

PRODUCT INFORMATION



GSK2269557

Item No. 17388

CAS Registry No.: 1254036-77-5

Formal Name: 6-(1H-indol-4-yl)-4-[[[4-(1-methylethyl)-1-piperazinyl]methyl]-2-oxazolyl]-1H-indazole, monohydrochloride

MF: C₂₆H₂₈N₆O • HCl

FW: 477.0

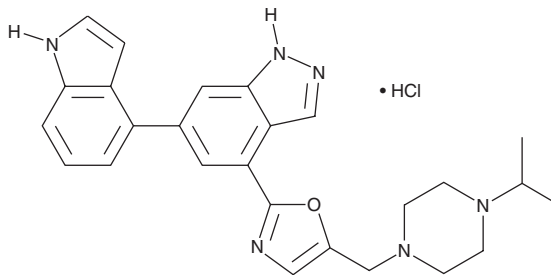
Purity: ≥98%

UV/Vis.: λ_{max}: 223, 303, 318 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

GSK2269557 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2269557 in the solvent of choice, which should be purged with an inert gas. GSK2269557 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GSK2269557 in these solvents is approximately 0.2, 20, and 14 mg/ml, respectively.

GSK2269557 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK2269557 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK2269557 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK2269557 is an inhibitor of PI3Kδ (pK_i = 9.9) that is >1,000-fold selective for PI3Kδ over PI3Kα, β, and γ isoforms.¹ It is reported to be active in a brown Norway rat acute ovalbumin-induced allergic asthma model of Th2-driven lung inflammation (ED₅₀ = 67 μg/kg).¹

Reference

1. Down, K., Amour, A., Baldwin, I.R., *et al.* Optimization of novel indazoles as highly potent and selective inhibitors of phosphoinositide 3-kinase δ for the treatment of respiratory disease. *J. Med. Chem.* **58**(18), 7381-7399 (2015).

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, 12/01/2022

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