

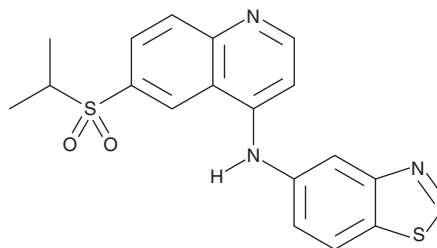
PRODUCT INFORMATION



GSK872

Item No. 23300

CAS Registry No.: 1346546-69-7
Formal Name: N-5-benzothiazolyl-6-[(1-methylethyl)sulfonyl]-4-quinolinamine
MF: C₁₉H₁₇N₃O₂S₂
FW: 383.5
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 252, 362 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

GSK872 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK872 in the solvent of choice, which should be purged with an inert gas. GSK872 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK872 in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

GSK872 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK872 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK872 has a solubility of approximately 0.15 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK872 is an inhibitor of receptor-interacting protein kinase 3 (RIPK3).¹ It is selective for RIPK3, having >1,000-fold selectivity over 300 kinases in a fluorescence polarization assay at a concentration of 1 μM. GSK872 inhibits TNF- and virus-induced necrosis in 3T3-SA fibroblasts. It also inhibits necrosis induced by toll-like receptor 3 (TLR3).

Reference

1. Kaiser, W.J., Sridharan, H., Huang, C., *et al.* Toll-like receptor 3-mediated necrosis via TRIF, RIP3, and MLKL. *J. Biol. Chem.* **288**(43), 31268-31279 (2013).

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, 10/20/2022

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