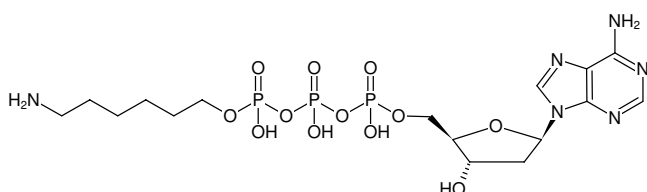




γ -(6-Aminoethyl)-dATP

γ -(6-Aminoethyl)-2'-deoxyadenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-864S	50 μ l (10 mM)
NU-864L	5 x 50 μ l (10 mM)



Structural formula of γ -(6-Aminoethyl)-dATP

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: C₁₆H₂₉N₆O₁₂P₃ (free acid)

Molecular Weight: 590.36 g/mol (free acid)

Exact Mass: 590.11 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{\max} 259 nm, ϵ 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)