# **PRODUCT** INFORMATION



## IWP-O1

Item No. 38418

CAS Registry No.:	2074607-48-8
Formal Name:	5-phenyl-N-(5-phenyl-2-
	pyridinyl)-4-(4-pyridinyl)-1H-
	1,2,3-triazole-1-acetamide
Synonym:	Inhibitor of Wnt Production-O1 $\sqrt{-1}$
MF:	$C_{26}H_{20}N_{6}O$
FW:	432.5
Purity:	≥98%
UV/Vis.:	$\lambda_{max}$ : 268 nm $\int \int \langle \cdot \rangle$
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

IWP-O1 is supplied as a solid. A stock solution may be made by dissolving the IWP-O1 in the solvent of choice, which should be purged with an inert gas. IWP-O1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IWP-O1 in these solvents is approximately 10 mg/ml.

#### Description

IWP-O1 is an inhibitor of porcupine (PORCN), an enzyme that catalyzes palmitoylation of Wnt proteins.<sup>1</sup> It inhibits Wnt signaling in L-Wnt-STF cells (IC<sub>50</sub> = 80 pM in a reporter assay) and phosphorylation of Dishevelled-2 (Dvl2) and Dvl3 in HeLa cells when used at a concentration of 1  $\mu$ M. IWP-O1 (1  $\mu$ M) enhances the cytotoxicity of the glycolysis inhibitors 2-deoxyglucose (2-DG) and lonidamine (Item No. 14640) in CAL 27 oral squamous cell carcinoma (OSCC) cells.<sup>2</sup>

### References

- 1. You, L., Zhang, C., Yarravarapu, N., et al. Development of a triazole class of highly potent Porcn inhibitors. Bioorg. Med. Chem. Lett. 26(24), 5891-5895 (2016).
- 2. Kleszcz, R. and Paluszczak, J. The Wnt signaling pathway inhibitors improve the therapeutic activity of glycolysis modulators against tongue cancer cells. Int. J. Mol. Sci. 23(3), 1248 (2022).

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

#### SAFETY DATA

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.

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, 06/05/2023

## 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