



γ -[(6-Aminohexyl)-imido]-dGTP

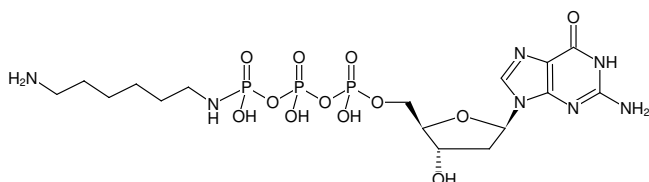
γ -[(6-Aminohexyl)-imido]-2'-deoxyguanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-849-5	5 mg
NU-849-25	25 mg

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

Selected References:

van Gijlswijk *et al.* (2002) Use of horseradish peroxidase- and fluorescein-modified cisplatin derivatives for simultaneous labeling of nucleic acids and proteins. *Clin Chem.* **48** (8):1352.



Structural formula of γ -[(6-Aminohexyl)-imido]-dGTP

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: C₁₆H₃₀N₇O₁₂P₃ (free acid)

Molecular Weight: 605.37 g/mol (free acid)

Exact Mass: 605.12 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 252 nm, ϵ 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)