

(±)-Verapamil . hydrochloride

Ca²⁺ channel inhibitor

Blocks slow voltage-sensitive L-type Ca²⁺ channels. Adrenergic antagonist. Inhibits platelet-activating factor and PMA-stimulated prostaglandin production in Kupffer cells. Induces apoptosis. *In vitro* resistance modifier in drug resistant human tumor cell lines.

Citations: 22

[View Online »](#)

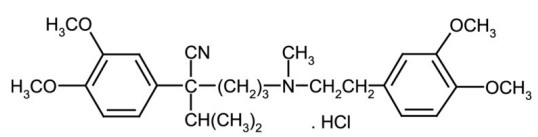
Ordering Information

[Order Online »](#)

ALX-550-306-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-[(3,4-Dimethoxyphenylethyl)methylamino]-2-(3,4-dimethoxyphenyl)-2-isopropyl-valeronitrile . HCl

Appearance White crystalline powder.

CAS 152-11-4

Couple Target Calcium channel

Couple Type Inhibitor

Formula $C_{27}H_{38}N_2O_4 \cdot HCl$

Identity Determined by IR.

MI 14: 9950

MW 454.6 . 36.5

Purity ≥99% (HPLC)

RTECS YV8320000

Solubility Soluble in 100% ethanol (26mg/ml), methanol or water (83mg/ml).

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