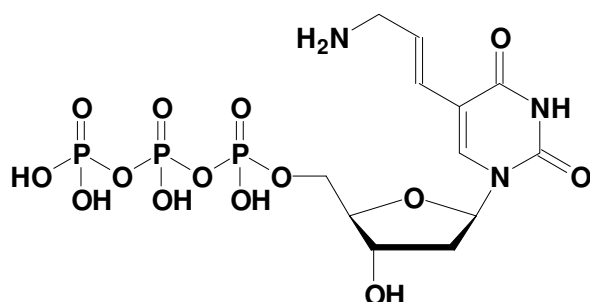


Aminoallyl-dUTP, lyophilized

5-(3-Aminoallyl)-2'-deoxy-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-803-1	1 mg
NU-803-5	5 mg



Cat. No.: NU-803

Molecular Formula: $C_{12}H_{20}N_3O_{14}P_3$ (free acid)

Molecular Weight: 523.22 (free acid)

Purity: > 95%, lyophilized solid

Spectroscopic Properties: λ_{max} 289nm;
 ϵ 7,100 $M^{-1}cm^{-1}$ Tris HCl 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!

Selected References:

Cherkasov *et al.* (2010) New Nucleotide Analogues with Enhanced Signal Properties. *Bioconjugate Chem.* **21** (1):122.

Kuwahara *et al.* (2006) Direct PCR amplification of various modified DNAs having amino acids: Convenient preparation of DNA libraries with high-potential activities for in vitro selection. *Bioorganic & Medicinal Chemistry* **14**:2518.

Nimmakayalu *et al.* (2000) Simple method for preparation of fluor/hapten-labeled dUTP. *Biotechniques* **28**:518.

Muhlegger *et al.* (1990) Non-radioactive labeling and detection of nucleic acids. IV. Synthesis and properties of digoxigeninmodified 2'-deoxyuridine-5'-triphosphates and a photoactivatable analog of digoxigenin (photodigoxigenin). *Biol. Chem. Hoppe Seyler* **371**:953.

Giaid *et al.* (1989) Non-isotopic RNA probes. Comparison between different labels and detection systems. *Histochemistry* **93**:191.