

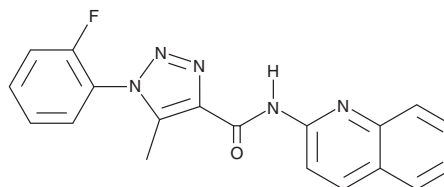
PRODUCT INFORMATION



YW1159

Item No. 27451

CAS Registry No.: 2131223-69-1
Formal Name: 1-(2-fluorophenyl)-5-methyl-N-2-quinolinyl-1H-1,2,3-triazole-4-carboxamide
Synonym: Compound 3d
MF: C₁₉H₁₄FN₅O
FW: 347.3
Purity: ≥98%
UV/Vis.: λ_{max}: 257, 318, 332 nm
Supplied as: A crystalline 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

YW1159 is supplied as a crystalline solid. A stock solution may be made by dissolving the YW1159 in the solvent of choice, which should be purged with an inert gas. YW1159 is soluble in organic solvents such as dimethyl formamide (DMF) and chloroform. The solubility of YW1159 in these solvents is approximately 5 and 30 mg/ml, respectively.

YW1159 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, YW1159 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. YW1159 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

YW1159 is an inhibitor of Wnt signaling with an IC₅₀ value of 1.2 nM in a reporter assay.^{1,2} It is not cytotoxic to HEK293 cells following incubations of up to 24 hours but cell viability is decreased in a concentration-dependent manner after 72 hours.²

References

1. Xue, F. and Shu, Y. Wnt signaling pathway inhibitors for treatments of disease. *University of Maryland, Baltimore*. **WO 2017/151786 A1** (2017).
2. 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, 07/10/2025

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