

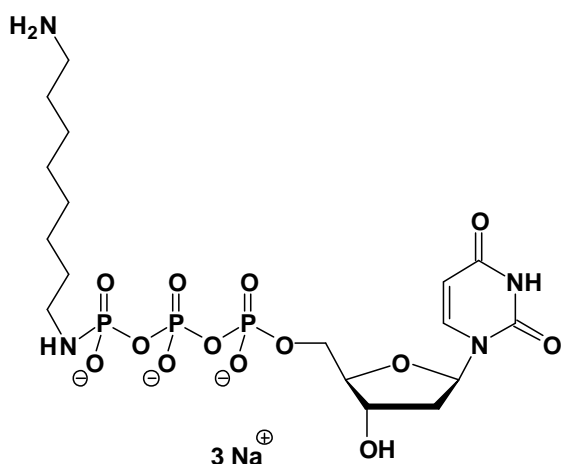
γ -[(8-Amino-octyl)-imido]-dUTP

γ -[(8-Amino-octyl)-imido]-2'-deoxy-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-818S	50 Units
NU-818L	250 Units

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.



Cat. No.: NU-818

Molecular Formula: C₁₇H₃₀N₄O₁₃P₃ (Anion)

Molecular Weight: 579.31 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution