

Tryptanthrin

Inhibitor of prostaglandin and leukotriene synthesis

Potent inhibitor of prostaglandin and leukotriene synthesis in various cell lines. Selective inhibitor of cyclooxygenase 2 (COX-2). Inhibitor of inducible nitric oxide synthase (iNOS; NOS II) expression. Inhibits the production of IFN- γ and IL-2 after stimulation of Peyer's patch lymphocytes with *staphylococcal* enterotoxin B (SEB). Inhibits P-glycoprotein (Pgp) (expressed by the MDR1 gene) and reverses doxorubicin (Prod. No. BML-GR319) resistance on breast cancer cells.

Tryptanthrin occurs as a natural product in various plants: *Isatis tinctoria*, *I. indigotica*, *Polygonum tinctorium*, *Couroupita guianensis*, *Wrightia tinctoria*, *Calanthe discolor*, *Phaius mishmensis*, *Strobilanthes cusia*.

Citations: 10

[View Online »](#)

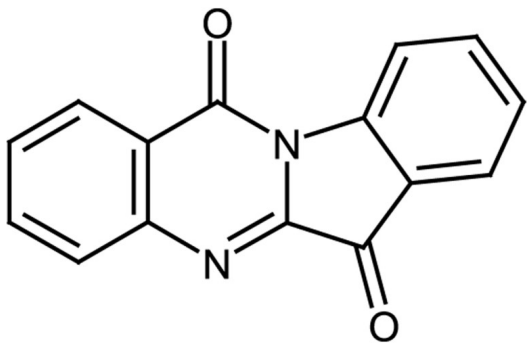
Ordering Information

[Order Online »](#)

ALX-270-360-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at +4°C in the dark. Stock solutions made in DMSO must be stored in the dark no longer than one month.
Handling	Protect from light and moisture.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Yellow needles.
CAS	13220-57-0
Couple Target	COX, Lipoxygenase
Couple Type	Inhibitor
Formula	$C_{15}H_8N_2O_2$
Identity	Determined by 1H-NMR, 13C-NMR, IR, MS and UV.
MW	248.2
Purity	≥98% (1H-NMR, HPLC)
Solubility	Soluble in DMSO (8mg/ml) or methanol (0.3mg/ml); sparingly soluble in water (2.5µg/ml).
Source	Synthetic.

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

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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