

# 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 »](#)

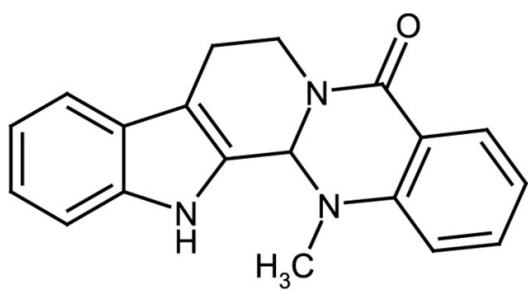
## Ordering Information

[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-  
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)