

# 3-Methyladenine

## PI3 kinase inhibitor

PI3 kinase inhibitor and a well-known inhibitor of autophagy. This compound has been shown to induce autophagy under nutrient-rich conditions as well as inhibiting starvation-induced autophagy due to differential effects on class I versus class III PI3 kinase.

Citations: 13

[View Online »](#)

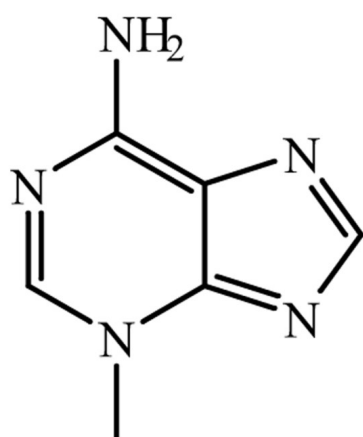
## Ordering Information

[Order Online »](#)

BML-AP502-0025	25mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions can be stored at -20°C for up to 3 months.
Handling	Solutions can be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	6-Amino-3-methylpurine
Appearance	White solid.
CAS	5142-23-4
Couple Target	PI3 kinase
Couple Type	Inhibitor
Formula	$C_6H_7N_5$
Identity	Determined by NMR.
MW	149.2
Purity	≥98% (HPLC)
RTECS	AU6520000
Solubility	Soluble in DMSO (10 mg/ml) or 100% ethanol (5 mg/ml, warm).



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-  
fr@enzolifesciences.com](mailto:info-fr@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)