

# Panepoxydone

NF-κB pathway inhibitor

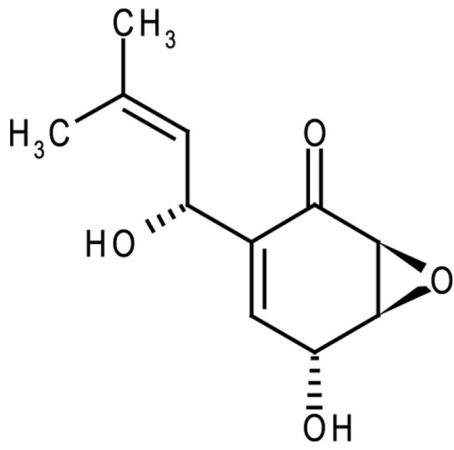
Antibiotic. Inhibitor of NF-κB activation by preventing the phosphorylation of IκB protein and thereby interrupts the signalling pathway. Strongly inhibits the expression of several NF-κB dependent pro-inflammatory genes.

**Citations:** 13

[View Online »](#)

**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-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-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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