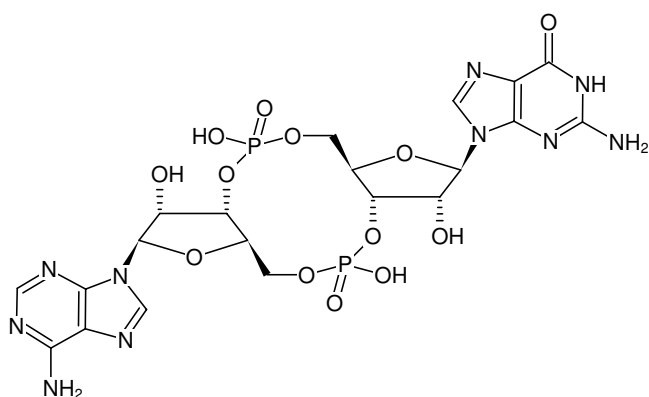


**3',3'-cGAMP**

3',3'-cyclic GMP-AMP

Adenosine-Guanosine-3',3'-cyclic monophosphate, Sodium salt

Cat. No.	Amount
NU-986S	1 μ mol
NU-986L	5 x 1 μ mol



Structural formula of 3',3'-cGAMP

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:** C₂₀H₂₄N₁₀O₁₃P₂ (free acid)**Molecular Weight:** 674.41 g/mol (free acid)**Exact Mass:** 674.10 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 260 nm, ϵ 25.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)