



γ -[(8-Aminoethyl)-imido]-dUTP

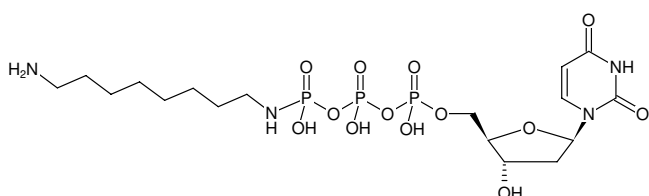
γ -[(8-Aminoethyl)-imido]-2'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-818-5	5 mg
NU-818-25	25 mg

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

Selected References:

Suto *et al.* (1998) Synthesis of γ -phosphate-linked nucleoside affinity chromatography resins for protein purification, including ribonucleoside triphosphate reductase. *Nucleosides and Nucleotides* **17** (8):1453.



Structural formula of γ -[(8-Aminoethyl)-imido]-dUTP

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: C₁₇H₃₃N₄O₁₃P₃ (free acid)

Molecular Weight: 594.39 g/mol (free acid)

Exact Mass: 594.13 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid