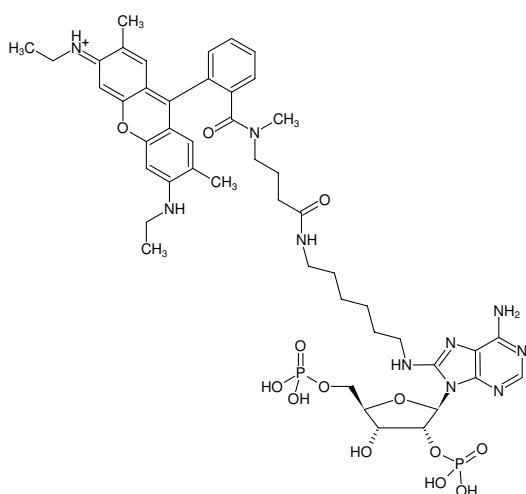




8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho6G

8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate, labeled with ATTO Rho6G, Triethylammonium salt

| Cat. No. | Amount |
|--------------|--------------|
| NU-812-RHO6G | 40 µl (1 mM) |



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho6G

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: C₄₇H₆₂N₁₀O₁₃P₂ (free acid)

Molecular Weight: 1037.02 g/mol (free acid)

Exact Mass: 1036.40 g/mol (free acid)

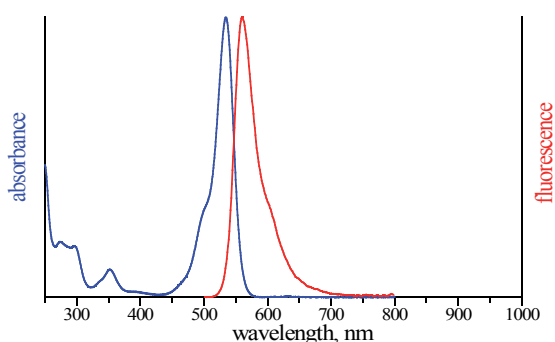
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 533 nm, λ_{em} 557 nm, ε 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Rho6G