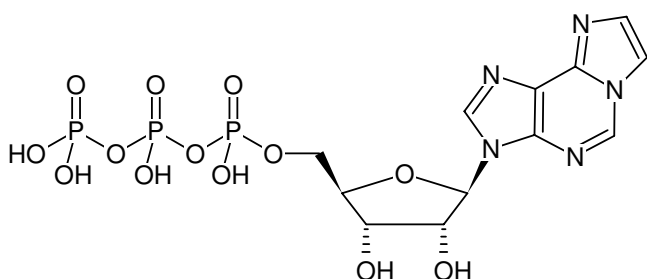


**Etheno-ATP (ϵ -ATP)**(1,N⁶-Etheno-ATP)1,N⁶-Etheno-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1103S	500 μ l (10 mM)
NU-1103L	5 x 500 μ l (10 mM)

Structural formula of Etheno-ATP (ϵ -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:** C₁₂H₁₆N₅O₁₃P₃ (free acid)**Molecular Weight:** 531.20 g/mol (free acid)**Exact Mass:** 531.00 g/mol (free acid)**CAS#:** 60777-99-3**Purity:** \geq 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 275 nm, ϵ 6.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 300 nm, λ_{em} 415 nm**Applications:**Fluorescence quenching^[1]Exchange rate from G-actin and wild type^[2]FRET-studies on conformational changes^[3]Formation in erythrocytes and endothelial cells^[4]Fluorescence anisotropy changes during protein binding^[5]Influence on centromeric binding protein E associated activity^[6]**Specific Ligands:**Yeast and muscle actins^[7]Cofilin and profilin^[8]P2Y receptor in cardiac endothelial cells^[9]**Selected References:**[1] Saratovskikh *et al.* (2010) Kinetic of transport of technogenic oxicans through model membranes. *J. Environmental Protection* **1**:410.[2] McKane *et al.* (2006) Effect of the substitution of muscle actin-specific subdomain 1 and 2 residues in yeast actin and actin function. *J. Biol. Chem.* **281**:29916.[3] Dedova *et al.* (2006) Thymosin β 4 induces a conformational change in actin monomers. *Biophysical J.* **90**:985.[4] Mattig *et al.* (2003) Modulation of adenine nucleotide concentrations in human plasma by erythrocytes and endothelial cells. *Thrombosis Res.* **110**:195.[5] Suarez *et al.* (202) Biochemical defects in retina-specific human ATP binding cassette transporter nucleotide binding domain 1 mutants associated with macular degeneration. *J. Biol. Chem.* **277**:21759.[6] Thrower *et al.* (1995) Mitotic HeLa cells contain a CENP-E associated minus end-directed microtubule motor. *EMBO Journal* **14**:918.[7] Wen *et al.* (2008) Control of the ability of profilin to bind and facilitate nucleotide exchange from G-actin. *J. Biol. Chem.* **283**:9444.[8] Kardos *et al.* (2009) The effects of ADF/cofilin and profilin on the conformation of the ATP-binding cleft of monomeric actin. *Biophysical J.* **96**:2335.[9] Yang *et al.* (1994) Purinergic axis in cardiac blood vessels: Agonist mediated release of ATP from cardiac endothelial cells. *Circ. Res.* **74** (3):401.Singh *et al.* (2011) Crystal structures explain functional differences in the two

**Etheno-ATP (ϵ -ATP)**(1,N⁶-Etheno-ATP)1,N⁶-Etheno-adenosine-5'-triphosphate, Sodium salt

actin depolymerization factors of the malaria parasite. *J. Biol. Chem.* **286** (32): 28256.

Aguilar *et al.* (2001) Ecto enzymatic breakdown of diadenosine polyphosphates by *Xenopus laevis* oocytes. *Eur. J. Biochem.* **268** (5):1289.

Churchich *et al.* (2000) A catalytic site of protein disulfide isomerase probed with adenosine-5'-triphosphate analogs. *BBA-Protein Struct. M.* **1479** (1):293.

Gualix *et al.* (1999) Studies of chromaffin granule functioning by flow cytometry: Transport of fluorescent epsilon-ATP and granular size increase induced by ATP. *Receptor Channel* **6** (6):449.

Franksskiba *et al.* (1994) Quenching of fluorescent nucleotides bound to myosin - a probe of the active-site conformation. *Biochemistry-US* **33** (42):12720.

Miki *et al.* (1994) Domain motion in actin observed by fluorescence resonance energy-transfer. *Biochemistry-US* **33** (33):10171.

Root *et al.* (1992) The accessibility of ethenonucleotides to collisional quenchers and the nucleotide cleft in G-Actin and F-Actin. *Protein Sci.* **1** (8):1014.

Conner *et al.* (1989) Sister chromatid exchange induced by etheno-ATP derivatives invitro. *Cancer Res.* **49** (14):3839.

Wang *et al.* (1981) Exchange of 1, N (6)-etheno-ATP with Actin-bound nucleotides as a tool for studying the steady-state exchange of subunits in F-Actin solutions. *P. Natl. Acad. Sci.-Biol.* **78** (9):5503.

Yanagida (1981) Angles of nucleotides bound to cross-bridges in glycerinated muscle-fiber at various concentrations of epsilon-ATP, epsilon-ADP and epsilon-AMPPNP detected by polarized fluorescence. *J. Mol. Biol.* **146** (4):539.

Burtnick *et al.* (1979) Circular-polarization of the fluorescence of Actin-bound epsilon-ATP - effects of binding DNase-I. *FEBS Lett.* **97** (1):166.

Kaplan *et al.* (1976) Mitochondrial ATPase activity and adenine-nucleotide transport with epsilon-ATP. *J. Cell Biol.* **70** (2):a414.

Hohne *et al.* (1975) New principle for activity measurement of ADP or ATP dependent enzymes - fluorescence quenching of epsilon-ADP and epsilon-ATP by divalent metal-ions. *Anal. Biochem.* **69** (2):607.