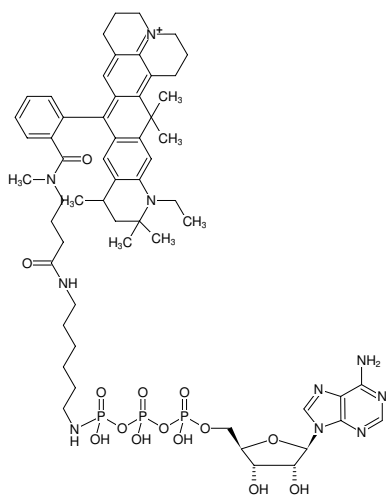




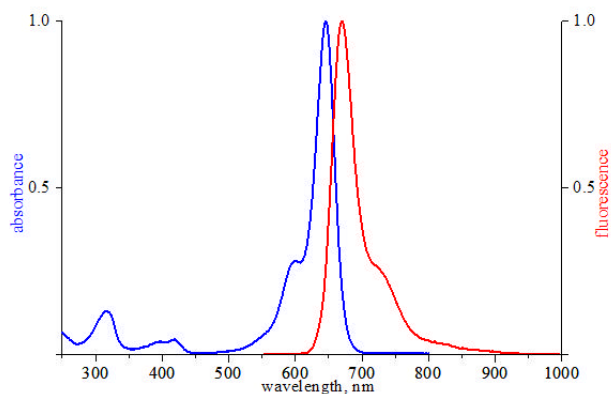
γ -[[(6-Aminoethyl)-imido]-ATP-ATTO-647N

γ -[[(6-Aminoethyl)-imido]-adenosine-5'-triphosphate, labeled with ATTO-647N, Triethylammonium salt

Cat. No.	Amount
NU-814-647N	120 μ l (0,5 mM)



Structural formula of γ -[[(6-Aminoethyl)-imido]-ATP-ATTO-647N



excitation and emission spectrum of ATTO 647N

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: C₅₈H₇₉N₁₀O₁₄P₃ (free acid)

Molecular Weight: 1233.23 g/mol (free acid)

Exact Mass: 1232.50 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: blue

Concentration: 0.50 mM - 0.55 mM

pH: 8.0 \pm 0.5

Spectroscopic Properties: λ_{exc} 646 nm, λ_{em} 664 nm, ϵ 150.0 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.

Selected References:

Götz *et al.* (2016) A Multicolor Single-Molecule FRET Approach to Study Protein Dynamics and Interactions Simultaneously. *Methods Enzymol.* **581**:487.