



γ -(6-Aminoethyl)-ATP-ATTO-Rho11

γ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-833-RHO11	80 μ l (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume \leq 2 ml).

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: C₅₁H₆₆N₉O₁₆P₃ (free acid)

Molecular Weight: 1154.06 g/mol (free acid)

Exact Mass: 1153.38 g/mol (free acid)

Purity: \geq 95 % (HPLC)

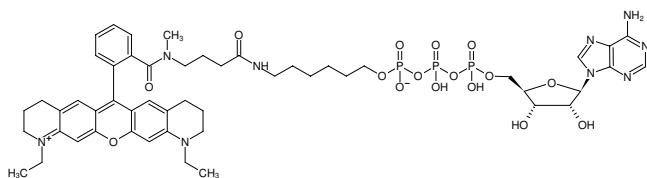
Form: solution in water

Color: violet

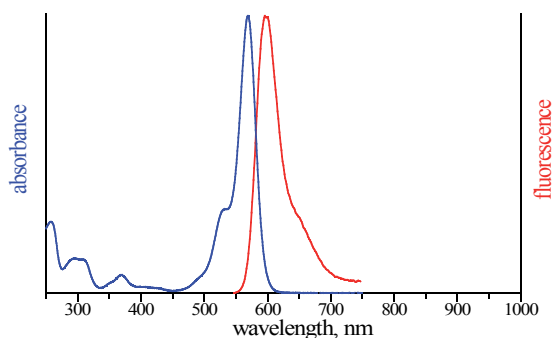
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 572 nm, λ_{em} 595 nm, ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of γ -(6-Aminoethyl)-ATP-ATTO-Rho11



excitation and emission spectrum of ATTO Rho11