

Dihydrocapsaicin

Dihydro-analog and congener of capsaicin (Prod. No. BML-EI125) in chili peppers (*Capsicum*). Like capsaicin it is irritant. Dihydrocapsaicin accounts for about 22% of the total capsaicinoids mixture and has about the same pungency as capsaicin. Antioxidant. Reduces oxidation of serum lipids. Mutagenic.

Citations: 5

[View Online »](#)

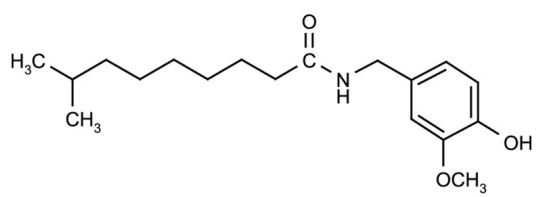
Ordering Information

[Order Online »](#)

ALX-350-052-M050	50mg
------------------	------

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 8-Methyl-N-vanillylnonanamide

Appearance White to off-white solid.

CAS 19408-84-5

Formula $C_{18}H_{29}NO_3$

MW 307.4

Purity ≥85% (HPLC)

RTECS RA5998000

Solubility Soluble in chloroform, DMSO or 100% ethanol; sparingly soluble in water.

Source Isolated from *Capsicum* fruit.

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