

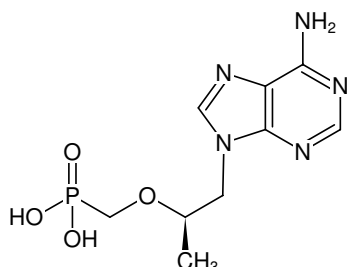
**Tenofovir**

PMPA

[(2R)-1-(6-aminopurin-9-yl)propan-2-yl]oxymethylphosphonic acid

(R)-9-(2-phosphonylmethoxypropyl)adenine

Cat. No.	Amount
NU-974	100 mg



Structural formula of Tenofovir

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:** C₉H₁₄N₅O₄P (free acid)**Molecular Weight:** 287.21 g/mol (free acid)**Exact Mass:** 287.08 g/mol (free acid)**CAS#:** 147127-20-6**Purity:** ≥ 98 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Krejčová *et al.* (2000) Phosphorylation of Purine (Phosphonomethoxy)alkyl Derivatives by Mitochondrial AMP Kinase (AK2 Type) from L1210 Cells. *Collect. Czech. Chem. Commun.* **65**:1653.

Robbins *et al.* (1998) Anti-human immunodeficiency virus activity and cellular metabolism of a potential prodrug of the acyclic nucleoside phosphonate 9-R-(2-phosphonomethoxypropyl)adenine (PMPA), Bis(isopropylloxymethylcarbonyl)PMPA. *Antimicrob. Agents Chemother.* **42**(3):612.

Robbins *et al.* (1995) Metabolic Pathways for Activation of the Antiviral Agent 9-(2-Phosphonylmethoxyethyl)Adenine in Human Lymphoid Cells. *Antimicrob. Agents Chemother.* **39**(10):2304.