



8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-ATTO-612Q

8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate, labeled with ATTO 612Q, Triethylammonium salt

Cat. No.	Amount
NU-811-612Q	40 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

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: C₁₆H₂₉N₇O₁₀P₂ - ATTO 612Q (free acid)

Molecular Weight: 1213.39 g/mol (free acid)

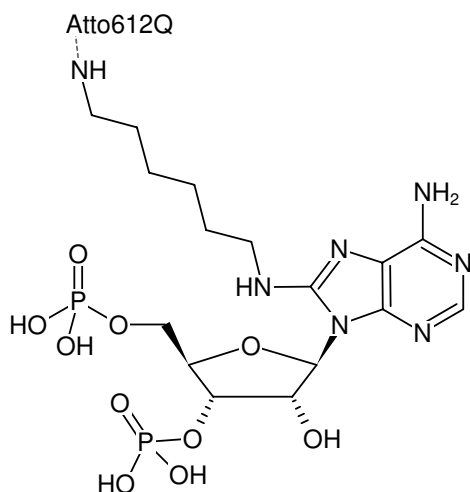
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

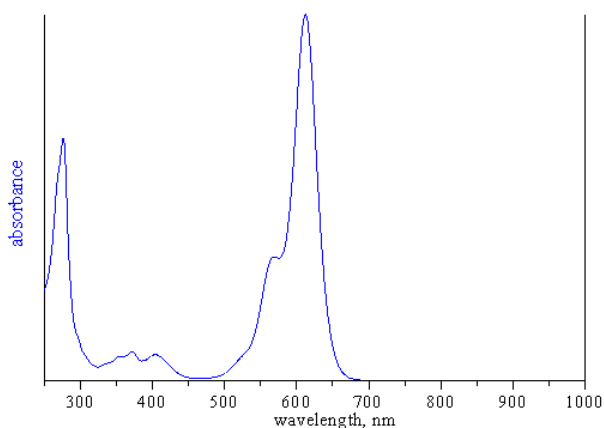
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 615 nm, ε 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-ATTO-612Q



absorption spectrum of ATTO 612Q