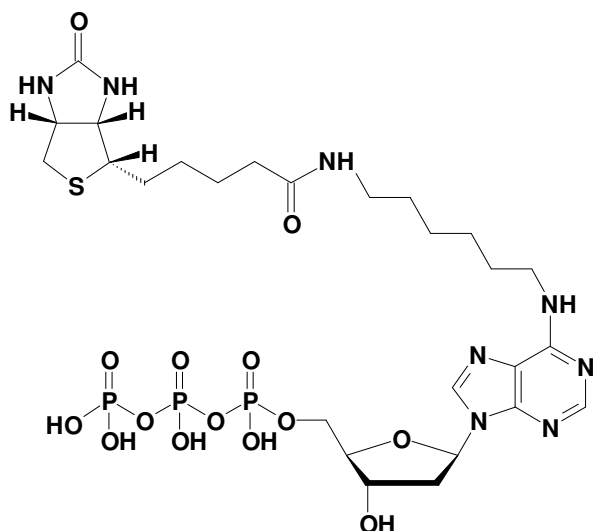


N⁶-(6-Amino)hexyl-dATP - Biotin

N⁶-(6-Amino)hexyl-2'-deoxy-adenosine-5'-triphosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-835-BIO	50 µl / 5 mM



Applications:

Incorporation by terminal deoxynucleotidyl transferase^[1]

Selected References:

[1] Flickinger *et al.* (1992) Differential incorporation of biotinylated nucleotides by terminal deoxynucleotidyl transferase. *Nucleic Acids Res.* **20**:2382.

Cat. No.: NU-835-BIO

Molecular Formula: C₂₆H₄₃N₈O₁₄P₃S (free acid)

Molecular Weight: 816.65 (free acid)

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 6 months after date of delivery.

For research use only!