

Necrostatin-5

Necroptosis inhibitor

Necroptosis inhibitor. In contrary to the mechanism of necrostatin-1 (Prod.No. BML-AP309) it was shown that necrostatin-5 inhibits RIP1 kinase indirectly. A potent inhibitor of immunoprecipitated, but not recombinant RIP1.

The necrostatins inhibit non-apoptotic cell death pathway by inhibition of the death domain receptor-associated adaptor kinase RIP1. There are three different mechanisms: T-loop dependent inhibition by necrostatin-1, partially T-loop independent inhibition by necrostatin-3 and indirect inhibition of RIP1 by necrostatin-5.

Citations: 3

[View Online »](#)

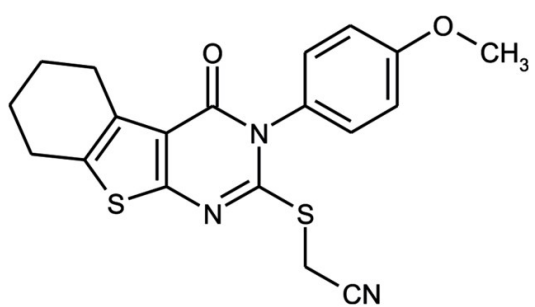
Ordering Information

[Order Online »](#)

ALX-430-169-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Nec-5, 3-p-Methoxyphenyl-5,6-tetramethylenothieno[2,3-d]pyrimidin-4-one-2-mercaptoethylcyanide

Appearance White to off-white solid.

CAS 337349-54-9

Couple Target RIP

Couple Type Inhibitor

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

MW 383.5

Purity ≥98% (HPLC, TLC)

Solubility Soluble in DMSO (>10mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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