

# 6-1MeBu-ATP

Specific triphosphate-source for chemically-genetically engineered protein kinases with selective sensitivity to N<sup>6</sup>-modified ATP analogs.

*BLG-M027-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

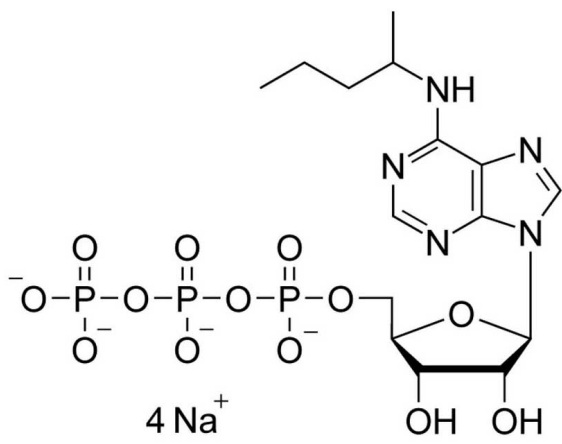
## Ordering Information

[Order Online »](#)

BLG-M027-05	5µmol
BLG-M027-25	5x5µmol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name N6-(1-Methylbutyl)-ATP. tetrasodium salt

Appearance 10 mM aqueous solution

CAS 252889-13-7

Couple Type Modified nucleotides

Formula  $C_{15}H_{26}N_5O_{13}P_3$  (free acid)

MW 577.3

Purity > 95% HPLC

Quantity 5µmol≈2.9mg

Solubility Soluble in water or buffer.

**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)