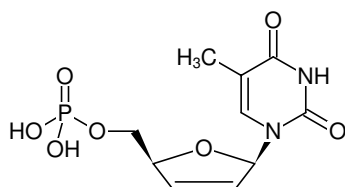


**d4TMP**

Stavudine monophosphate, Sodium Salt

2',3'-Didehydro-2',3'-dideoxythymidine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1603S	20 µl (10 mM)
NU-1603L	5 x 20 µl (10 mM)



Structural formula of d4TMP

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:** C₁₀H₁₃N₂O₇P (free acid)**Molecular Weight:** 304.19 g/mol (free acid)**Exact Mass:** 304.05 g/mol (free acid)**CAS#:** 3715-64-8**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 266 nm, ε 10.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

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.

De Clercq (2002) Highlights in the development of new antiviral agents. *Mini Rev. Med. Chem.* **2**:163.

Mas *et al.* (2002) Multidrug-resistant HIV-1 reverse transcriptase: involvement of ribonucleotide-dependent phosphorolysis in cross-resistance to nucleoside analogue inhibitors. *J. Mol. Biol.* **267**:181.

Selmi *et al.* (2001) The valine-to-threonine 75 substitution in human immunodeficiency virus type 1 reverse transcriptase and its relation with stavudine resistance. *J. Biol. Chem.* **276**:13965.

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.