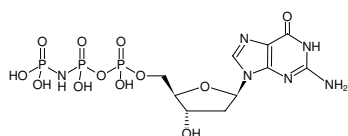


**dGppNHp**

(dGMPPNP)

2'-Deoxyguanosine-5'-[(β,γ)-imido]triphosphate, Lithium salt

Cat. No.	Amount
NU-258	5 mg



Structural formula of dGppNHp

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: 6 months after date of delivery**Molecular Formula:** C₁₀H₁₇N₆O₁₂P₃ (free acid)**Molecular Weight:** 506.20 g/mol (free acid)**Exact Mass:** 506.01 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 252 nm, ϵ 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)