

Sauchinone

Cytoprotective and antioxidant

Diastereomeric lignan with cytoprotective and antioxidant activities in cultured hepatocytes. Inhibitor of LPS-inducible iNOS (NOS II), TNF- α and COX-2 expression in macrophages through suppression of I κ B α phosphorylation and p65 nuclear translocation and of C/EBP and/or AP-1 activation, which may have constitutive anti-inflammatory effects. Suppressor of NF- κ B by inhibiting transactivation activity of the RelA subunit. Inhibits staurosporine (Prod. No. ALX-380-014)-induced apoptosis. Inhibits RANKL-induced osteoclastogenesis by reducing ROS generation.

Citations: 6

[View Online »](#)

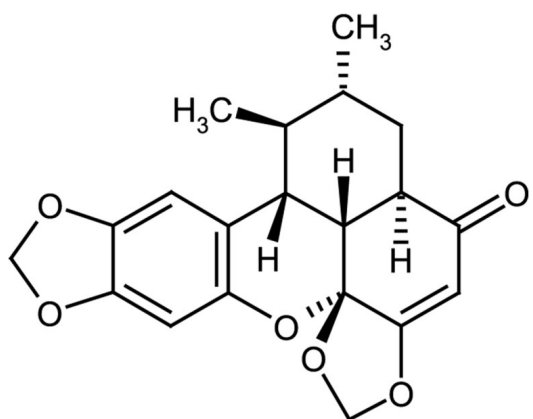
Ordering Information

[Order Online »](#)

ALX-350-116-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white powder.

CAS 177931-17-8

Couple Target COX, NOS, TNF

Couple Type Inhibitor

Formula $C_{20}H_{20}O_6$

MW 356.1

Purity ≥95% (HPLC)

Solubility Soluble in DMSO or chloroform.

Source Isolated from *Saururus chinensis*.



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)