



Tenofovir-diphosphate

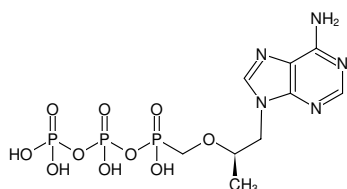
PMPApp

Tenofovir-DP, Sodium salt

[2-(6-Amino-9H-purin-9-yl)-1-methylethoxy]methyl-triphosphate;

[(1R)-2-(6-aminopurin-9-yl)-1-methyl-ethoxy]methyl-[hydroxy(phosphonoxy)phosphoryl]oxy-phosphinic acid

Cat. No.	Amount
NU-975	1 mg



Structural formula of Tenofovir-diphosphate

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: C₉H₁₆N₅O₁₀P₃ (free acid)

Molecular Weight: 447.17 g/mol (free acid)

Exact Mass: 447.01 g/mol (free acid)

CAS#: 166403-66-3

Purity: ≥ 95 % (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:

Zheng *et al.* (2016) Application of an intracellular assay for determination of Tenofovir-diphosphate and Emtricitabine-triphosphate from erythrocytes using Dried Blood Spots. *J. Pharm. Biomed. Anal.* **122**:16.

Dickinson *et al.* (2015) Plasma Tenofovir, Emtricitabine, and Rilpivirine and Intracellular Tenofovir Diphosphate and Emtricitabine Triphosphate Pharmacokinetics following Drug Intake Cessation. *Antimicrob. Agents Chemother.* **59** (10):6080.