



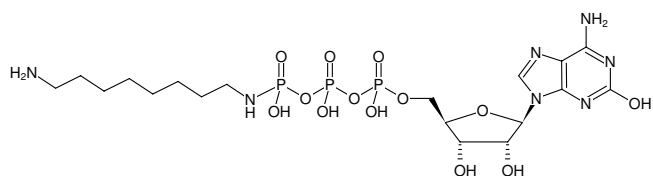
γ -[(8-Aminoethyl)-imido]-2-Hydroxy-ATP

γ -[(8-Aminoethyl)-imido]-2-hydroxy-adenosine-5'-triphosphate, Sodium salt

γ -[(8-Aminoethyl)-imido]-iso-GTP

Cat. No.	Amount
NU-843S	1 mg
NU-843L	5 mg

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.



Structural formula of γ -[(8-Aminoethyl)-imido]-2-Hydroxy-ATP

Please note: This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.

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: C₁₈H₃₄N₇O₁₃P₃ (free acid)

Molecular Weight: 649.43 g/mol (free acid)

Exact Mass: 649.14 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{\max} 292 nm, ϵ 11.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)