

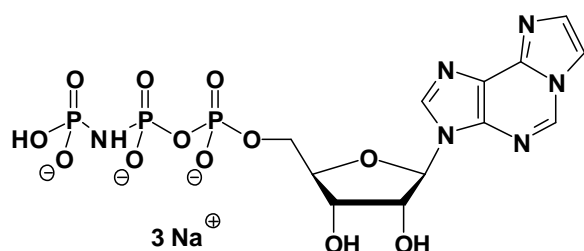
## $\epsilon$ -AppNHp ( $\epsilon$ -AMPPNP)

1,N<sup>6</sup>-Etheno-adenosine-5'-[( $\beta,\gamma$ )-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-1104S	5 Units
NU-1104L	25 Units

### Selected References:

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.



**Cat. No.:** NU-1104

**Molecular Formula:** C<sub>12</sub>H<sub>14</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 527.19 (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 6 months after date of delivery.

### For research use only!

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