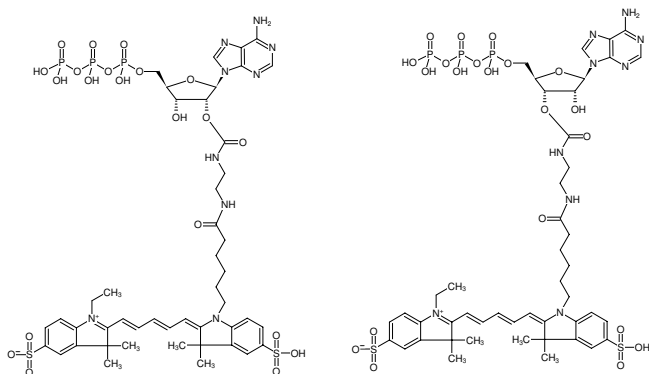
Applications:

Single molecule studies^[1]

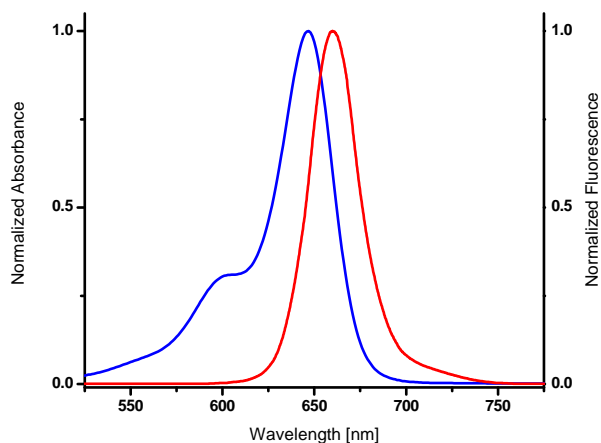
Fluorescence microscopy^[1]

Selected References:

[1] Oiwa *et al.* (2000) Comparative Single-Molecule and Ensemble Myosin Enzymology: Sulfoindocyanine ATP and ADP Derivatives. *Biophys. J.* **78**:3048.



Structural formula of EDA-ATP-Cy5



excitation and emission spectrum of Cy5