

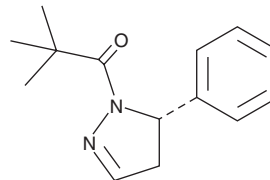
PRODUCT INFORMATION



GSK963

Item No. 39900

CAS Registry No.: 2049868-46-2
Formal Name: 1-[(5S)-4,5-dihydro-5-phenyl-1H-pyrazol-1-yl]-2,2-dimethyl-1-propanone
MF: C₁₄H₁₈N₂O
FW: 230.3
Purity: ≥95%
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

GSK963 is supplied as a solid. A stock solution may be made by dissolving the GSK963 in the solvent of choice, which should be purged with an inert gas. GSK963 is soluble in acetonitrile and DMSO.

Description

GSK963 is an inhibitor of receptor-interacting serine/threonine kinase 1 (RIPK1).¹ It is selective for RIPK1 with less than 50% inhibition of 339 kinases in a panel at 10 μM. Unlike the RIPK1 inhibitor necrostatin-1 (Item No. 11658), GSK963 is inactive against indolamine-2,3-dioxygenase (IDO). GSK963 inhibits necroptosis induced by TNF in the presence of the caspase inhibitor zVAD in L-929 mouse and U937 human cells (IC₅₀s = 1 and 4 nM, respectively). It also inhibits RIPK1-mediated lethal hypothermia induced by TNF with zVAD in mice when administered at a dose of 2 mg/kg in a model of sterile shock.

Reference

- Berger, S.B., Harris, P., Nagilla, R., *et al.* Characterization of GSK'963: A structurally distinct, potent and selective inhibitor of RIP1 kinase. *Cell Death Discov.* **1**, 15009 (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, 12/19/2023

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