

(+)-Egenine

First phthalideisoquinoline hemiacetal from a natural source. Inhibits Ca^{2+} currents and contractile responses *in vitro*.

Citations: 6

[View Online »](#)

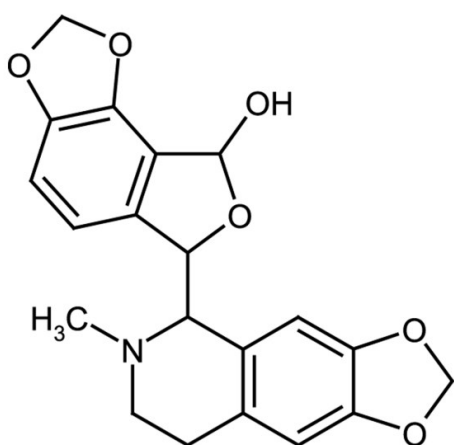
Ordering Information

[Order Online »](#)

ALX-350-362-M025	25mg
ALX-350-362-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 6-(6-Methyl-5,6,7,8-tetrahydro[1,3]dioxolo[4,5-g]isoquinolin-5-yl)-6,8-dihydrofuro[3,4-E][1,3]benzodioxol-8-ol, (+)-Decumbensine

Appearance White to off-white crystalline solid.

CAS 6883-44-9

Formula $C_{20}H_{19}NO_6$

MW 369.4

Purity ≥98% (HPLC)

Solubility Soluble in ethyl acetate or chloroform; sparingly soluble in 100% ethanol or ether; practically insoluble in water.

Source Isolated from *Fumaria vaillantii* or *Corydalis decumbens*.

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)