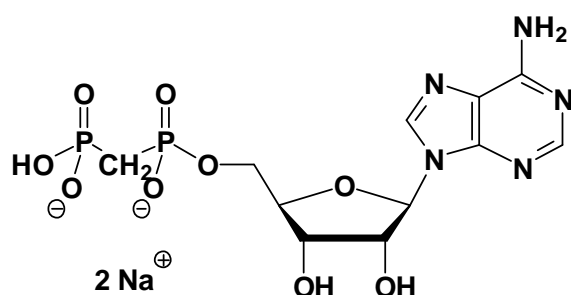


ApCp (AMPCP)

Adenosine-5'-[(α,β)-methyleno]diphosphate, Sodium salt

Cat. No.	Amount
NU-420S	150 Units
NU-420L	750 Units



Cat. No.: NU-420

Molecular Formula: C₁₁H₁₅N₅O₉P₂ (Anion)

Molecular Weight: 423.03 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Cometti *et. al.* (2003) Oviduct cells express the cyclic AMP-adenosine pathway. *Biol. Reprod.* **69** (3):868.

Park *et. al.* (1999) Differential effects of adenine nucleotide analogues on shape change and aggregation induced by adenosine 5-diphosphate (ADP) in human platelets. *Br. J. Pharmacol.* **127** (6):1359.