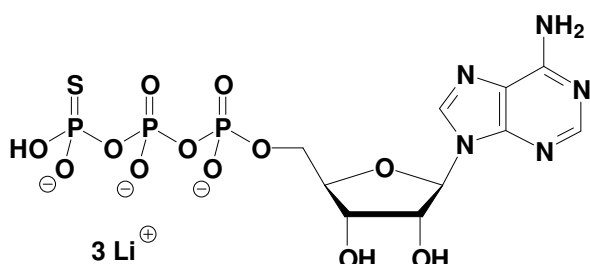


## ATP $\gamma$ S, lyophilized

Adenosine-5'-( $\gamma$ -thio)-triphosphate, Lithium salt

Cat. No.	Amount
NU-406-25	25 mg
NU-406-50	50 mg



**Cat. No.:** NU-406

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub>S (Anion)

**Molecular Weight:** 520.22 (Anion)

**Purity:** > 90%, lyophilized powder

### 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!**

### Selected References:

Menendez-Arias *et al.* (2005) Suppression of Multidrug-resistant HIV-1 Reverse Transcriptase Primer Unblocking Activity by  $\alpha$ -Phosphate-modified Thymidine Analogues. *J. Mol. Biol.* **349** (3):451

Yengo *et al.* (2001) Myosin V is weakly bound to actin in the presence of ATP $\gamma$ S. *Biophys. J.* **80** (1):2609, Part 2.

Buchbinder *et al.* (1993) Stereochemistry of metal-ion coordination to the terminal thiophosphoryl group of adenosine 5'- $\alpha$ -(3-thiotriphosphate) at the active-site of pyruvate-kinase. *Biochemistry-US* **32** (51):14111.

Wang *et al.* (1993) Mg(2+) and ATP or adenosine 5'-[gamma-thio]-triphosphate (ATP-gamma-S) enhances intrinsic fluorescence and induces aggregation which increases the activity of spinach rubisco activase. *Biochim. Biophys. Acta* **1202** (1):47.

Schwer *et al.* (1992) A conformational rearrangement in the spliceosome is dependent on prp16 and atp hydrolysis. *EMBO J.* **11** (13):5033.

Shrimankar *et al.* (1992) Purification and characterization of a mutant dnab protein specifically defective in ATP hydrolysis. *J. Bacteriol.* **174** (23):7689.

Chidiac *et al.* (1992) Effects of adenylnucleotides and carbachol on cooperative interactions among G-proteins. *Biochemistry-US* **31** (44):10908.