

Capsazepine

Capsaicin antagonist

Analog of capsaicin (Prod. No. BML-EI125-0200) that acts as a capsaicin antagonist. Inhibits resiniferatoxin and capsaicin-induced contractions of guinea pig trachea. Inhibits C-fiber-smooth muscle contractions induced by capsaicin, bradykinin and neurokinin A. Increases intracellular Ca^{2+} levels.

Citations: 14

[View Online »](#)

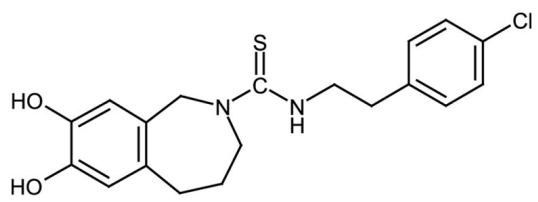
Ordering Information

[Order Online »](#)

ALX-550-145-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-[2-(4-Chlorophenyl)ethyl]-1,3,4,5-tetrahydro-7,8-dihydroxy-2H-2-benzazepine-2-carbothioamide

Appearance White to off-white powder.

CAS 138977-28-3

Formula $C_{19}H_{21}ClN_2O_2S$

MW 376.9

Purity ≥97% (HPLC, TLC)

Solubility Soluble in methanol, 100% ethanol (25mg/ml) or DMSO (25mg/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)