

Triptolide

Immunosuppressant

Immunosuppressive agent with anti-inflammatory and antitumor properties. More effective in preventing T cell proliferation and interferon- γ production than FK506 (Prod. No. ALX-380-008). Induces apoptosis in T cells by activating DEVD cleaving caspases. Blocks TNF- α mediated induction of c-IAP1 and c-IAP2 and NF- κ B activation. Inhibits induction of MMP-3, MMP-13, and ADAMTS-4 in cell culture.

Citations: 11

[View Online »](#)

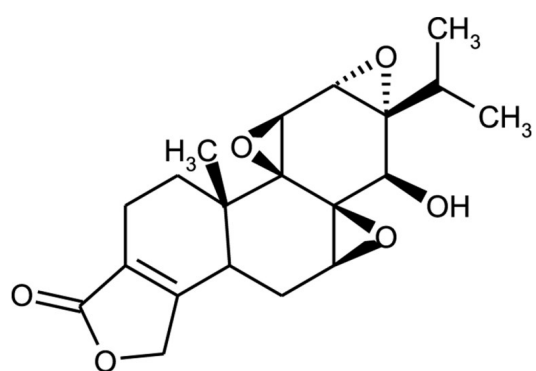
Ordering Information

[Order Online »](#)

ALX-350-259-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PG490
Appearance	White to off-white solid.
CAS	38748-32-2
Couple Target	Caspase
Couple Type	Activator
Formula	$C_{20}H_{24}O_6$
MW	360.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (1mg/ml) (1h at 37°C).
Source	Isolated from <i>Tripterygium wilfordii</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com