



8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho101

8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate, labeled with ATTO Rho101, Triethylammonium salt

Cat. No.	Amount
NU-812-RHO101	40 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤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: C₅₃H₆₆N₁₀O₁₃P₂ (free acid)

Molecular Weight: 1113.12 g/mol (free acid)

Exact Mass: 1112.43 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

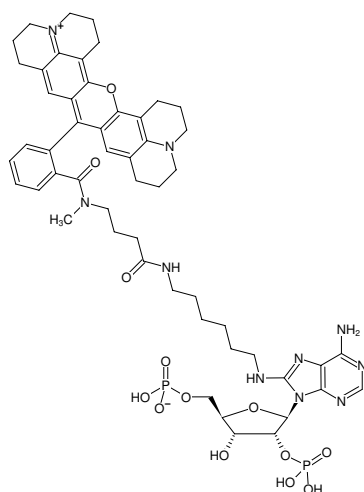
Form: solution in water

Color: violet

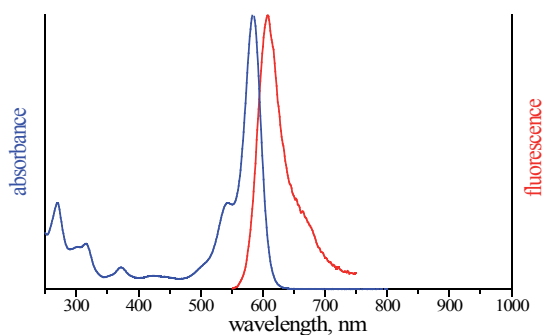
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 587 nm, λ_{em} 609 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho101



excitation and emission spectrum of ATTO Rho101