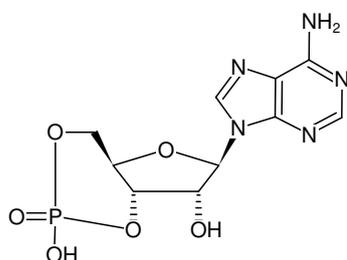


**cAMP**

Adenosine-3',5'-cyclic monophosphate, free acid

Cat. No.	Amount
NU-1503S	500 mg
NU-1503L	5 x 500 mg



Structural formula of cAMP

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:** C₁₀H₁₂N₅O₆P (free acid)**Molecular Weight:** 329.21 g/mol (free acid)**Exact Mass:** 329.05 g/mol (free acid)**CAS#:** 60-92-4**Purity:** ≥ 99 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Chen *et al.* (1998) Up-regulation of the cAMP/PKA pathway inhibits proliferation, induces differentiation, and leads to apoptosis in malignant gliomas. *Lab. Invest.* **78**:165.Daniel *et al.* (1998) Cyclic AMP signaling and gene regulation. *Annu. Rev. Nutr.* **18**:353.Spada *et al.* (1997) Cyclic AMP and calcium in the transduction of hypothalamic neurohormone action in human pituitary tumors. *Horm. Res.* **47**:235.