



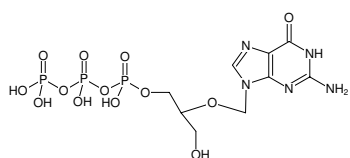
Ganciclovir-triphosphate

Sodium salt

GCV-TP

9-(1,3-Dihydroxy-2-propoxymethyl)guanine-5'-triphosphate

Cat. No.	Amount
NU-275S	10 µl (100 mM)
NU-275L	5 x 10 µl (100 mM)



Structural formula of Ganciclovir-triphosphate

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: C₉H₁₆N₅O₁₃P₃ (free acid)

Molecular Weight: 495.17 g/mol (free acid)

Exact Mass: 495.00 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 252 nm, ε 13.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Davis *et al.* (2007) Hypoxia Enhances the Phosphorylation and Cytotoxicity of Ganciclovir and Zidovudine in Kaposi's Sarcoma-Associated Herpesvirus-Infected Cells. *Cancer Res.* **67**:7003.

Zhu *et al.* (2005) Potent Antiviral Activity of North-Methanocarbothymidine against Kaposi's Sarcoma-Associated Herpesvirus. *Antimicrob. Agents Chemother.* **49**:4965.

Rubsam *et al.* (1999) Cytotoxicity and accumulation of ganciclovir triphosphate in bystander cells cocultured with herpes simplex virus type 1 thymidine kinase-expressing human glioblastoma cells. *Cancer Res.* **59**:669.