



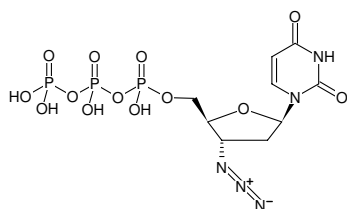
3'-Azido-2',3'-ddUTP

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 research use only!

Shipping: shipped on blue ice

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: C₉H₁₄N₅O₁₃P₃ (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_{max} 262 nm, ε 10.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)