

Torin 1

mTOR inhibitor

Catalog code: inh-tor1, inh-tor1-5

<https://www.invivogen.com/torin1>

For research use only

Version 23L08-MM

PRODUCT INFORMATION

Contents Torin 1 is available in two quantities:

- **inh-tor1:** 10 mg Torin 1
- **inh-tor1-5:** 5 x 10 mg Torin 1

Storage and stability

- Torin 1 is shipped at room temperature. Upon receipt, store at -20 °C.

- Upon resuspension, prepare aliquots of Torin 1 and store at -20 °C. Resuspended Torin 1 is stable for 6 months when properly stored. Avoid repeated freeze-thaw cycles.

Quality control

- Purity ≥97% (UHPLC)

- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) is confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Torin 1 is a potent and selective ATP-competitive inhibitor of mTOR (mammalian target of rapamycin) kinase¹, the catalytic subunit of two functionally distinct complexes (mTORc1 and mTORc2) that coordinately promote cell growth, proliferation, and survival. Unlike classical mTOR inhibitors such as rapamycin, Torin 1 is able to effectively block phosphorylation of mTORc1 and mTORc2². Torin 1 is an effective inducer of autophagy, as inhibition of mTOR mimics cellular starvation by blocking signals required for cell growth and proliferation^{3,4}. Interestingly, Torin 1 has also been shown to have cytotoxic and cytostatic effects in rapamycin-resistant cells⁵.

1. Liu Q. *et al.*, 2010. Discovery of 1-(4-(4-propionylpiperazin-1-yl)-3-(trifluoromethyl) phenyl)-9-(quinolin-3-yl)benzo[h][1,6]naphthyridin-2(1H)-one as a highly potent, selective Mammalian Target of Rapamycin (mTOR) inhibitor for the treatment of cancer. *Med Chem.* 53(19):7146-55. 2. Thoreen C.C. *et al.*, 2009. An ATP-competitive mammalian target of rapamycin inhibitor reveals rapamycin-resistant functions of mTORC1. *J Biol Chem.* 284(12):8023-32. 3. Jung C.H. *et al.*, 2010. mTOR regulation of autophagy. *FEBS Lett.* 584(7):1287-95. 4. Wang, R.C. *et al.*, 2012. Akt-mediated regulation of autophagy and tumorigenesis through Beclin 1 phosphorylation. *Science* 338(6109):956-9. 5. Park S. *et al.*, 2016. Rapamycin-resistant and torin-sensitive mTOR signaling promotes the survival and proliferation of leukemic cells. *BMB Rep.* 49(1):63-68. 6. 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.

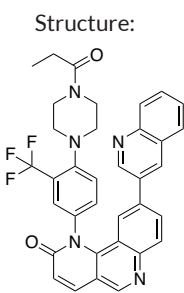
CHEMICAL PROPERTIES

Solubility: 2 mg/ml (3.29 mM) in DMSO

CAS number: 1222998-36-8

Formula: C₃₅H₂₈F₃N₅O₂

Molecular weight: 607.6 g/mol



METHODS

Preparation of 3 mM (1.823 mg/ml) stock solution

1. Weigh 2 mg of Torin 1.
2. Add 1.097 ml of DMSO to the 2 mg of Torin 1. Mix by vortexing.
3. Prepare dilutions using sterile, endotoxin-free water or aqueous buffer.

Working concentration: 250 nM (152 ng/ml) to 1 μM (607.6 ng/ml) for cell culture assays

Autophagy reporter assay:

Described below is a protocol to study the effects of Torin 1 in HeLa-Difluo™ hLC3b cells, an autophagy reporter cell line derived from the human epithelial carcinoma HeLa cell line. These 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. For more information, visit www.invivogen.com/hela-difluo-hlc3.

Day 1

1. Prepare a HeLa-Difluo™ hLC3b cell suspension at ~100,000 cells/ml.
2. Add 500 μl of cell suspension 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 μl of Torin 1 at a final concentration of 250 nM to 1 μM and incubate at 37 °C for 30 min to 1 hour.
2. Perform image-based analysis using a fluorescent microscope.

PROTOCOLS

For reference only; as described in the indicated publications.

Cell Culture Assay⁴

Cells: HeLa expressing GFP-LC3

Working concentration: 250 nM

Incubation time: 2-4 hours

Cell Culture Assay⁶

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

Working concentration: 1 μM

Incubation time: 3 hours

Animal Study¹

Animal model: Mice bearing U87MG xenografts

Dose: 20 mg/kg daily for 10 days

Administration: Intraperitoneal (IP)

RELATED PRODUCTS

Product	Description	Cat. Code
HeLa-Difluo™ hLC3 Cells	Autophagy reporter cells	heldf-hlc3b
Rapamycin	mTOR inhibitor	tlrl-rap
RAW-Difluo™ hLC3 Cells	Autophagy reporter cells	rawdf-mlc3b

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 3622-3480

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com