

PRODUCT INFORMATION



IWP-12

Item No. 22729

CAS Registry No.: 688353-45-9
Formal Name: N-(6-methyl-2-benzothiazolyl)-2-
[(3,4,6,7-tetrahydro-3,6-dimethyl-4-
oxothieno[3,2-d]pyrimidin-2-yl)thio]-
acetamide

Synonym: Inhibitor of Wnt Production-12

MF: C₁₈H₁₈N₄O₂S₃

FW: 418.6

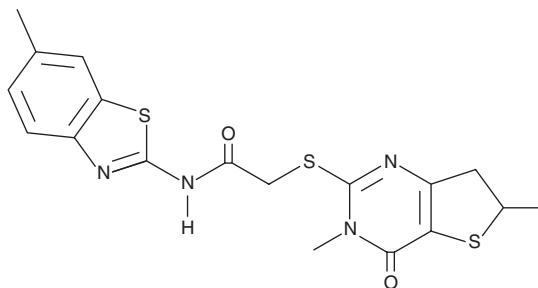
Purity: ≥98%

UV/Vis.: λ_{max}: 278, 302, 335 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

IWP-12 is supplied as a crystalline solid. A stock solution may be made by dissolving the IWP-12 in the solvent of choice. IWP-12 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of IWP-12 in these solvents is approximately 2 and 5 mg/ml, respectively.

IWP-12 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IWP-12 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. IWP-12 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

IWP-12 is a small molecule inhibitor of porcupine (PORCN), which is an acyltransferase that palmitoylates and activates Wnt proteins.¹ IWP-12 inhibits cell-autonomous Wnt signaling with an IC₅₀ value of 15 nM.

Reference

1. Wang, X., Moon, J., Dodge, M.E., *et al.* The development of highly potent inhibitors for porcupine. *J. Med. Chem.* **56**(6), 2700-2704 (2013).

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, 09/12/2017

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