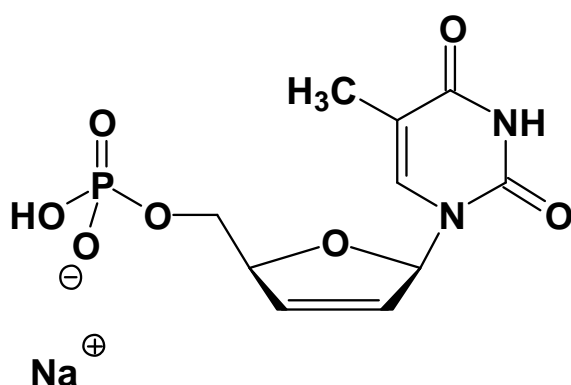


## d4TMP

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

Cat. No.	Amount
NU-1603S	20 Units
NU-1603L	100 Units



**Cat. No.:** NU-1603

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>2</sub>O<sub>7</sub>P (Anion)

**Molecular Weight:** 303.18 (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:

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.