

Betulinic acid (high purity)

Antitumor and anti-HIV agent

Antitumor and anti-HIV agent. Induces apoptosis by activating mitochondrial permeability transition (MPT). Inhibits NF- κ B activation and NF- κ B-regulated gene expression induced by carcinogens and inflammatory stimuli. Decreases expression of Bcl-2 and cyclin D1. Potent proteasome activator. Triggers CD95- and p53-independent apoptosis via activation of caspases. It induces mitochondrial permeability pore opening and the release of cytochrome c and apoptosis-inducing factor. Displays anti-HIV activity. EC₅₀ for inhibition of HIV-1 replicaton in H9 lymphocytes: 1.4 μ M.

Citations: 22

[View Online »](#)

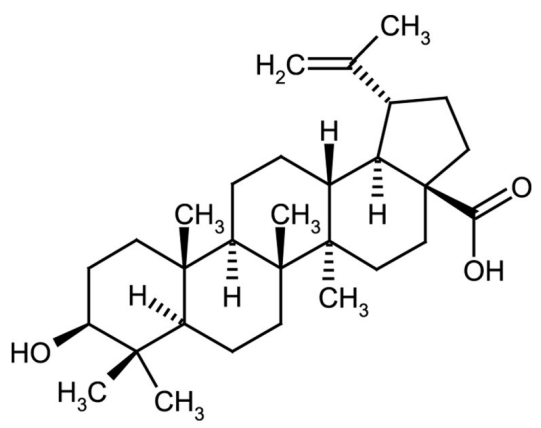
Ordering Information

[Order Online »](#)

ALX-350-277-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep cool and dry.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3β-Hydroxy-20(29)-lupaene-28-oic acid
Appearance	White to off-white powder.
CAS	472-15-1
Formula	$C_{30}H_{48}O_3$
MW	456.7
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO.
Source	Isolated from <i>Platanus acerifolia</i> (plane) tree bark.

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-
be@enzolifesciences.com](mailto:info-be@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)