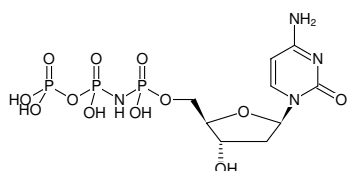


**dCpNHpp**

(dCMPNPP)

2'-Deoxycytidine-5'-[(α,β)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-439-1	1 mg



Structural formula of dCpNHpp

For research use only!**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:** C₉H₁₇N₄O₁₂P₃ (free acid)**Molecular Weight:** 466.17 g/mol (free acid)**Exact Mass:** 466.01 g/mol (free acid)**CAS#:** 791761-70-1 (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 271 nm, ϵ 9.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Moscato *et al.* (2016) Induced Fit in the Selection of Correct versus Incorrect Nucleotides by DNA Polymerase β . *Biochemistry* **55** (2):382.Patra *et al.* (2015) Structural Basis for Error-Free Bypass of the 5-N-Methylformamidopyrimidine-dG Lesion by Human DNA Polymerase η and *Sulfolobus solfataricus* P2 Polymerase IV. *J. Am. Chem. Soc.* **137** (22):7011.