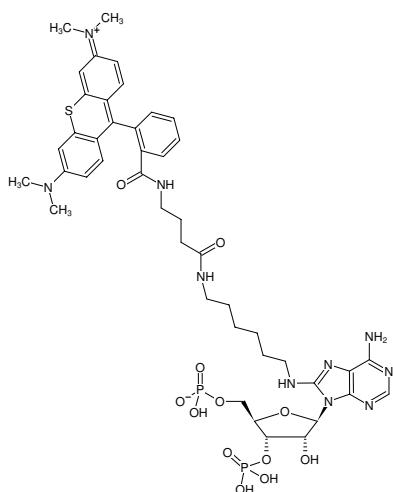




8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-ATTO-Thio12

8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-811-THIO12	40 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-ATTO-Thio12

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: C₄₅H₅₈N₁₀O₁₂P₂S (free acid)

Molecular Weight: 1025.02 g/mol (free acid)

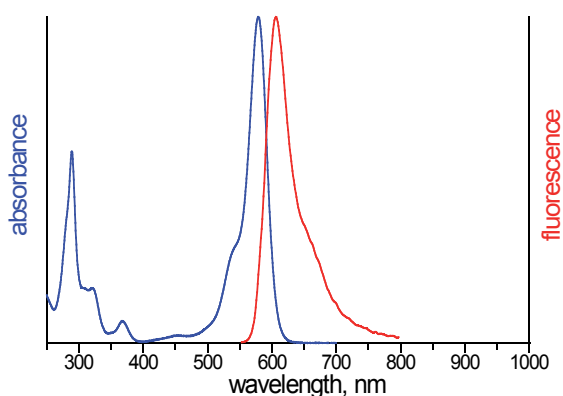
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 582 nm, λ_{em} 607 nm, ε 110.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Thio12