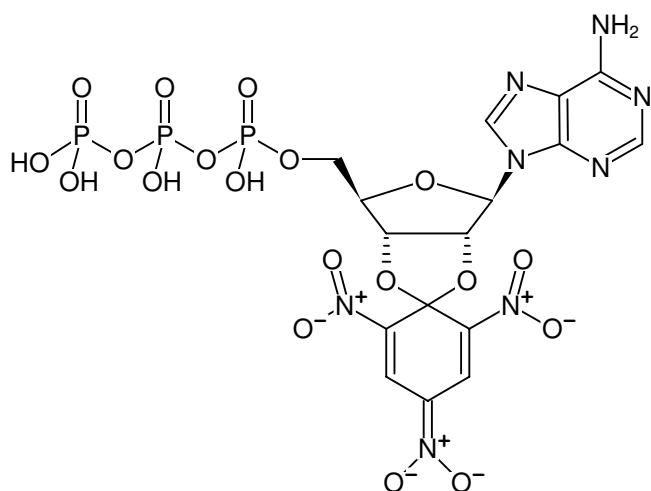


**TNP-ATP**

2',3'-O-Trinitrophenyl-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-221S	200 µl (10 mM)
NU-221L	5 x 200 µl (10 mM)



Structural formula of TNP-ATP

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:** C₁₆H₁₆N₈O₁₉P₃**Molecular Weight:** 718.27 g/mol**Exact Mass:** 717.98 g/mol**CAS#:** 120360-48-7**Purity:** ≥ 95 % (HPLC)**Form:** orange solution in water**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 259/408/470 nm,
 ε 25.0/26.4/18.5 L mol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 408/470 nm,
 λ_{em} 552 nm

Applications:

Agonistic ligand, mainly for nucleoside receptor A₁
 Nucleoside-triphosphates can be converted by different membrane-bound phosphatases into nucleosides acting as nucleoside receptor ligands. The ester form is protected during uptake and transport and can be well-directed released through activation.

Specific Ligands:Ligand for purinergic receptors:P2X₁, P2X₂^[1]Antagonist for purinergic receptors:P2X₁^[2,3], P2X₃^[3,6], P2X₄^[4], P2X₅^[5,7]**Selected References:**

- [1] Linan-Rico *et al.* (2015) Neuropharmacology of purinergic receptors in human sub- mucous plexus: Involvement of P2X₁, P2X₂, P2X₃ channels, P2Y and A₃ metabotropic receptors in neurotransmission. *Neuropharmacology* **95**:83.
- [2] Kur *et al.* (2014) Purinergic control of vascular tone in retina. *J. Physiol.* **592** (3):491.
- [3] Alkayed *et al.* (2012) P2Y₁₁ purinoceptor mediates the ATP-enhanced chemotactic response of neutrophils. *J. Pharmacol. Sci.* **120** (4):288.
- [4] Manohar *et al.* (2012) ATP release and autocrine signaling Through P2X₄ receptors regulate ?? T cell activation. *J. Leukoc. Biol.* **92** (4):787.
- [5] Kamai *et al.* (2006) Involvement of ionotropic purinergic receptors in the histamine-induced enhancement of the cough reflex sensitivity in guinea pigs. *Eur. J. Pharmacol.* **547** (1-3):160.
- [6] Jarvis *et al.* (2001) Modulation of BzATP and formalin induced nociception: attenuation by the P2X receptor antagonist, TNP-ATP and enhancement by the P2X₃ allosteric modulator, cibacon blue. *Br. J. Pharmacol.* **132**:259.
- [7] Ruan *et al.* (2004) Identification of P2X receptors in cultured mouse and rat parasympathetic otic ganglion neurones including P2X knockout studies. *Neuropharmacology* **46**:1039.
- Seifert *et al.* (2012) Inhibitors of membranous adenylyl cyclases. *Trends Pharmacol. Sci.* **33** (2):64.
- Adina-Zada *et al.* (2011) Probing the allosteric activation of pyruvate carboxylase using 2',3'-O- (2,4,6-trinitrophenyl) adenosine 5'-triphosphate as a fluorescent mimic of the allosteric activator acetyl CoA. *Arch. Biochem. Biophys.* **509** (2):117.
- 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.

**TNP-ATP**

2',3'-O-Trinitrophenyl-adenosine-5'-triphosphate, Triethylammonium salt

Goettle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.

Sprang *et al.* (2006) Broad Specificity of Mammalian Adenylyl Cyclase for Interaction with 2',3'-Substituted Purine- and Pyrimidine Nucleotide Inhibitors. *Mol. Pharmacol.* **70**:878.

Berman *et al.* (2003) Interaction of an aromatic dibromoisothiuronium derivative with the Ca (2+)-ATPase of skeletal muscle sarcoplasmic reticulum. *Biochemistry* **42**:3556.

Milgrom *et al.* (1998) Bi-site activation occurs with the native and nucleotide-depleted mitochondrial F1-ATPase. *Biochem J.* **330**:1037.

Faller *et al.* (1990) Binding of the fluorescent substrate analogue 2',3'-O-(2,4,6-trinitrophenylcyclohexadienylidene)-adenosine 5'-triphosphate to the gastric H⁺,K⁺ (+)-ATPase: evidence for cofactor-induced conformational changes in the enzyme. *Biochemistry* **29**:3179.

Hiratsuka *et al.* (1982) Biological activities and spectroscopic properties of chromophoric and fluorescent analogs of adenine nucleoside and nucleotides, 2',3'-O- (2,4,6-trinitrocyclohexadienylidene) adenosine derivatives. *Biochim Biophys Acta.* **719**:509.