

GSK872

Necroptosis inhibitor / RIP3 inhibitor

GSK872 is a potent and selective inhibitor of receptor interacting protein 3 (RIP3). It is able to prevent necroptosis, virus-induced necrosis, and TLR3-induced necrosis. GSK872 can induce Caspase8-mediated apoptosis at higher concentrations.

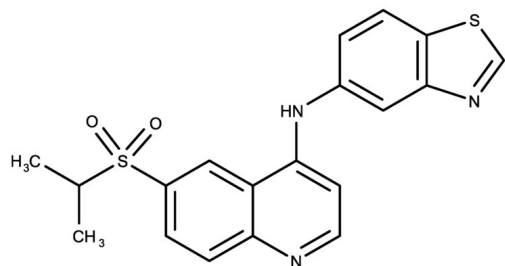
Ordering Information

[Order Online »](#)

ENZ-CHM499-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-(6-Propan-2-ylsulfonylquinolin-4-yl)-1,3-benzothiazol-5-amine

Appearance Off-white or pale yellow solid.

CAS 1346546-69-7

Couple Target RIP

Couple Type Inhibitor

Formula $C_{19}H_{17}N_3O_2S_2$

Identity Determined by NMR.

MW 383.49

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (>25 mg/ml).

Last modified: February 27, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com