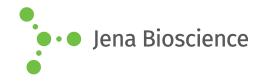
DATA SHEET

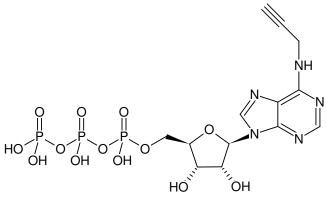




■ N⁶-Propargyl-ATP (N⁶pATP)

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

Cat. No.	Amount
CLK-NU-001-1	1 mg
CLK-NU-001-5	5 mg



Structural formula of N⁶-Propargyl-ATP (N⁶pATP)

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: C₁₃H₁₈N₅O₁₃P₃ (free acid)
Molecular Weight: 545.23 g/mol (free acid)
Exact Mass: 545.01 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Solubility: 10 mM Tris-HCl pH 7.5

Spectroscopic Properties: λ_{max} 262 nm, ϵ 18.0 L mmol⁻¹ cm⁻¹ (Tris-HCl

pH 7.5)

Applications:

in vitro AMPylation of proteins^[1,2]

in vitro polyadenylation of RNA^[3]

The resulting alkyne-functionalized protein $^{[1,2]}$ or RNA $^{[3]}$ can subsequently be processed via Cu(I)-catalyzed (azide-alkyne) click chemistry that offers the choice

- to introduce a Biotin group for subsequent purification tasks (via Azides of Biotin)
- to introduce fluorescent group for subsequent microscopic imaging (via Azides of fluorescent dyes)
- to crosslink the RNA to azide-functionalized biomolecules e.g.proteins

Presolski *et al.*^[4] and Hong *et al.*^[5] provide a general protocol for Cu(I)-catalyzed click chemistry reactions that may be used as a starting point for the set up and optimization of individual assays.

Agonistic ligand, mainly for nucleoside receptor A₁

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.

Please note: This compound contains a phosphoramide linkage which is hydrolyzed at pH <7.0.

For preparation of a 10 mM solution use 100 mM buffer (for example: bicarbonate, borate, phosphate and Tris) to prevent degradation at acidic pH.

Related Products:

Copper (II)-Sulphate (CuSO₄), #CLK-MI004 Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010 Sodium Ascorbate (Na-Ascorbate), #CLK-MI005

Selected References:

[1] Grammel et al. (2011) A Chemical Reporter for Protein AMPylation. J. Am. Chem. Soc. 133:17103.

[2] Broncel *et al.* (2012) A New Chemical Handle for Protein AMPylation at the Host-Pathogen Interface. *ChemBioChem* **13**:183.

[3] Grammel et al. (2012) Chemical Reporter for Monitoring RNA Synthesis and Poly (A) Tail Dynamics. ChemBioChem 13:1112.

[4] Presolski *et al.* (2011) Copper-Catalyzed Azide-Alkyne Click Chemistry for Bioconjugation. *Current Protocols in Chemical Biology* **3**:153.

[5] Hong et al. (2011) Analysis and Optimization of Copper-Catalyzed Azide-Alkyne Cycloaddition for Bioconjugation. Angew. Chem. Int. Ed. 48:9879.

Sirci et al. (2012) Ligand-, structure- and pharmacophore-based molecular fingerprints: a case study on adenosine A1, A2A, A2B, and A3 receptor

DATA SHEET





■ N⁶-Propargyl-ATP (N⁶pATP)

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

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.

Volpini et al. (2002) N6-Alkyl-2-alkynyl derivatives of adenosine as potent and selective agonists of the human adenosine A3 receptor and starting point for searching A2B ligands. J. Med. Chem. 45 (15):3271.

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.

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