

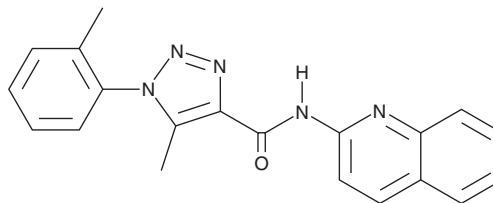
PRODUCT INFORMATION



YW1128

Item No. 27450

CAS Registry No.: 2131223-64-6
Formal Name: 5-methyl-1-(2-methylphenyl)-N-2-quinolinyl-1H-1,2,3-triazole-4-carboxamide
MF: C₂₀H₁₇N₅O
FW: 343.4
Purity: ≥98%
UV/Vis.: λ_{max}: 257, 318, 332 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

YW1128 is supplied as a crystalline solid. A stock solution may be made by dissolving the YW1128 in the solvent of choice, which should be purged with an inert gas. YW1128 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of YW1128 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of YW1128 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of YW1128 in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

YW1128 is an inhibitor of Wnt/β-catenin signaling with an IC₅₀ value of 4.1 nM in a reporter assay.¹ It decreases protein levels of β-catenin in the presence of the GSK3β inhibitor lithium chloride and increases protein levels of Axin1 in HEK293 cells. YW1128 decreases lipid accumulation and the expression of gluconeogenic and lipogenic genes in Huh7 cells. It decreases the hepatic expression of Wnt target genes, improves glucose tolerance, and prevents body weight increases and hepatic lipid accumulation in mice fed a high-fat diet, but not mice fed normal chow, when administered at a dose of 40 mg/kg every other day for 11 weeks.

Reference

1. Obianom, O.N., Ai, Y., Li, Y., et al. Triazole-based inhibitors of the Wnt/β-catenin signaling pathway improve glucose and lipid metabolism in diet-induced obese mice. *J. Med. Chem.* **62**(2), 727-741 (2019).

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, 01/30/2020

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