



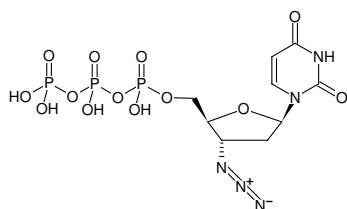
### 3'-Azido-2',3'-ddUTP

3'-Azido-2',3'-dideoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-251S	10 µl (100 mM)
NU-251L	5 x 10 µl (100 mM)

#### Selected References:

Eriksson *et al.* (1989) Phosphorylation of 3'-azido-2',3'-dideoxyuridine and preferential inhibition of human and simian immunodeficiency virus reverse transcriptases by its 5'-triphosphate. *Antimicrob. Agents Chemother.* **33** (10):1729.



Structural formula of 3'-Azido-2',3'-ddUTP

**For general laboratory use.**

**Shipping:** shipped on gel packs

**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:** C<sub>9</sub>H<sub>14</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 493.15 g/mol (free acid)

**Exact Mass:** 492.98 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 100 mM - 110 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 10.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)