

N6-Propargyl-ATP (N6pATP) (1 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-NU-001-1	1mg
------------------	-----

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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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