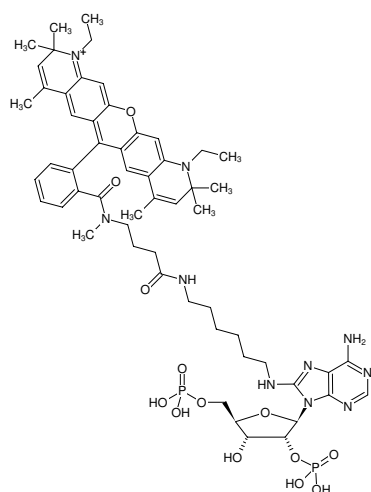




8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho13

8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate, labeled with ATTO Rho13, Triethylammonium salt

Cat. No.	Amount
NU-812-RHO13	40 μ l (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho13

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: C₅₇H₇₄N₁₀O₁₃P₂ (free acid)

Molecular Weight: 1169.22 g/mol (free acid)

Exact Mass: 1168.49 g/mol (free acid)

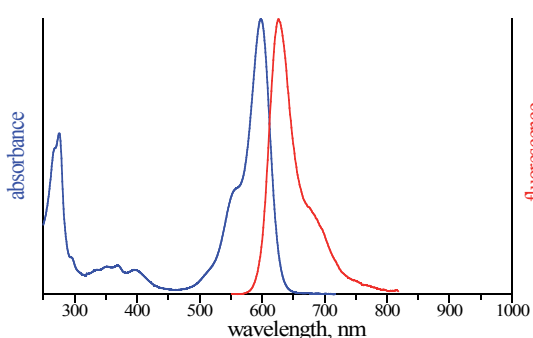
Purity: \geq 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 627 nm, ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Rho13