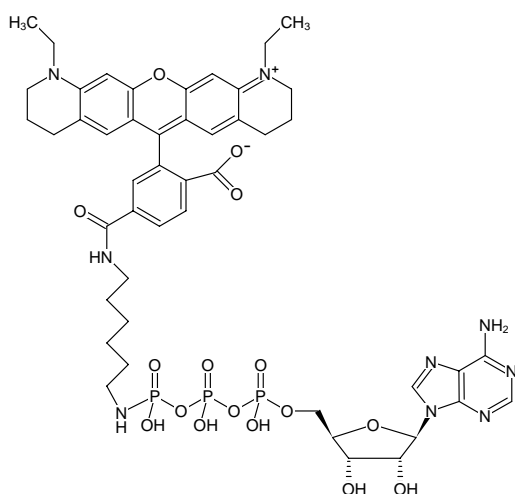




γ -[(6-Aminoethyl)-imido]-ATP-ATTO-565

γ -[(6-Aminoethyl)-imido]-adenosine-5'-triphosphate, labeled with ATTO-565, Triethylammonium salt

Cat. No.	Amount
NU-814-565	120 μ l (1 mM)



Structural formula of γ -[(6-Aminoethyl)-imido]-ATP-ATTO-565

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: C₄₇H₅₈N₉O₁₆P₃ (free acid)

Molecular Weight: 1097.95 g/mol (free acid)

Exact Mass: 1097.32 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

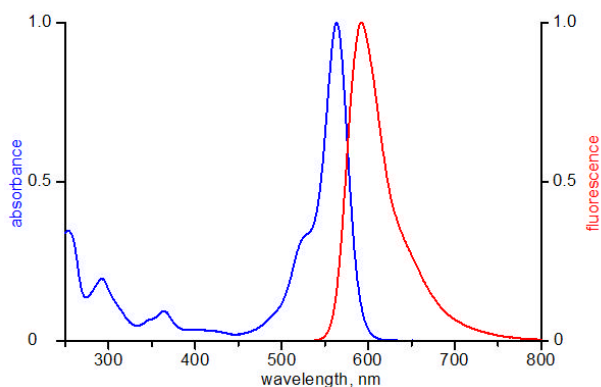
Color: red

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 564 nm, λ_{em} 590 nm, ϵ 120.0 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.



excitation and emission spectrum of ATTO 565