

PRODUCT INFORMATION

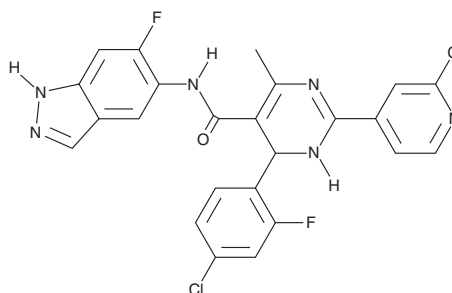


GSK25

Item No. 39163

CAS Registry No.: 874119-56-9
Formal Name: 4-(4-chloro-2-fluorophenyl)-2-(2-chloro-4-pyridinyl)-N-(6-fluoro-1H-indazol-5-yl)-1,4-dihydro-6-methyl-5-pyrimidinecarboxamide

MF: C₂₄H₁₆Cl₂F₂N₆O
FW: 513.3
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

GSK25 is supplied as a solid. A stock solution may be made by dissolving the GSK25 in the solvent of choice, which should be purged with an inert gas. GSK25 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK25 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

GSK25 is an inhibitor of Rho-associated kinase 1 (ROCK1; IC₅₀ = 7 nM for the human kinase).¹ It is selective for ROCK1 over p90 ribosomal S6 kinase 1 (RSK1) and p70 ribosomal S6 kinase (p70S6K; IC₅₀s = 398 and 1,000 nM, respectively) and is greater than 100-fold selective over a panel of 31 additional kinases. GSK25 induces relaxation in precontracted isolated rat aortic rings (EC₅₀ = 256 nM). *In vivo*, GSK25 (30 mg/kg) decreases mean arterial blood pressure in spontaneously hypertensive rats.

Reference

1. Sehon, C.A., Wang, G.Z., Viet, A.Q., *et al.* Potent, selective and orally bioavailable dihydropyrimidine inhibitors of Rho kinase (ROCK1) as potential therapeutic agents for cardiovascular diseases. *J. Med. Chem.* **51(21)**, 6631-6634 (2008).

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, 10/26/2023

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