

# N6-Propargyl-ATP (N6pATP) (5 mg)

## Ordering Information

[Order Online »](#)

JBS-CLK-NU-001-5	5mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N6-Propargyl-adenosine-5'-triphosphate, Sodium salt

**Appearance** white to off-white solid

**Application Notes** in vitro AMPylation of proteins, in vitro polyadenylation of RNA  
The resulting alkyne-functionalized protein or RNA can subsequently be processed via Cu(I)-catalyzed (azide-alkyne) click chemistry that offers the choiceto 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

**Couple Type** Modified nucleotides

**Formula**  $C_{13}H_{18}N_5O_{13}P_3$  (free acid)

**MW** 545.23 (free acid)

**Purity** ≥ 95 % (HPLC)

**Solubility** 10 mM Tris-HCl pH 7.5

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)