

# PRODUCT INFORMATION

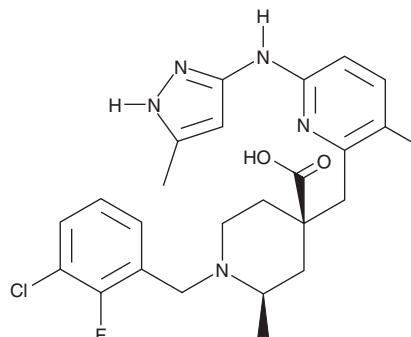
**LY3295668**

Item No. 27914



**CAS Registry No.:** 1919888-06-4  
**Formal Name:** (2R,4R)-1-[(3-chloro-2-fluorophenyl)methyl]-4-[[[3-fluoro-6-[(5-methyl-1H-pyrazol-3-yl)amino]-2-pyridinyl)methyl]-2-methyl-4-piperidinecarboxylic acid

**Synonym:** AK-01  
**MF:** C<sub>24</sub>H<sub>26</sub>ClF<sub>2</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 490.0  
**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

LY3295668 is supplied as a solid. A stock solution may be made by dissolving the LY3295668 in the solvent of choice, which should be purged with an inert gas. LY3295668 is sparingly soluble (1-10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

LY3295668 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

## Description

LY3295668 is an inhibitor of Aurora A kinase (IC<sub>50</sub> = <1 nM).<sup>1</sup> It is selective for Aurora A kinase over a panel of 498 other kinases at 1 μM. LY3295668 (200 nM) induces apoptosis in HeLa cervical cancer cells. It induces cell cycle arrest at the G<sub>2</sub> phase in HeLa cells (IC<sub>50</sub> = 108 nM). *In vivo*, LY3295668 (40 mg/kg per day) decreases tumor volume in a patient-derived xenograft (PDX) mouse model of small cell lung cancer.

## Reference

1. Du, J., Yan, L., Torres, R., *et al.* Aurora A-selective inhibitor LY3295668 leads to dominant mitotic arrest, apoptosis in cancer cells, and shows potent preclinical antitumor efficacy. *Mol. Cancer Ther.* **18**(12), 2207-2219 (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, 10/08/2024

## 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