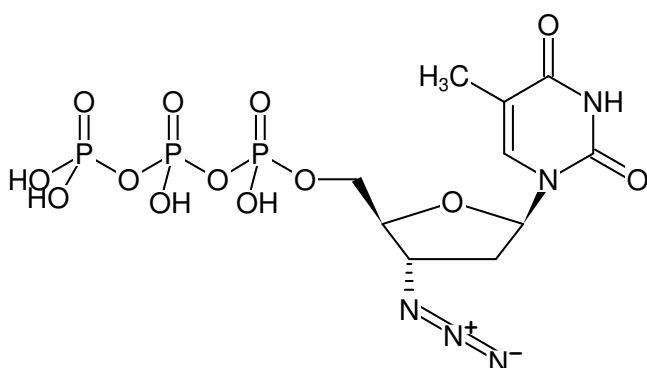


**AzTTP**

Zidovudine triphosphate

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

Cat. No.	Amount
NU-989S	10 µl (100 mM)
NU-989L	5 x 10 µl (100 mM)



Structural formula of AzTTP

**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:** C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 507.18 g/mol (free acid)**Exact Mass:** 507.00 g/mol (free acid)**CAS#:** 906479-25-2**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>max</sub> 267 nm, ε 10.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**Suppression of HIV-type 1 replication<sup>[1]</sup>Removal of iron from transferrin<sup>[2]</sup>Inhibition of telomerase activity<sup>[3]</sup>Inhibition of thymidine phosphorylation<sup>[4]</sup>**Specific Ligands:**High binding affinity to HIV-1 mutants<sup>[5]</sup>**Selected References:**

[1] Sayed *et al.* (2009) AZT 5'-triphosphate nanoformulation suppresses human immunodeficiency virus type 1 replication in peripheral blood mononuclear cells. *J. Neurobiology* **15** (4):343.

[2] D'Andrea *et al.* (2008) AZT: an old drug with new perspectives. *Curr. Clin. Pharmacol.* **3** (1):20.

[3] Liu *et al.* (2007) 3'-Azido-2',3'-dideoxynucleoside 5'-triphosphates inhibit telomerase activity in vitro, and the corresponding nucleosides cause telomere shortening in human HL60 cells. *Nucleic Acid Res.* **35** (21):7140.

[4] Lynx *et al.* (2006) 3'-Azido-3'-deoxythymidine (AZT) inhibits thymidine phosphorylation in isolated rat liver mitochondria: A possible mechanism of AZT hepatotoxicity. *Biochem. Pharmacol.* **71** (9):1342.

[5] Jamburuthugoda *et al.* (2005) Kinetic evidence for interaction of HIV type 1 reverse transcriptase with the 3'-OH of the incoming dTTP substrate. *Biochemistry* **44** (31):10635.

Coulier *et al.* (2008) Simultaneous determination of endogenous deoxynucleotides and phosphorylated nucleoside reverse transcriptase inhibitors in peripheral blood mononuclear cells using ion-pair liquid chromatography coupled to mass spectrometry. *Proteomics Clin. Appl.* **2**:1557.

Cruchaga *et al.* (2005) Inhibition of Phosphorolysis Catalyzed by HIV-1 Reverse Transcriptase Is Responsible for the Synergy 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.

**AzTTP**

Zidovudine triphosphate

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

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.