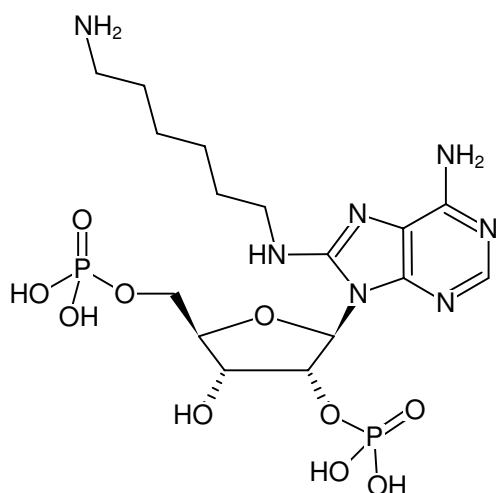




8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate

8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-812S	30 µl (10 mM)
NU-812L	5 x 30 µl (10 mM)



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

For research use only!

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: C₁₆H₂₉N₇O₁₀P₂ (free acid)

Molecular Weight: 541.39 g/mol (free acid)

Exact Mass: 541.15 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 273 nm, ε 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)