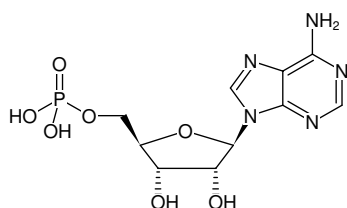


**AMP - Solid**

Adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1025-5G	5 g
NU-1025-25G	25 g



Structural formula of AMP - Solid

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:** C₁₀H₁₄N₅O₇P (free acid)**Molecular Weight:** 347.22 g/mol (free acid)**Exact Mass:** 347.06 g/mol (free acid)**CAS#:** 4578-31-8**Purity:** ≥ 95 % (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)**Applications:**

Agonistic ligand, mainly for nucleoside receptor A_{2A} and less to A₃. After dephosphorylation by different membrane-bound phosphatases the formed adenosine can directly activate the receptor.

Selected References:

Volonte *et al.* (2009) Membrane components and purinergic signalling: the purinome, a complex interplay among ligands, degrading enzymes, receptors and transporters. *FEBS J.* **276**:318.

Yegutkin (2008) Nucleotide and nucleoside converting enzymes: Important modulators of purinergic signalling cascade. *Biochim. Biophys. Acta* **1783**:673.

Van Galen *et al.* (1994) A binding site model and structure-activity relationships for rat A₃ adenosine receptor. *Molecular Pharmacology* **45**:1101.

Williams *et al.* (1986) Effects of purine nucleotides on the binding of [³H]cyclopentyladenosine to adenosine A₁-receptors in rat brain membranes. *J. Neurochem.* **47** (1):88.