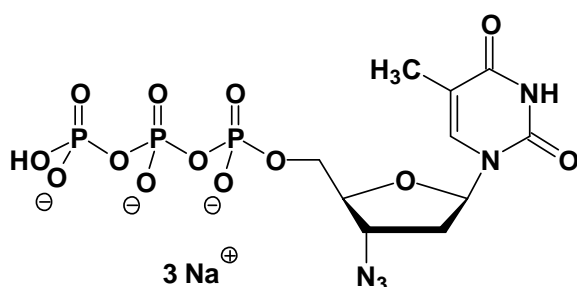


AzTTP

3'-Azido-2',3'-dideoxy-thymidine-5'-triphosphate, Sodium salt
(Zidovudine triphosphate)

Cat. No.	Amount
NU-1602S	10 Units
NU-1602L	50 Units



Cat. No.: NU-1602

Molecular Formula: C₁₀H₁₃N₅O₁₃P₃ (Anion)

Molecular Weight: 504.16 (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:

Cruchaga *et al.* (2005) Inhibition of Phosphorolysis Catalyzed by HIV-1 Reverse Transcriptase Is Responsible for the Energy Found in Combinations of 3'-Azido-3'-deoxythymidine with Nonnucleoside Inhibitors *Biochemistry*. **44** (9):3535.

Akeb *et al.* (2001) The production and evaluation of antibodies for enzyme immunoassay of AZTTP. *Nucleosides Nucleotides Nucleic Acids*. **20**:243.

Fletcher *et al.* (2000) Zidovudine triphosphate and lamivudine triphosphate concentration-response relationships in HIV-infected persons. *AIDS* **14**:2137.

Faraj *et al.* (2000) Effects of beta-L-3'-azido-3'-deoxythymidine 5'-triphosphate on host and viral DNA polymerases. *Antiviral Res.* **47**:97.

Font *et al.* (1999) Determination of zidovudine triphosphate intracellular concentrations in peripheral blood mononuclear cells from human immunodeficiency virus-infected individuals by tandem mass spectrometry. *Antimicrob. Agents Chemother.* **43**:2964.

Elimadi *et al.* (1997) Differential effects of zidovudine and zidovudine triphosphate on mitochondrial permeability transition and oxidative phosphorylation. *Br. J. Pharmacol.* **121**:1295.

Jaju *et al.* (1995) Human immunodeficiency virus type 1 reverse transcriptase. 3'-Azidodeoxythymidine 5'-triphosphate inhibition indicates two-step binding for template-primer. *J. Biol. Chem.* **270**:9740.

Tornevik *et al.* (1995) Cytotoxicity of 3'-azido-3'-deoxythymidine correlates with 3'-azidothymidine-5'-monophosphate (AZTMP) levels, whereas anti-human immunodeficiency virus (HIV) activity correlates with 3'-azidothymidine-5'-triphosphate (AZTTP) levels in cultured CEM T-lymphoblastoid cells. *Biochem. Pharmacol.* **49**:829.