

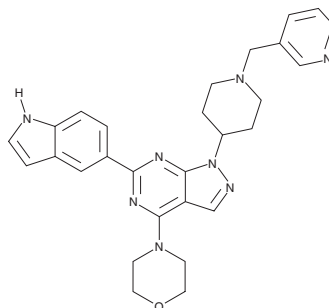
# PRODUCT INFORMATION



## WAY-600

Item No. 29214

**CAS Registry No.:** 1062159-35-6  
**Formal Name:** 6-(1H-indol-5-yl)-4-(4-morpholinyl)-1-[1-(3-pyridinylmethyl)-4-piperidinyl]-1H-pyrazolo[3,4-d]pyrimidine  
**MF:** C<sub>28</sub>H<sub>30</sub>N<sub>8</sub>O  
**FW:** 494.6  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

WAY-600 is supplied as a solid. A stock solution may be made by dissolving the WAY-600 in the solvent of choice, which should be purged with an inert gas. WAY-600 is soluble in the organic DMSO.

### Description

WAY-600 is an inhibitor of mammalian target of rapamycin (mTOR; IC<sub>50</sub> = 0.009 μM).<sup>1</sup> It selectively inhibits mTOR over PI3Kα and PI3Kγ (IC<sub>50</sub>s = 1.96 and 8.45 μM, respectively) and 24 additional lipid and protein kinases in a panel (IC<sub>50</sub>s = >50 μM for all). WAY-600 decreases phosphorylation of the mTORC1 and mTORC2 substrates Akt and S6 kinase (S6K1), respectively, in a cell-free assay when used at concentrations of 0.2, 1, and 5 μM. It decreases proliferation of nine cancer cell lines, including breast, prostate, glioma, kidney, and colorectal cancer cells, with IC<sub>50</sub> values ranging from 0.6 to 2.5 μM. WAY-600 reduces tumor growth in a HepG2 mouse xenograft model when administered at a dose of 10 mg/kg for 14 days.<sup>2</sup>

### References

1. Yu, K., Toral-Barza, L., Shi, C., *et al.* Biochemical, cellular, and *in vivo* activity of novel ATP-competitive and selective inhibitors of the mammalian target of rapamycin. *Cancer Res.* **69(15)**, 6232-6240 (2009).
2. Wang, K., Fan, Y., Chen, S., *et al.* MEK-ERK inhibition potentiates WAY-600-induced anti-cancer efficiency in preclinical hepatocellular carcinoma (HCC) models. *Biochem. Biophys. Res. Commun.* **474(2)**, 330-337 (2016).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/03/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM