

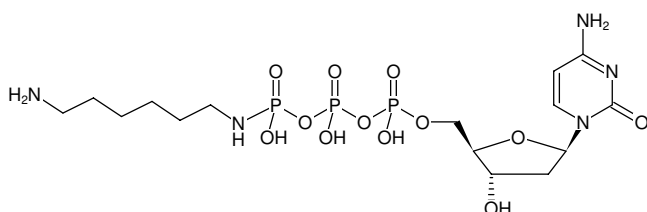


γ -[(6-Aminoethyl)-imido]-dCTP

γ -[(6-Aminoethyl)-imido]-2'-deoxycytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-862-5	5 mg
NU-862-25	25 mg

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



Structural formula of γ -[(6-Aminoethyl)-imido]-dCTP

For research use only!

Shipping: shipped on blue ice

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: C₁₅H₃₀N₅O₁₂P₃ (free acid)

Molecular Weight: 565.35 g/mol (free acid)

Exact Mass: 565.11 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Spectroscopic Properties: λ_{\max} 271 nm, ϵ 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)