

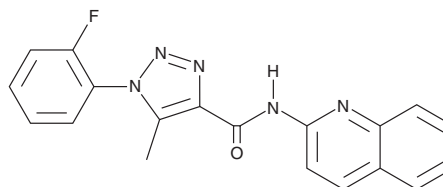
# PRODUCT INFORMATION



**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:** C<sub>19</sub>H<sub>14</sub>FN<sub>5</sub>O  
**FW:** 347.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 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

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 IC<sub>50</sub> value of 1.2 nM in a reporter assay.<sup>1,2</sup> 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.<sup>2</sup>

## 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, 12/05/2019

## 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](http://WWW.CAYMANCHEM.COM)