

Betulinic acid (~95%)

Proteasome activator

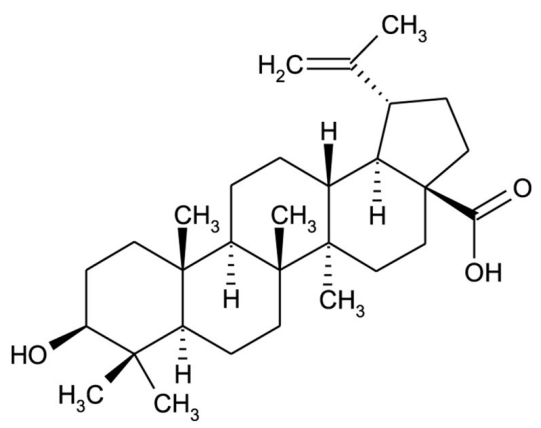
Antitumor and anti-HIV agent. Induces apoptosis by activating mitochondrial permeability transition (MPT). Inhibits NF- κ B activation and NF- κ B-regulated gene expression induced by carcinogens and inflammatory stimuli. Decreases expression of Bcl-2 and cyclin D1. Potent proteasome activator.

Citations: 19

[View Online »](#)

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 3β-Hydroxy-20(29)-lupaene-28-oic acid

Appearance White to off-white powder.

CAS 472-15-1

Couple Target Proteasome

Couple Type Activator

Formula $C_{30}H_{48}O_3$

MW 456.7

Purity ≥90% (HPLC)

Solubility Soluble in DMSO.

Source Isolated from *Platanus acerifolia* (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)