

Pseudolaric acid B

Tubulin ligand

Activates PPAR α and the phospholipase C (PLC) signalling pathway. Stimulates peroxisomal fatty acyl-CoA oxidase activity. These effects can be blocked by staurosporine (Prod. No. ALX-380-014). Exhibits significant cytotoxic activities against numerous tumor cell lines. Binds tubulin. Inhibits angiogenesis. Reduces HIF-1 α protein levels. Induces apoptosis. Antifungal.

Citations: 16

[View Online »](#)

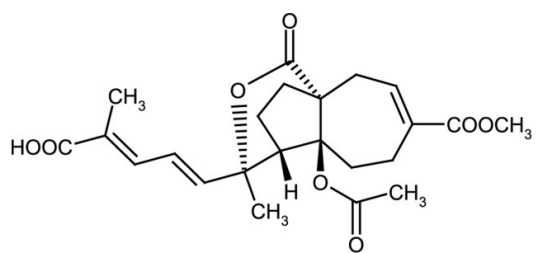
Ordering Information

[Order Online »](#)

ALX-350-108-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Pseudolarix acid B, PLAB, PAB
Appearance	Off-white to light brown solid.
CAS	82508-31-4
Couple Target	Tubulin
Couple Type	Ligand
Formula	$C_{23}H_{28}O_8$
MW	432.5
Purity	≥95% (HPLC)
RTECS	KJ4620000
Solubility	Soluble in DMSO, 100% ethanol or chloroform.
Source	Isolated from <i>Pseudolarix kaempferi</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-
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)