

Necrostatin-1

Inhibits necroptosis

Necrostatin-1 inhibits necroptosis, a non-apoptotic cell death pathway. Inhibits the loss of mitochondrial membrane potential in TNF α -treated Jurkat cells (EC_{50} =490 nM). Does not inhibit FAS-induced apoptosis and has no effect on apoptotic morphology. It displays a pronounced protective effect in a mouse model of ischemic brain injury and inhibits myocardial cell death. Inhibits RIP1 kinase the key upstream kinase involved in the activation of necroptosis (EC_{50} =180nM).

Citations: 65

[View Online »](#)

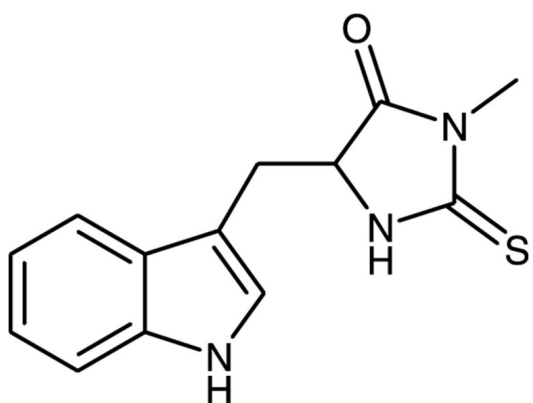
Ordering Information

[Order Online »](#)

BML-AP309-0020	20mg
BML-AP309-0100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Methylthiohydantoin-DL-tryptophan
Appearance	Yellow solid.
CAS	4311-88-0
Couple Target	RIP
Couple Type	Inhibitor
Formula	$C_{13}H_{13}N_3OS$
MW	259.3
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (20mg/ml) or 100% ethanol (5mg/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-
be@enzolifesciences.com](mailto:info-be@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)