

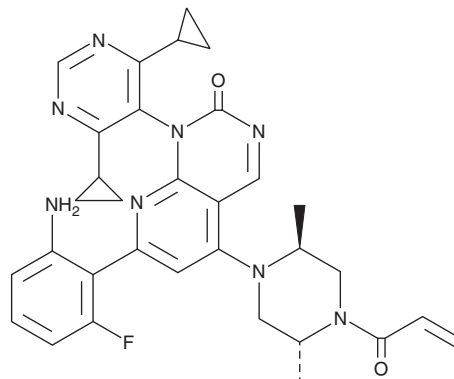
PRODUCT INFORMATION



Garsorasib

Item No. 44250

CAS Registry No.: 2559761-14-5
Formal Name: 7-(2-amino-6-fluorophenyl)-1-(4,6-dicyclopropyl-5-pyrimidinyl)-4-[(2S,5R)-2,5-dimethyl-4-(1-oxo-2-propen-1-yl)-1-piperazinyl]-6-fluoro-pyrido[2,3-d]pyrimidin-2(1H)-one
Synonym: D-1553
MF: C₃₂H₃₂F₂N₈O₂
FW: 598.7
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

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

Description

Garsorasib is an inhibitor of K-Ras containing the activating glycine-to-cysteine mutation at position 12 (K-Ras^{G12C}; IC₅₀ = 18 nM).¹ It selectively binds to GDP-bound K-Ras^{G12C} over GDP-bound wild-type K-Ras in thermal shift assays and selectively inhibits ERK phosphorylation in NCI H358 cells, which express K-Ras^{G12C}, over NCI H1993 cells, which express wild-type K-Ras (IC₅₀s = 6.9 and >1,000 nM, respectively). Garsorasib inhibits proliferation in 11 K-Ras^{G12C}-expressing cancer cell lines (IC₅₀s = 4.6-33 nM) but not in K-Ras^{G12D}- or wild-type-expressing cancer cells (IC₅₀s = >10,000 nM). *In vivo*, garsorasib (30 mg/kg) induces tumor regression in NCI H358 non-small cell lung cancer (NSCLC) and MIA PaCa-2 pancreatic cancer mouse xenograft models.

Reference

1. Shi, Z., Weng, J., Niu, H., *et al.* D-1553: A novel KRAS^{G12C} inhibitor with potent and selective cellular and *in vivo* antitumor activity. *Cancer Sci.* **114**(7), 2951-2960 (2023).

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, 10/03/2025

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