

Lupeol

Anti-inflammatory, antioxidant and antimutagenic compound

Anti-inflammatory, antioxidant and antimutagenic compound. Shows preventive effects on DMBA induced DNA alkylation damage in mouse skin. Modulator of NF- κ B and PI(3)K/Akt (protein kinase B; PKB) pathways. Induces Fas-mediated apoptosis via inhibition of Ras signalling. Inhibits tumor growth in mouse *in vivo*.

Citations: 11

[View Online »](#)

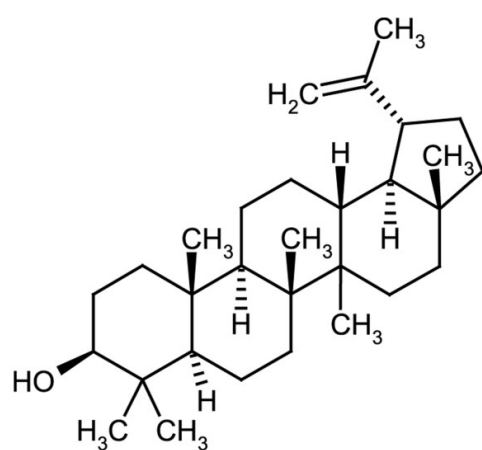
Ordering Information

[Order Online »](#)

ALX-350-355-M010	10mg
ALX-350-355-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Fagarasterol, 20(29)-Lupen-3β-ol, 3β-Hydroxy-20(29)-lupene

Appearance White to off-white solid.

CAS 545-47-1

Formula $C_{30}H_{50}O$

MI 14: 5608

MW 426.7

Purity ≥98% (HPLC)

RTECS OK5763000

Solubility Soluble in ether or warm 100% ethanol; insoluble in water

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)