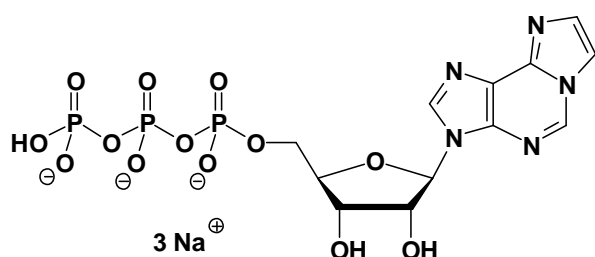


## ε-ATP

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

Cat. No.	Amount
NU-1103S	20 Units
NU-1103L	100 Units



**Cat. No.:** NU-1103

**Molecular Formula:** C<sub>12</sub>H<sub>13</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 528.18 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

#### Spectroscopic properties:

$\lambda_{\max}$  275 nm;  $\epsilon$  6000;  $\lambda_{\text{exc}}$  300 nm;  $\lambda_{\text{em}}$  415 nm

#### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

#### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution

#### Selected References:

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** (1-2):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.