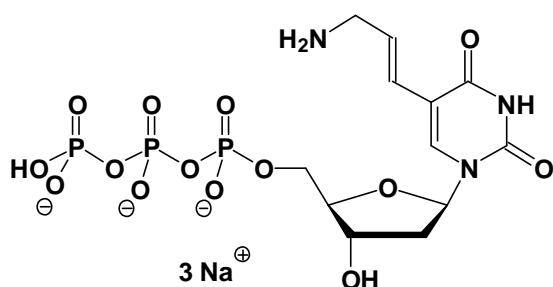


## Aminoallyl-dUTP

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

Cat. No.	Amount
NU-803S	50 Units
NU-803L	250 Units



**Cat. No.:** NU-803

**Molecular Formula:** C<sub>12</sub>H<sub>17</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 520.19 (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 µl of a 10 mM solution

### Selected References:

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.