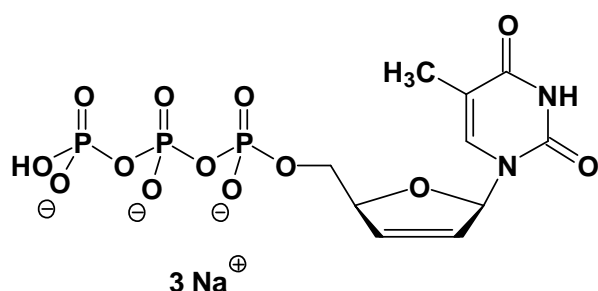


d4TTP

**2',3'-Didehydro-2',3'-dideoxy-thymidine-5'-triphosphate, Sodium salt
(Stavudine triphosphate)**

Cat. No.	Amount
NU-1604S	10 Units
NU-1604L	50 Units



Cat. No.: NU-1604

Molecular Formula: $\text{C}_{10}\text{H}_{12}\text{N}_2\text{O}_{13}\text{P}_3$ (Anion)

Molecular Weight: 461.13 (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:

Ray *et al.* (2002) Insights into the molecular mechanism of inhibition and drug resistance for HIV-1 RT with carbovir triphosphate. *Biochemistry* **41**:5150.

Hoggard *et al.* (2000) Correlation between intracellular pharmacological activation of nucleoside analogues and HIV suppression in vitro. *Antivir. Chem. Chemother.* **11**:353.

Vaccaro *et al.* (2000) Mechanism of inhibition of the human immunodeficiency virus type 1 reverse transcriptase by d4TTP: an equivalent incorporation efficiency relative to the natural substrate dTTP. *Antimicrob. Agents. Chemother.* **44**:217.

Ueno *et al.* (1997) Comparative enzymatic study of HIV-1 reverse transcriptase resistant to 2',3'-dideoxynucleotide analogs using the single-nucleotide incorporation assay. *Biochemistry* **36**:1092.

Im *et al.* (1993) Identification of the amino acid in the human immunodeficiency virus type 1 reverse transcriptase involved in the pyrophosphate binding of antiviral nucleoside triphosphate analogs and phosphonoformate. Implications for multiple drug resistance. *Biochem. Pharmacol.* **46**:2307.

Huang *et al.* (1992) Selective action of 2',3'-didehydro-2',3'-dideoxythymidine triphosphate on human immunodeficiency virus reverse transcriptase and human DNA polymerases. *J. Biol. Chem.* **267**:2817.