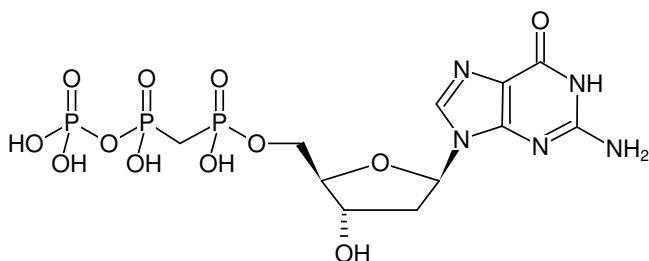


**dGpCpp**

(dGMPCPP)

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

Cat. No.	Amount
NU-431S	50 μ l (10 mM)
NU-431L	5 x 50 μ l (10 mM)



Structural formula of dGpCpp

For research use only!**Shipping:** shipped on blue ice**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**CAS#:** 941569-34-2**Purity:** \geq 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 252 nm, ϵ 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)