

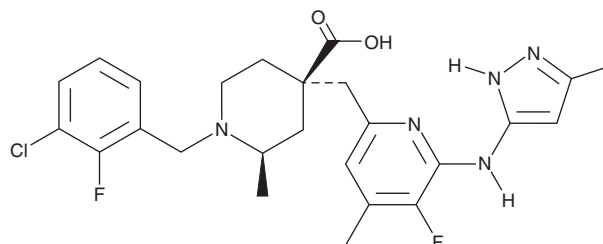
PRODUCT INFORMATION



JAB-2485

Item No. 44291

CAS Registry No.: 2899209-55-1
Formal Name: 1-[(3-chloro-2-fluorophenyl)methyl]-4-[[5-fluoro-4-methyl-6-[(3-methyl-1H-pyrazol-5-yl)amino]-2-pyridinyl)methyl]-2-methyl-4-piperidinecarboxylic acid
MF: C₂₅H₂₈ClF₂N₅O₂
FW: 504.0
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

JAB-2485 is supplied as a solid. A stock solution may be made by dissolving the JAB-2485 in the solvent of choice, which should be purged with an inert gas. JAB-2485 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

JAB-2485 is an inhibitor of Aurora A kinase (IC₅₀ = 0.327 nM).¹ It is selective for Aurora A over Aurora B and Aurora C kinases at 2 μM. JAB-2485 decreases levels of phosphorylated Aurora A kinase in NCI H446 small-cell lung cancer cells (IC₅₀ = 2.08 nM) and reduces the viability of 11 lung cancer cell lines (IC₅₀s = 29.7-982 nM). JAB-2485 (0.3, 1, and 3 μM) induces apoptosis and cell cycle arrest at the G₂/M phase in NCI H446 cells. It decreases tumor volume without affecting body weight in an NCI H446 mouse xenograft model when administered at doses of 10 or 20 mg/kg per day.

Reference

1. Yang, G., Lin, Y., Sun, X., *et al.* Preclinical evaluation of JAB-2485, a potent AURKA inhibitor with high selectivity and favorable pharmacokinetic properties. *ACS Omega* **9(19)**, 21416-21425 (2024).

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, 09/30/2025

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