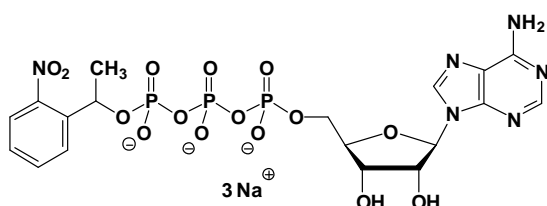


## NPE-caged-ATP

Adenosine-5'-triphosphate, P<sup>3</sup>-(1-(2-nitrophenyl)-ethyl)-ester, Sodium salt

Cat. No.	Amount
NU-301S	150 Units
NU-301L	750 Units



**Cat. No.:** NU-301

**Molecular Formula:** C<sub>18</sub>H<sub>20</sub>N<sub>6</sub>O<sub>15</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 653.30 (Anion)

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

### 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. Store in a dark place and protect from light!

### For research use only!

1 unit = 1 µl of a 10 mM solution

### Selected References:

Zscherp *et al.* (2001) Reaction-induced infrared difference spectroscopy for the study of protein reaction mechanisms. *Biochemistry* **40**:1875.

Scheirlinckx *et al.* (2001) Monitoring of secondary and tertiary structure changes in the gastric H<sup>+</sup>/K<sup>+</sup>-ATPase by infrared spectroscopy *Eur. J. Biochem.* **268 (13)**:3644.

Barth *et al.* (2000) Substrate binding and enzyme function investigated by infrared spectroscopy. *FEBS Lett.* **477**:151.

Hess (1999) Light-Activated (Caged) Biological Ligands. *Encyclopedia of Molecular Biology*, Vol. 3, T.E. Creighton (Ed.), pp. 1385-1391.

Broustovetsky *et al.* (1997) Biochemical and physical parameters of the electrical currents measured with the ADP/ATP carrier by photolysis of caged ADP and ATP. *Biochemistry* **36**:13865.

Higuchi *et al.* (1997) Kinetics of force generation by single kinesin molecules activated by laser photolysis of caged ATP. *Proc. Natl. Acad. Sci. USA* **94**:4395.

Emoto *et al.* (1995) Tension relaxation induced by pulse photolysis of caged ATP in partially crosslinked fibers from rabbit psoas muscle. *Proc. Natl. Acad. Sci. USA* **92**:1461.

Barth *et al.* (1995) Photochemical Release of ATP from "Caged ATP" Studied by Time-Resolved Infrared Spectroscopy. *J. Am. Chem. Soc.* **117**:10311.

Hyman *et al.* (1992) Microtubule-motor activity of a yeast centromere-binding protein complex. *Nature* **359**:533.

Fajer *et al.* (1990) Myosin heads have a broad orientational distribution during isometric muscle contraction: time-resolved EPR studies using caged ATP. *Proc. Natl. Acad. Sci. USA* **87**:5538.

Walker *et al.* (1988) Photolabile 1-(2-Nitrophenyl)ethyl Phosphate Esters of Adenine Nucleotide Analogues. Synthesis and Mechanism of Photolysis. *J. Am. Chem. Soc.* **110**:7170.

Goldman *et al.* (1982) Relaxation of muscle-fibers by photolysis of caged ATP. *Nature* **300**:701.