

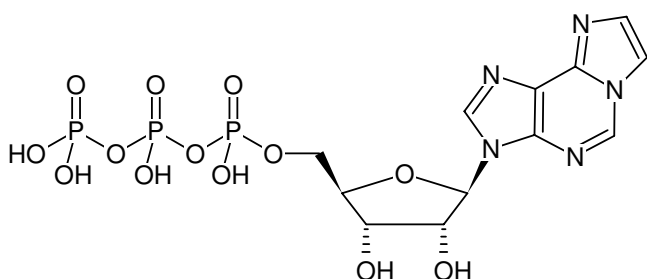


## Etheno-ATP ( $\epsilon$ -ATP)

(1,N<sup>6</sup>-Etheno-ATP)

1,N<sup>6</sup>-Etheno-adenosine-5'-triphosphate, Sodium salt

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



Structural formula of Etheno-ATP ( $\epsilon$ -ATP)

### 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>12</sub>H<sub>16</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (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:**  $\lambda_{max}$  275 nm,  $\epsilon$  6.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5),  $\lambda_{exc}$  300 nm,  $\lambda_{em}$  415 nm

### Applications:

Fluorescence quenching<sup>[1]</sup>

Exchange rate from G-actin and wild type<sup>[2]</sup>

FRET-studies on conformational changes<sup>[3]</sup>

Formation in erythrocytes and endothelial cells<sup>[4]</sup>

Fluorescence anisotropy changes during protein binding<sup>[5]</sup>

Influence on centromeric binding protein E associated activity<sup>[6]</sup>

### Specific Ligands:

Yeast and muscle actins<sup>[7]</sup>

Cofilin and profilin<sup>[8]</sup>

P2Y receptor in cardiac endothelial cells<sup>[9]</sup>

### 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  $\beta$ 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 ( $\epsilon$ -ATP)**(1,N<sup>6</sup>-Etheno-ATP)1,N<sup>6</sup>-Etheno-adenosine-5'-triphosphate, Sodium salt

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

Aguilar *et al.* (2001) Ectoenzymatic 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.