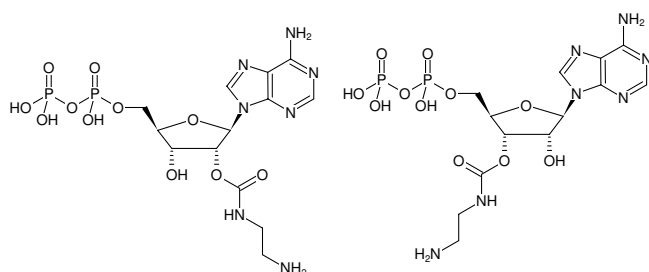


**EDA-ADP**

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

Cat. No.	Amount
NU-802S	30 µl (10 mM)
NU-802L	5 x 30 µl (10 mM)



Structural formula of EDA-ADP

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₁₃H₂₁N₇O₁₁P₂ (free acid)**Molecular Weight:** 513.30 g/mol (free acid)**Exact Mass:** 513.08 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**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.Oiwa *et al.* (1997) Microscopic kinetic measurements of single Cy3-EDA-ADP molecules interacting with Myosin filaments in vitro. *Biophys. J.* **72** (2):TU240 Part 2.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.