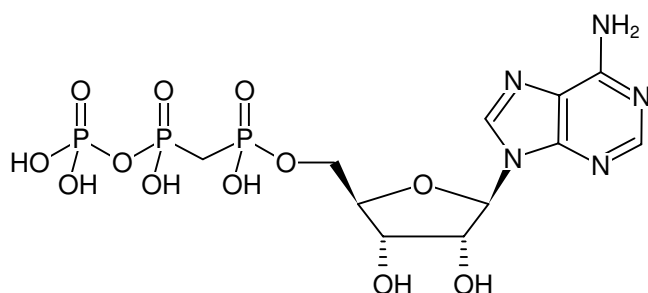


**ApCpp**

(AMPCPP)

 $\alpha,\beta$ -Methyleneadenosine-5'-triphosphateAdenosine-5'-[( $\alpha,\beta$ )-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-421-2	2 mg
NU-421-10	10 mg
NU-421-25	25 mg



Structural formula of ApCpp

**For research use only!****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:** C<sub>11</sub>H<sub>18</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 505.21 g/mol (free acid)**Exact Mass:** 505.02 g/mol (free acid)**CAS#:** 7292-42-4 (acid), 1343364-54-4 (Trisodium salt)**Purity:** ≥ 98 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:**  $\lambda_{\max}$  259 nm,  $\epsilon$  15.4 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Specific Ligands:**Nonhydrolyzable ligand for P2 purinoreceptor<sup>[1]</sup>**Selected References:**[1] Burnstock *et al.* (1985) Is there a basis for distinguishing two types of P2-purinoreceptor? *Gen. Pharmacol.* **16**:433.Zebisch *et al.* (2012) Crystallographic Evidence for a Domain Motion in Rat Nucleoside Triphosphate Diphosphohydrolase (NTPDase) 1. *Journal of molecular biology* **415** (2):288.Riballo *et al.* (2009) XLF-Cernunos promotes DNA ligase IV $\beta$ XRCC4 re-adenylation following ligation. *Nucleic Acids Research* **37** (2):482.Bivalacqua *et al.* (2002) Comparative responses to alpha,beta-methylene-ATP in cat pulmonary, mesenteric, and hindquarter vascular beds. *Appl. Physiol.* **93** (4):1287.Trezise (1994) The use of antagonists to characterize the receptors mediating depolarization of the rat isolated vagus nerve by  $\alpha,\beta$ -methylene adenosine 5'-triphosphate. *Br. J. Pharmacol.* **112**:282.Bo (1995) Characterization and autoradiographic localization of [<sup>3</sup>H] $\alpha,\beta$ -methylene adenosine 5'-triphosphate binding sites in human urinary bladder. *Br. J. Urol.* **76**:297.