

3-Methyladenine

Autophagy Inhibitor; PI3K inhibitor

Catalog code: inh-3ma-2

<https://www.invivogen.com/3-methyladenine>

For research use only

Version 23D21-MM

PRODUCT INFORMATION

Contents

- 2 x 50 mg of 3-Methyladenine

Storage and stability:

- 3-Methyladenine is shipped at room temperature. Upon receipt, store at 15-25°C (room temperature).
- Upon resuspension in DMSO, prepare aliquots of 3-Methyladenine and store at -20°C. 3-Methyladenine resuspended in DMSO is stable for 3 months when properly stored. Avoid repeated freeze-thaw cycles.

Quality control:

- ≥90% (UHPLC).
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

3-Methyladenine (also known as 3-MA) is an inhibitor of phosphatidylinositol 3-kinases (PI3K). PI3K plays an important role in many biological processes, including controlling the activation of mTOR, a key regulator of autophagy. 3-MA inhibits autophagy by blocking autophagosome formation via the inhibition of class III PI3K¹. 3-Methyladenine plays a dual role in autophagy. Prolonged treatment with 3-Methyladenine promotes autophagy under nutrient-rich conditions, whereas 3-Methyladenine inhibits starvation-induced autophagy¹. These results are attributed to differential effects on class I versus class III PI3K.

In addition to its role in autophagy, 3-Methyladenine has been implicated in cancer therapy². It has been revealed that 3-Methyladenine suppresses the invasion of highly metastatic cancer cells through the inhibition of class I and II PI3K³. Further studies demonstrated that 3-Methyladenine can induce caspase-dependent cell death that is independent of autophagy inhibition⁴.

1. **Wu Y. et al., 2010.** Dual role of 3-methyladenine in modulation of autophagy via different temporal patterns of inhibition on class I and III phosphoinositide 3-kinase. *J Biol Chem.* 285(14):10850-61. 2. **Cheong H et al., 2012.** Therapeutic targets in cancer cell metabolism and autophagy. *Nature Biotechnology* 30:671-678. 3. **Ito S. et al., 2007.** 3-Methyladenine suppresses cell migration and invasion of HT1080 fibrosarcoma cells through inhibiting phosphoinositide 3-kinases independently of autophagy inhibition. *Int J Oncol* 31 261-268. 4. **Hou H. et al., 2012.** Inhibitors of phosphatidylinositol 3'-kinases promote mitotic cell death in HeLa cells. *PLoS One.* 7(4):e35665. 5. **Jing C. et al., 2012.** Autophagy activation is associated with neuroprotection against apoptosis via a mitochondrial pathway in a rat model of subarachnoid hemorrhage. *Neuroscience.* 213:144-53.

CHEMICAL PROPERTIES

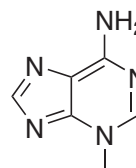
CAS number: 5142-23-4

Formula: C₆H₇N₅

Molecular weight: 149.2 g/mol

Solubility: 7.5 mg/ml (50 mM) in DMSO heated for 5 minutes at 55°C

Structure:



METHODS

Preparation of stock solution at 7.5 mg/ml (50 mM)

1. Weigh 5 mg of 3-Methyladenine.
2. Add 670 µl of DMSO.
3. Heat to 55°C for 5 minutes.
4. Once 3-Methyladenine has been dissolved in DMSO, it can be used immediately or stored at -20°C for 3 months.
5. Prepare further dilutions using sterile culture medium. We do not recommend storing the aqueous solution for more than one day.

PROTOCOLS

For reference only; as described in the indicated publications.

Cell Culture Assay¹

Cell type: Mouse embryonic fibroblasts (MEFs)

Working concentration: 5 mM

Pre-incubation time: 1-9 hours

Method: Confocal microscopy observation of LC3 autophagy

Animal Study⁵

Animal model: Adult male Sprague-Dawley rats

Dose: 400 nM

Administration: Intracerebral ventricular

Solubility: 30% propylene glycol, 5% Tween 80, 65% D5W, 30 mg/ml

RELATED PRODUCTS

Product	Description	Cat. Code
Rapamycin	mTOR inhibitor	tlrl-rap
Bafilomycin A1	Autophagy inhibitor	tlrl-baf1

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-34-80

E-mail: info@invivogen.com