

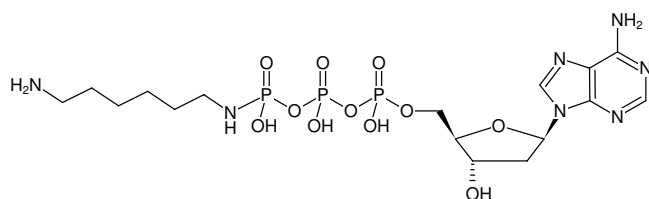


γ -[(6-Aminohexyl)-imido]-dATP

γ -[(6-Aminohexyl)-imido]-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-861-5	5 mg
NU-861-25	25 mg

Please note: This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.



Structural formula of γ -[(6-Aminohexyl)-imido]-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: C₁₆H₃₀N₇O₁₁P₃ (free acid)

Molecular Weight: 589.38 g/mol (free acid)

Exact Mass: 589.12 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{\max} 259 nm, ϵ 15.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)