

(-)-Eburnamone

Antihypoxic. Vasodilator.

Cerebral metabolic stimulant; shows antihypoxic effects. Vasodilator.

Citations: 4

[View Online »](#)

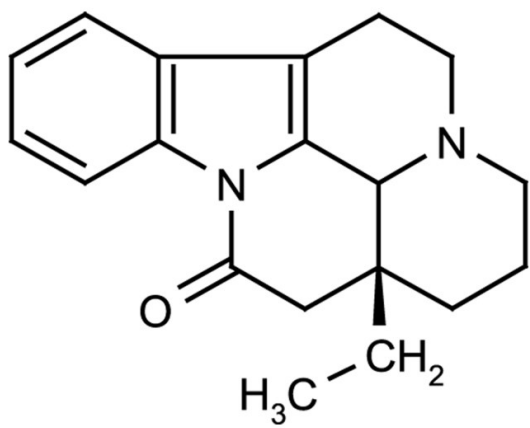
Ordering Information

[Order Online »](#)

ALX-350-216-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (3 α ,16 α)-Eburnamenin-14(15H)-one, Vincamone, Vinburnine

Appearance White to slightly yellow crystalline powder.

CAS 4880-88-0

Formula C₁₉H₂₂N₂O

MI 14: 3487

MW 294.4

Purity ≥99% (Assay)

RTECS YY8575570

Solubility Soluble in chloroform.

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)