

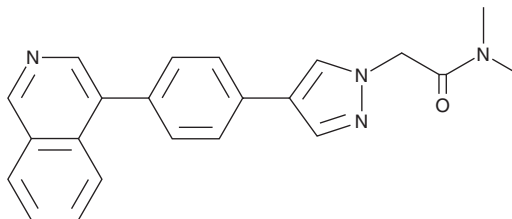
# PRODUCT INFORMATION



**BI-1347**

Item No. 43667

**CAS Registry No.:** 2163056-91-3  
**Formal Name:** 4-[4-(4-isoquinolinyl)phenyl]-N,N-dimethyl-1H-pyrazole-1-acetamide  
**MF:** C<sub>22</sub>H<sub>20</sub>N<sub>4</sub>O  
**FW:** 356.4  
**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

BI-1347 is supplied as a solid. A stock solution may be made by dissolving the BI-1347 in the solvent of choice, which should be purged with an inert gas. BI-1347 is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

## Description

BI-1347 is an inhibitor of Cdk8 and Cdk19 (IC<sub>50</sub>s = 1.4 and 1.7 nM, respectively).<sup>1</sup> It is selective for Cdk8 and Cdk19 over Cdk1, -2, -4, -6, -7, and -9 and a panel of 318 additional kinases at 1 μM. *In vivo*, BI-1347 (10 mg/kg) reduces tumor volume in an MV4-11 mouse xenograft model.

## Reference

1. Hofmann, M.H., Mani, R., Engelhardt, H., *et al.* Selective and potent CDK8/19 inhibitors enhance NK-cell activity and promote tumor surveillance.. *Mol. Cancer Ther.* **19**(4), 1018-1030 (2020).

### 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, 08/08/2025

## 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