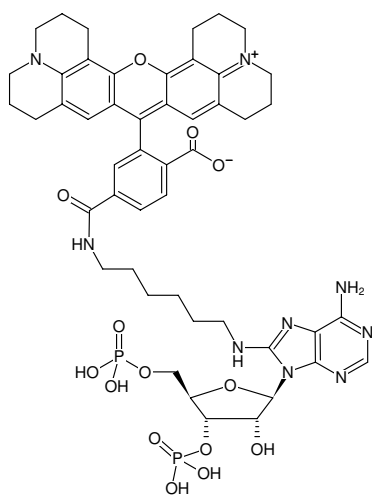




8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-6-ROX

8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate, labeled with 6-ROX, Triethylammonium salt

Cat. No.	Amount
NU-811-ROX	160 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-6-ROX

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: C₄₉H₅₇N₉O₁₄P₂ (free acid)

Molecular Weight: 1057.99 g/mol (free acid)

Exact Mass: 1057.35 g/mol (free acid)

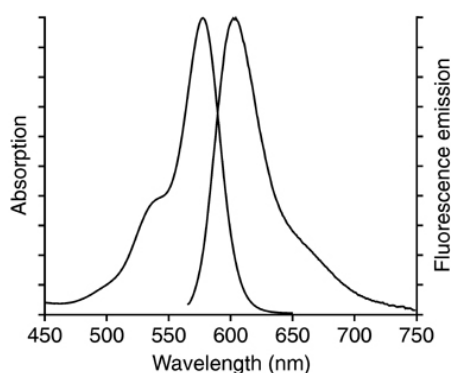
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 575 nm, λ_{em} 600 nm, ε 82.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of 6-ROX