

Panepoxydone

NF- κ B pathway inhibitor

Antibiotic. Inhibitor of NF- κ B activation by preventing the phosphorylation of I κ B protein and thereby interrupts the signalling pathway. Strongly inhibits the expression of several NF- κ B dependent pro-inflammatory genes.

Citations: 13

[View Online »](#)

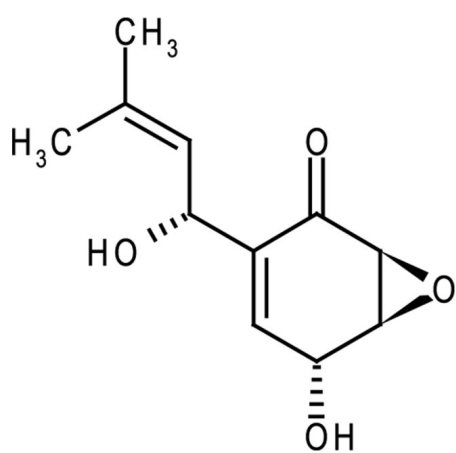
Ordering Information

[Order Online »](#)

ALX-350-109-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep cool and dry. After reconstitution, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow oil.
CAS	31298-54-1
Couple Target	IkappaB, NF-kappaB
Couple Type	Inhibitor
Formula	$C_{11}H_{14}O_4$
MW	210.2
Purity	≥95% (HPLC, MS)
Solubility	Soluble in DMSO, 100% ethanol or methanol.
Source	Isolated from <i>Lentinus crinitus</i> .

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)