

# Celastrol

## Anti-inflammatory and immunosuppressive

Cell permeable triterpenoid; anti-inflammatory and immunosuppressive compound. Antioxidant. Suppresses LPS-induced cytokine release in macrophages and monocytes ( $IC_{50}$ =40nM for IL-1 $\beta$  and IL-1 $\alpha$ , 80nM for IL-6, 110nM for prostaglandin E<sub>2</sub> and 210nM for IL-8 and TNF- $\alpha$ ). Suppresses nitric oxide (NO) production ( $IC_{50}$ =230nM) and LPS-induced NF- $\kappa$ B activation ( $IC_{50}$ =270nM). Inhibits chymotrypsin-like activity of 20S proteasome ( $IC_{50}$ =2.5 $\mu$ M). Inhibits lipid peroxidation induced by ADP and Fe<sup>2+</sup> in rat liver mitochondria ( $IC_{50}$ =7 $\mu$ M). Induces expression of a wider set of heat shock proteins. Inhibits topoisomerase II.

Citations: 20

[View Online »](#)

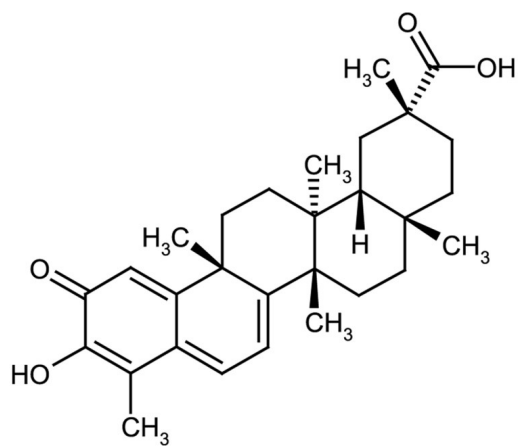
## Ordering Information

[Order Online »](#)

ALX-350-332-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	Tripterine, 3-Hydroxy-24-nor-2-oxo-1(10),3,5,7-friedelatetraen-29-oic acid
Appearance	Red cubic crystals.
CAS	34157-83-0
Couple Target	Proteasome, Topoisomerase
Couple Type	Inhibitor
Formula	$C_{29}H_{38}O_4$
MI	14: 10370
MW	450.6
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol, DMSO or dimethyl formamide; insoluble in water.
Source	Isolated from the root of <i>Tripterygium</i> sp.

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)