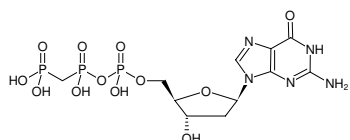


**dGppCp**

(dGMPPCP)

2'-Deoxyguanosine-5'-[(β,γ)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-260	5 mg



Structural formula of dGppCp

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:** C₁₁H₁₈N₅O₁₂P₃ (free acid)**Molecular Weight:** 505.21 g/mol (free acid)**Exact Mass:** 505.02 g/mol (free acid)**Purity:** ≥ 95 % (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)