

Pristimerin

Antitumor agent

Triterpenoid. Antitumor agent. Induces caspase-dependent apoptosis by inhibiting the chymotrypsin-like activity of the proteasome. Inhibits NF- κ B activation via inhibition of IKK α or IKK β . Inhibits induction of inducible nitric oxide synthase (iNOS; NOS II). Shows *in vitro* antimalarial and insecticidal activity. Induces mitochondrial cell death. Has antiproliferative, antifungal and antiviral properties. Potent and reversible monoacylglycerol lipase (MGL) inhibitor.

Citations: 17

[View Online »](#)

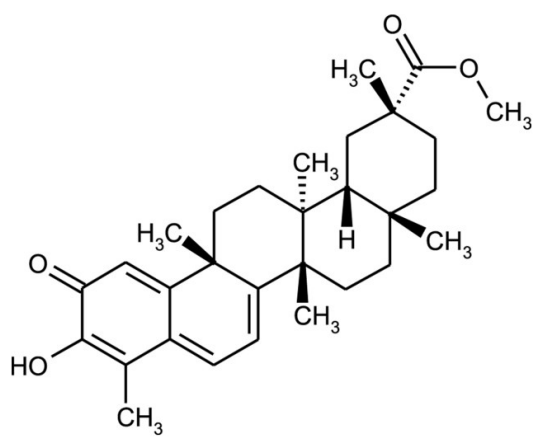
Ordering Information

[Order Online »](#)

ALX-350-411-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Celastrol methyl ester
Appearance	Orange crystalline solid.
CAS	1258-84-0
Couple Target	Monoacylglycerol lipase, NOS
Couple Type	Inhibitor
Formula	$C_{30}H_{40}O_4$
Identity	Determined by IR.
MI	14: 10370
MW	464.6
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO or ethyl acetate; insoluble in water.
Source	Isolated from <i>Tripterygium wilfordii</i> Hook.F.

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)