

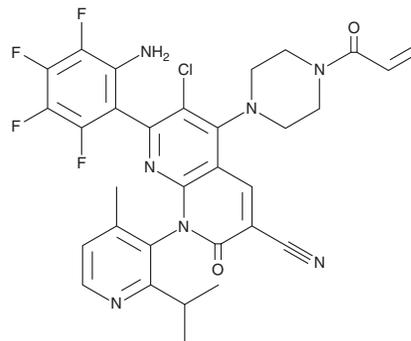
PRODUCT INFORMATION



Gleclirasib

Item No. 45041

CAS Registry No.: 2657613-87-9
Formal Name: 7-(2-amino-3,4,5,6-tetrafluorophenyl)-6-chloro-1,2-dihydro-1-[4-methyl-2-(1-methylethyl)-3-pyridinyl]-2-oxo-4-[4-(1-oxo-2-propen-1-yl)-1-piperazinyl]-1,8-naphthyridine-3-carbonitrile
MF: C₃₁H₂₆ClF₄N₇O₂
FW: 640.0
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

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

Description

Gleclirasib is a covalent inhibitor of K-Ras containing the activating glycine-to-cysteine mutation at position 12 (K-Ras^{G12C}; IC₅₀ = 1.6 nM).¹ It inhibits ERK phosphorylation in K-Ras^{G12C}-expressing NCI H358 non-small cell lung cancer (NSCLC) cells (IC₅₀ = 7.3 nM) and reduces their viability (IC₅₀ = 23 nM). *In vivo*, gleclirasib (10 mg/kg) inhibits tumor growth in an NCI H1373 NSCLC mouse xenograft model.

Reference

1. Li, A., Li, S., Wang, P., *et al.* Design, structure optimization, and preclinical characterization of JAB-21822, a covalent inhibitor of KRAS^{G12C}. *J. Med. Chem.* **68(3)**, 2422-2436 (2025).

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, 02/12/2026

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