

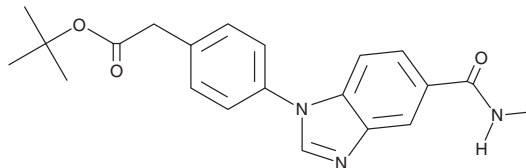
PRODUCT INFORMATION



GSK840

Item No. 35661

CAS Registry No.: 2361146-30-5
Formal Name: 4-[5-[(methylamino)carbonyl]-1H-benzimidazol-1-yl]-benzeneacetic acid, 1,1-dimethylethyl ester
MF: C₂₁H₂₃N₃O₃
FW: 365.4
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

GSK840 is supplied as a solid. A stock solution may be made by dissolving the GSK840 in the solvent of choice, which should be purged with an inert gas. GSK840 is soluble in methanol.

Description

GSK840 is a receptor-interacting serine/threonine kinase 3 (RIPK3) inhibitor (IC₅₀ = 0.3 nM).¹ It is selective for RIPK3 over a panel of 300 additional kinases at 1 μM. GSK840 (0.04-1 μM) inhibits TNF-α-induced necroptosis in primary human neutrophils isolated from human whole blood. It also inhibits LPS-induced neutrophil accumulation in *D. rerio* larvae.²

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. Wen, W., Chen, J., Zhou, Y., *et al.* Loss of Ripk3 attenuated neutrophil accumulation in a lipopolysaccharide-induced zebrafish inflammatory model. *Cell Death Discov.* **8**(1), 88 (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, 11/02/2023

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