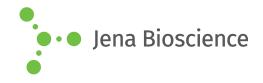
DATA SHEET





■ N⁶-Benzyl-ATP

N⁶-Benzyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1196S	300 μl (10 mM)
NU-1196L	5 x 300 μl (10 mM)

Structural formula of N⁶-Benzyl-ATP

For general laboratory use.

Shipping: shipped on gel packs **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:** C₁₇H₂₂N₅O₁₃P₃ (free acid) **Molecular Weight:** 597.30 g/mol (free acid)

Exact Mass: 597.04 g/mol (free acid)

CAS#: 40922-97-2
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} 269 nm, ϵ 20.5 L mmol⁻¹ cm⁻¹ (Tris-HCl

pH 7.5)

Applications:

Agonistic ligand, mainly for nucleoside receptor A_1 , with less affinity to $A_{2\Delta}$ and A_3

Nucleoside-triphosphates can be converted by different membranebound phosphatases into nucleosides acting as nucleoside receptor ligands. In some cases nucleoside phosphates act also directly on nucleoside receptors.

Selected References:

Sirci et al. (2012) Ligand-, structure- and pharmacophore-based molecular fingerprints: a case study on adenosine A1, A2A, A2B, and A3 receptor antagonists. *J. Comput. Aided Mol. Des.* **26**:1247.

Volonte *et al.* (2009) Membrane components and purinergic signalling: the purinome, a complex interplay among ligands, degrading enzymes, receptors and transporters. *FEBS J.* **276**:318.

Yegutkin (2008) Nucleotide and nucleoside converting enzymes: Important modulators of purinergic signalling cascade. *Biochim. Biophys. Acta* 1783:673.

Joshi et al. (2005) Purine derivatives as ligands for A3 adenosine receptors. Current Topics in Medicinal Chemistry **5**:1275.

Zhou et al. (2005) High affinity ATP/ADP analogues as new tools for studying CFTR gating. J. Physiol. **569 (2)**:447.

Hess (2001) Recent advantages in adenosine receptor antagonist research. Expert Opin. Ther. Patents 11 (10):1533.

Jacobson (2001) Probing adenosine and P2 receptors: design of novel purines and nonpurines as selective ligands. *Drug Development Res.* **52**:178.

Jacobson et al. (2001) Ribose modified nucleosides and nucleotides as ligands for purine receptors. Nucleosides, Nucleotides & Nucleic Acids 20 (4):333.

Gillespie et al. (1999) Engineering of the myosin-ibeta nucleotide-binding pocket to create selective sensitivity to N (6)-modified ADP analogs. *J. Biol. Chem.* **274 (44)**:31373.

Van Galen et al. (1994) A binding site model and structure-activity relationships for rat A3 adenosine receptor. Molecular Pharmacology 45:1101.