

MHY-1485

mTOR activator

Catalog # F17-901B

Lot # D3370-17

CAS # 326914-06-1

Product Description

Molecular Formula: C₁₇H₂₁N₇O₄

Appearance: off white powder

Molecular Weight: 387.40

Purity: >98% (TLC); NMR (conforms)

Solubilization: May be dissolved in DMSO (20 mg/ml); or DMF (10 mg/ml)

Alias

4,6-Di-4-morpholinyl-N-(4-nitrophenyl)-1,3,5-triazin-2-amine

Specific Activity

Pan-kinase inhibitor

Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to two months.

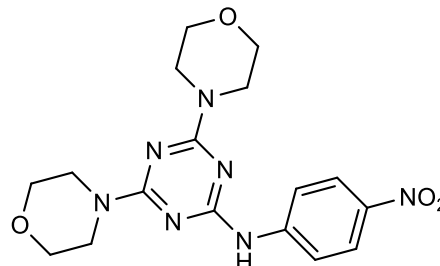
Scientific Background

mTOR activator suppresses autophagy by inhibiting the fusion of autophagosomes with lysosomes leading to the accumulation of enlarged autophagosomes¹. Increases ovarian follicle development². Displays cellular protective effects from cytotoxic doses of dexamethasone³ or UV radiation⁴. A useful tool for probing the cellular role of mTOR⁵. Cell permeable.

References

1. YJ Choy *et al.* *PLoS One* 2012 7(8):e43418
2. Y Cheng *et al.* *PLoS One* 2015 10(2):e0117769
3. S Zhao *et al.* *Biochem. Biophys. Res. Commun.* 2016 481:212
4. B Yang *et al.* *Oncotarget* 2017 481:212 Jan 4, Epub
5. C Li and HM Siragy *Am. J. Physiol. Endocrinol. Metab.* 2015 309:E302

Molecular Structure



MHY-1485

mTOR activator

Catalog #	F17-901B
CAS #	326914-06-1
Lot #	D3370-17
Purity	>98% (HPLC); NMR (conforms)
Format	white powder
Solubilization	May be dissolved in DMSO (20 mg/ml); or DMF (10 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to two months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.