

Evodiamine

TRPV1 antagonist

A cell permeable, non-pungent TRPV1 antagonist ($K_i=5.95\mu\text{M}$ in rat) that induces apoptosis in U937 cells. Inhibits nitric oxide (NO) production. Also inhibits tumor cell migration *in vitro*. Is effective against multidrug-resistant cancer cells. Inhibits angiogenesis.

Citations: 9

[View Online »](#)

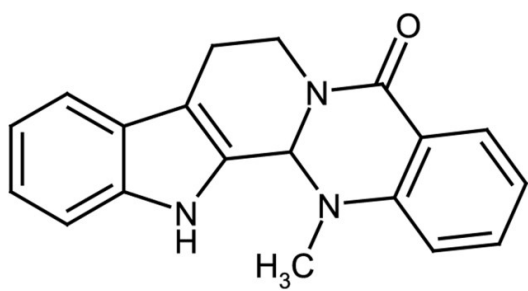
Ordering Information

[Order Online »](#)

ALX-350-330-M050	50mg
------------------	------

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 8,13,13b,14-Tetrahydro-14-methylindolo[2'3'-3,4]pyrido[2,1-b]quinazolin-5-[7H]-one

Appearance Off-white to light yellow solid.

CAS 518-17-2

Couple Target TRP channel

Couple Type Ligand

Formula $C_{19}H_{17}N_3O$

MI 14: 3908

MW 303.4

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (5mg/ml; warm to 40°C).

Source Synthetic. Originally isolated from *Evodia rutaecarpa*.

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)