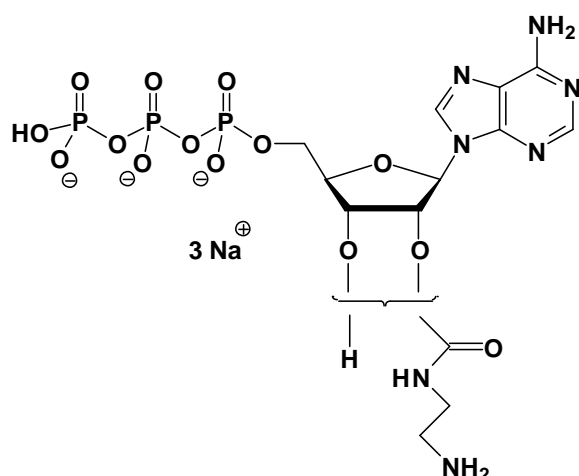


EDA-ATP

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-808S	30 Units
NU-808L	150 Units



Cat. No.: NU-808

Molecular Formula: C₁₃H₁₉N₇O₁₄P₃ (Anion)

Molecular Weight: 590.25 (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.

For research use only!

1 unit = 1 µl of a 10 mM solution

Selected References:

Oiwa *et al.* (2000) Comparative Single-Molecule and Ensemble Myosin Enzymology: Sulfoindocyanine ATP and ADP Derivatives. *Biophys. J.* **78**:3048.

Jameson *et al.* (1997) Fluorescent analogs: Synthesis and Applications. *Methods in Enzymology* **278**:363.

Eccleston *et al.* (1996) Ribose-linked sulfoindocyanine conjugates of ATP: Cy3-EDA-ATP and Cy5-EDA-ATP. *Biophys. J.* **70** (2): MPO29 Part 2.

Oiwa *et al.* (1996) Microscopic observations of Cy3-EDA-ATP and Cy5-EDA-ATP binding to myosin filaments in vitro. *Biophys. J.* **70** (2):MPO30 Part 2.

Conibear *et al.* (1996) Measurement of nucleotide exchange kinetics with isolated synthetic myosin filaments using flash photolysis. *FEBS Lett.* **380** (1-2):13.

Conibear *et al.* (1996) Kinetic and spectroscopic characterization of fluorescent ribose-modified ATP analogs upon interaction with skeletal muscle Myosin subfragment 1. *Biochemistry* **35** (7):2299.

Watson *et al.* (1995) Macromolecular arrangement in the aminoacyl-transfer-RNA-elongation factor-gtp ternary complex - a fluorescence energy-transfer study. *Biochemistry* **34** (24):7904.

Braxton *et al.* (1988) The synthesis of a novel class of ribose-modified nucleotide analogs. 1. affinity purification of skeletal Myosin subfragment-1. *Biophys. J.* **53** (2):A178 Part 2.