

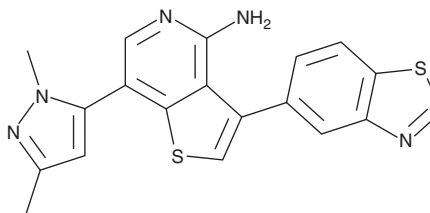
PRODUCT INFORMATION



GSK843

Item No. 39835

CAS Registry No.: 1601496-05-2
Formal Name: 3-(5-benzothiazolyl)-7-(1,3-dimethyl-thieno[3,2-c]pyridin-4-amine
MF: C₁₉H₁₅N₅S₂
FW: 377.5
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

GSK843 is supplied as a solid. A stock solution may be made by dissolving the GSK843 in the solvent of choice, which should be purged with an inert gas. GSK843 is soluble in DMSO. GSK843 is slightly soluble in chloroform and DMSO.

Description

GSK843 is an inhibitor of receptor-interacting serine/threonine kinase 3 (RIPK3; IC₅₀ = 6.5 nM).¹ It is selective for RIPK3 over a panel of 300 additional kinases at 1 μM. GSK843 (10 μM) induces apoptosis in L-929 fibroblasts. It inhibits TNF-α-induced necroptosis in primary human neutrophils in a concentration-dependent manner. GSK843 (3 μM) also inhibits necrosis induced by the toll-like receptor 3 (TLR3) agonist poly (I:C) in IFN-β-primed 3T3-SA fibroblasts.² Intrathecal administration of GSK843 (5 μl of a 13 mM solution) increases the paw withdrawal threshold and latency in a mouse model of inflammatory pain induced by complete Freund's adjuvant.³

References

1. Mandal, P., Berger, S.B., Pillay, S., *et al.* RIP3 induces apoptosis independent of pronecrotic kinase activity. *Mol. Cell* **56**(4), 481-495 (2014).
2. 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).
3. Zhang, Y., Wang, Y., Zhao, W., *et al.* Role of spinal RIP3 in inflammatory pain and electroacupuncture-mediated analgesic effect in mice. *Life Sci.* **306**, 120839 (2022).

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, 02/28/2024

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