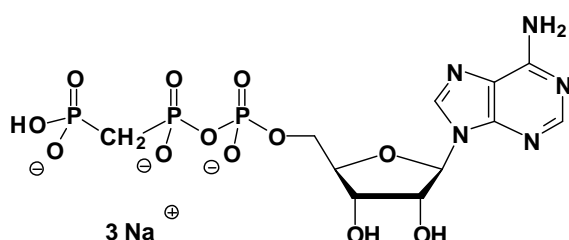


AppCp (AMPPCP)

Adenosine-5'-[(β,γ)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-422S	100 Units
NU-422L	500 Units



Cat. No.: NU-422

Molecular Formula: C₁₁H₁₅N₅O₁₂P₃ (Anion)

Molecular Weight: 501.99 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5
Spectroscopic properties: λ_{\max} 259 nm; ϵ 15400

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:

Toyoshima *et al.* (2004) Crystal structure of the calcium pump with a bound ATP analogue. *Nature* **430**:529.

Trezise *et al.* (1995) The selective P2X purinoceptor agonist, beta,gamma-methylene-L-adenosine 5'-triphosphate, discriminates between smooth muscle and neuronal P2X purinoceptors. *Arch. Pharmacol.* **351** (6):603.

Mayer *et al.* (1991) Re-evaluation of the specificity of adenylyl (beta,gamma-methylene)diphosphonate as a substrate for adenylyl cyclase. *Histochem J.* **23**(2):100.

O'Connor *et al.* (1990) Characterization of P2x-receptors in rabbit isolated ear artery. *Br. J. Pharmacol.* **101** (3):640.

Mayer *et al.* (1985) Specificity of cytochemical demonstration of adenylyl cyclase in liver using adenylyl-(beta, gamma-methylene) diphosphate as substrate. *Histochemistry.* **82**(2):135.