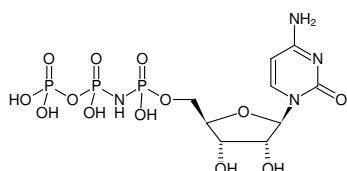


**CpNHpp**

(CMPNPP)

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

Cat. No.	Amount
NU-264	1 mg



Structural formula of CpNHpp

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:** C₉H₁₇N₄O₁₃P₃ (free acid)**Molecular Weight:** 482.17 g/mol (free acid)**Exact Mass:** 482.00 g/mol (free acid)**CAS#:** 497064-76-3 (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)