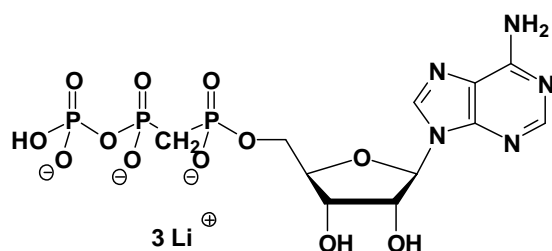


ApCpp (AMPCPP)

Adenosine-5'-[(α,β)-methylene]triphosphate, Lithium salt

Cat. No.	Amount
NU-421S	100 Units
NU-421L	500 Units



Cat. No.: NU-421

Molecular Formula: C₁₁H₁₅N₅O₁₂P₃ (Anion)

Molecular Weight: 501.99 (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:

Bivalacqua *et al.* (2002) Comparative responses to alpha,beta-methylene-ATP in cat pulmonary, mesenteric, and hindquarter vascular beds. *Appl. Physiol.* **93** (4):1287.

Treize (1994) The use of antagonists to characterize the receptors mediating depolarization of the rat isolated vagus nerve by α,β -methylene adenosine 5'-triphosphate. *Br. J. Pharmacol.* **112**:282.

Burnstock *et al.* (1985) Is there a basis for distinguishing two types of P2-purinoceptor? *Gen. Pharmacol.* **16**:433.

Bo (1995) Characterization and autoradiographic localization of [³H] α,β -methylene adenosine 5'-triphosphate binding sites in human urinary bladder. *Br. J. Urol.* **76**:297.