

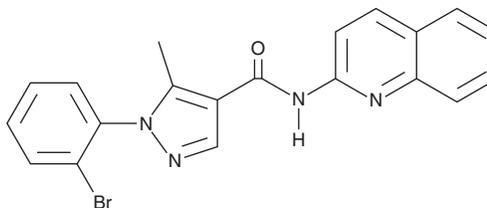
# PRODUCT INFORMATION



**YW2065**

Item No. 42959

**CAS Registry No.:** 2131223-85-1  
**Formal Name:** 1-(2-bromophenyl)-5-methyl-N-2-quinolinyl-1H-pyrazole-4-carboxamide  
**MF:** C<sub>20</sub>H<sub>15</sub>BrN<sub>4</sub>O  
**FW:** 407.3  
**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

YW2065 is supplied as a solid. A stock solution may be made by dissolving the YW2065 in the solvent of choice, which should be purged with an inert gas. YW2065 is slightly soluble (0.1-1 mg/ml) in DMSO.

## Description

YW2065 is an inhibitor of Wnt/ $\beta$ -catenin signaling and an activator of AMP-activated protein kinase (AMPK).<sup>1</sup> It induces degradation of  $\beta$ -catenin and increases the levels of Axin1 and phosphorylated AMPK and acetyl-CoA carboxylase (ACC), which is a downstream substrate of AMPK, in SW620 colorectal cancer cells when used at a concentration of 3  $\mu$ M. YW2065 inhibits recombinant human Wnt3a-induced reporter activation in HEK293 cells (IC<sub>50</sub> = 2.3 nM) and the growth of SW480 colorectal cancer cells (GI<sub>50</sub> = 2.4 nM). *In vivo*, YW2065 (50 mg/kg) decreases tumor volume and weight in an SW620 mouse xenograft model.

## Reference

1. Yang, W., Li, Y., Ai, Y., et al. Pyrazole-4-carboxamide (YW2065): A therapeutic candidate for colorectal cancer via dual activities of Wnt/ $\beta$ -catenin signaling inhibition and AMP-activated protein kinase (AMPK) activation. *J. Med. Chem.* **62**(24), 11151-11164 (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, 07/14/2025

## 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