

Deguelin

COX-2 inhibitor

Inhibitor of Akt (protein kinase B; PKB) in an *in vitro* lung carcinogenesis progression model. Inhibits cell proliferation - cells accumulate in the G2-M phase of the cell cycle. Premalignant and malignant human bronchial epithelial cells are highly sensitive to deguelin compared to normal cells. Induces apoptosis. Exhibits a marked inhibitory effect on mouse skin tumor promotion in an *in vivo* two-stage carcinogenesis test. Inhibits cyclooxygenase-2 (COX-2) expression.

Citations: 9

[View Online »](#)

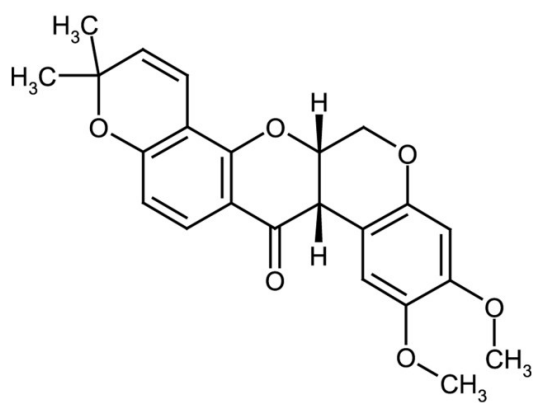
Ordering Information

[Order Online »](#)

ALX-350-118-M005	5mg
ALX-350-118-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name [-]-Deguelin

Appearance White to yellow powder.

CAS 522-17-8

Couple Target COX

Couple Type Inhibitor

Formula $C_{23}H_{22}O_6$

MI 14: 2863

MW 394.4

Purity ≥95% (HPLC)

RTECS DX1500000

Solubility Soluble in acetone, acetonitrile, DMSO (25mg/ml) or 100% ethanol (25mg/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-
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)