

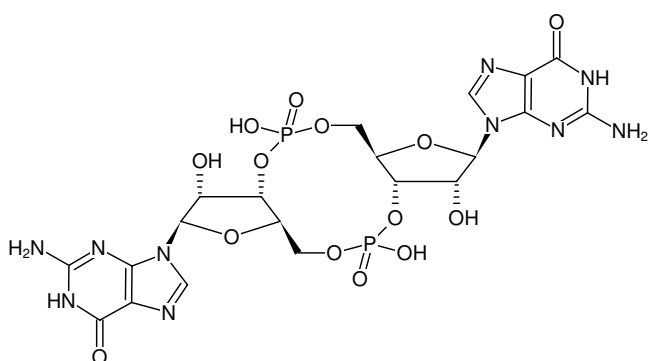
**cyclic-di-GMP**

(c-di-GMP)

cyclic diguanosine monophosphate, Sodium salt

cyclic bis(3'→5')diguanylic acid

Cat. No.	Amount
NU-951S	1 µmol
NU-951L	5 x 1 µmol



Structural formula of cyclic-di-GMP

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:** C₂₀H₂₄N₁₀O₁₄P₂ (free acid)**Molecular Weight:** 690.41 g/mol (free acid)**Exact Mass:** 690.09 g/mol (free acid)**CAS#:** 61093-23-0 (free acid)**Purity:** ≥ 98 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 253 nm, ε 23.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)