

**3',3',3'-cAAG**

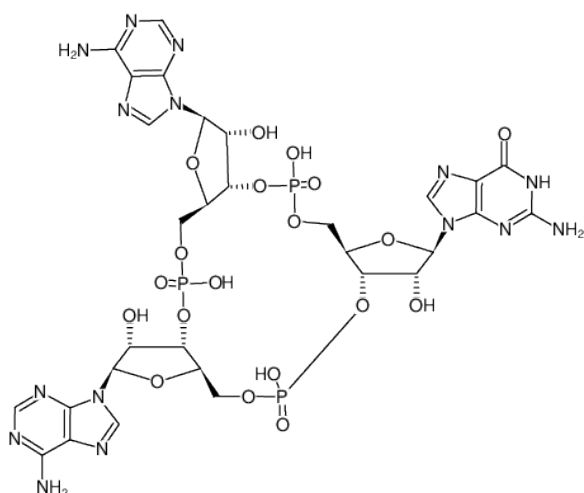
3',3',3'-cyclic AMP-AMP-GMP

3',3',3'-Cyclic adenosine monophosphate-adenosine monophosphate-guanosine monophosphate, Sodium salt

Cat. No.	Amount
NU-456S	1 μ mol
NU-456L	5 x 1 μ mol

Selected References:

Wenzl *et al.* (2024) How enzyme-centered approaches are advancing research on cyclic oligo-nucleotides. *FEBS Lett.* **598**(8): 839.



Structural formula of 3',3',3'-cAAG

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:** C₃₀H₃₆N₁₅O₁₉P₃ (free acid)**Molecular Weight:** 1003.63 g/mol (free acid)**Exact Mass:** 1003.15 g/mol (free acid)**CAS#:** 2365165-54-2**Purity:** \geq 98 % (HPLC) / \leq 2.5 EU/ml Endotoxin (LAL test)**Form:** solid (lyophilized from filtered (10 kDa) solution in water)**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 257 nm, ϵ 39.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)