

PRODUCT INFORMATION

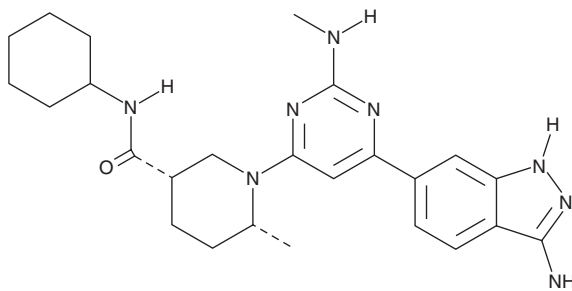


GSK2334470

Item No. 18095

CAS Registry No.: 1227911-45-6
Formal Name: (3S,6R)-1-[6-(3-amino-1H-indazol-6-yl)-2-(methylamino)-4-pyrimidinyl]-N-cyclohexyl-6-methyl-3-piperidinecarboxamide

MF: C₂₅H₃₄N₈O
FW: 462.6
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 258, 328 nm
Supplied as: A crystalline 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

GSK2334470 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2334470 in the solvent of choice, which should be purged with an inert gas. GSK2334470 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GSK2334470 in these solvents is approximately 20 mg/ml.

GSK2334470 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK2334470 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. GSK2334470 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3-Phosphoinositide-dependent protein kinase 1 (PDK1) is a serine-threonine kinase that plays an important role in growth factor signaling by phosphorylating and activating a group of kinases, including protein kinase (PK)B, PKA, and certain isoforms of PKC. GSK2334470 is an inhibitor of PDK1 (IC₅₀ = ~10 nM) that even at 100-fold higher concentrations does not affect the activity of 93 other protein kinases, including that of 13 closely related AGC kinase family members.¹ In HEK293 cells, 30 nM - 1 μM GSK2334470 suppresses T-loop phosphorylation and activation of SGK, S6K1, and RSK, as well as the activation of Akt.¹

Reference

1. Najafov, A., Sommer, E.M., Axten, J.M., *et al.* Characterization of GSK2334470, a novel and highly specific inhibitor of PDK1. *Biochem. J.* **433**(2), 357-369 (2011).

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, 12/05/2022

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