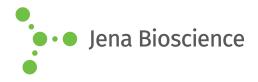
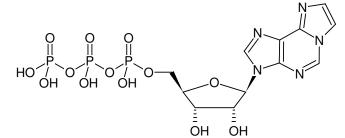
DATA SHEET



Είλου Είλουο-ΑΤΡ (ε-ΑΤΡ)

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

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



Structural formula of Etheno-ATP (ε-ATP)

For general laboratory use.

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: C₁₂H₁₆N₅O₁₃P₃ (free acid)

Molecular Weight: 531.20 g/mol (free acid)

Exact Mass: 531.00 g/mol (free acid)

CAS#: 60777-99-3

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±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]

Coflin and profilin^[8]

P2Y receptor in cardiac endothelial cells^[9]

Selected References:

[1] Saratovskikh *et al.* (2010) Kinetic of transport of technogenic oxicants 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.

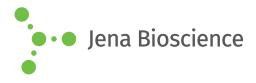
[8] Kardos *et al.* (2009) The effects of ADF/cofilin and profilin on the conformation of the ATP-binding cleft of monomeric actin. *Biophysical J.* **96**:2335.

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









Etheno-ATP (ε-ATP)

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

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.

