

pp242

mTOR inhibitor

Catalog # inh-pp242

For research use only

Version # 16113-MM

PRODUCT INFORMATION

Contents:

- 5 mg pp242

Storage and stability:

- pp242 is provided lyophilized and shipped at room temperature. Store at -20 °C. Lyophilized pp242 is stable for 2 years when properly stored.
- Upon resuspension, prepare aliquots of pp242 and store at -20 °C. Resuspended pp242 is stable for 6 months when properly stored.

Quality control:

- Purity $\geq 97\%$ (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) is confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

pp242 (also known as Torkinib) is a potent, selective and ATP-competitive inhibitor of mTOR (mammalian target of rapamycin) kinase, a key protein kinase of the PI3K/AKT signaling pathway that plays a central role in controlling cell growth, proliferation, and survival. mTOR is a major negative regulatory axis of autophagy. Thus, pp242 is an effective inducer of autophagy¹. Specifically, pp242 is able to induce autophagy by inhibiting the two mTOR-containing complexes (mTORC1 and mTORC2), neither of which is fully blocked by classical mTOR inhibitors such as rapamycin^{2,4}.

1. Zhou J. et al., 2013. Activation of lysosomal function in the course of autophagy via mTORC1 suppression and autophagosome-lysosome fusion. *Cell Res.* 23(4):508-23.
2. Feldman, ME. et al., 2009. Active-site inhibitors of mTOR target rapamycin-resistant outputs of mTORC1 and mTORC2. *PLoS Biol.* 7:371-83.
3. Shao H. et al., 2009. Dual targeting of mTORC1/C2 complexes enhances histone deacetylase inhibitor-mediated anti-tumor efficacy in primary HCC cancer in vitro and in vivo. *J. Hepatol.* 56(1):176-83.
4. Janes MR. et al., 2010. Effective and selective targeting of Ph+ leukemia cells using a TORC1/2 kinase inhibitor. *Nat Med.* 16(2):205-213.

CHEMICAL PROPERTIES

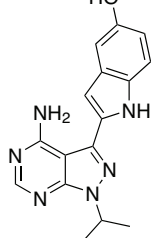
Solubility: 62 mg/ml (201.1 mM) in DMSO
18 mg/ml (58.4 mM) in ethanol

CAS number: 1092351-67-1

Formula: C₁₆H₁₆N₆O

Molecular weight: 308.3

Structure:



METHODS

Preparation of 20 mM (6.2 mg/ml) stock solution

- Add 810 μ l of DMSO to 5 mg pp242. Mix by vortexing.
- Prepare further dilutions using sterile, endotoxin-free water or an aqueous buffer.

Working concentration: 250 nM (77.1 ng/ml) to 1 μ M (308.3 ng/ml) for cell culture assays

Autophagy reporter assay:

Described below is a protocol to study the effects of pp242 in HeLa-Difluo™ hLC3b cells, autophagy reporter cells derived from the human epithelial carcinoma HeLa cell line. HeLa-Difluo™ hLC3b cells express two fluorescent reporter genes (RFP and GFP) fused to the N-terminal of the LC3 protein. The expression of this fusion protein enables the monitoring of autophagic flux in real time.

Day 1

1. Prepare a HeLa-Difluo™ hLC3b cell suspension at $\sim 100,000$ cells/ml.
2. Add 500 μ l of cell suspension ($\sim 50,000$ cells) per well of a 24-well plate.
3. Leave to incubate overnight at 37 °C in a 5% CO₂ incubator.

Day 2

1. Add 20 μ l of pp242 at a final concentration of 250 nM to 1 μ M and incubate at 37 °C for 30 min to 1 hour.
2. Perform image-based analysis for autophagy using a fluorescent microscope.

PROTOCOLS

For reference only; as described in the indicated publications.

Cell Culture Assay¹

Cells: MEF (Murine Embryonic Fibroblasts) expressing GFP-LC3

Working concentration: 1 μ M

Incubation time: 3 hours

Method: Immunoblotting and GFP-LC3 puncta

Cell Culture Assay⁴

Cells: Fibroblasts and solid tumor cell lines (SKOV3, PC-3, 786-O, U87)

Working concentration: 50-400 nM

Incubation time: 1.5-3 hours

Method: Immunoblotting and immunofluorescence analysis

Animal Study⁴

Animal model: Balb/cJ mice

Dose: 60 mg/kg daily for 9 days

Administration: Oral gavage

RELATED PRODUCTS

Product	Description	Cat. Code
HeLa-Difluo™ hLC3 cells	Autophagy reporter cells	heldf-hlc3b
Rapamycin	mTOR inhibitor	trl1-rap
RAW-Difluo™ hLC3 cells	Autophagy reporter cells	rawd-f-mlc3b
Torin 1	mTOR inhibitor	inh-tor1

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 Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com