

Flunarizine . dihydrochloride

Ca²⁺ channel inhibitor

Selectively blocks the low threshold (T-type) calcium channel in rat hippocampal CA1 pyramidal neurons. Inhibits Ca²⁺-influx-induced vasoconstriction of peripheral and cerebral blood vessels with little effect on myogenic activity and Ca²⁺-dependent myocardial contractility.

Citations: 3

[View Online »](#)

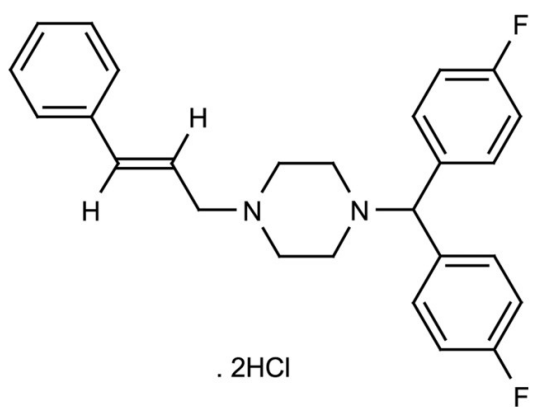
Ordering Information

[Order Online »](#)

ALX-550-268-G001	1g
------------------	----

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 1-[bis(4-Fluorophenyl)methyl]-4-(3-phenyl-2-propenyl)piperazine . 2HCl

Appearance White solid.

CAS 30484-77-6

Couple Target Calcium channel

Couple Type Inhibitor

Formula $C_{26}H_{26}F_2N_2 \cdot 2HCl$

Identity Determined by IR.

MI **14: 4144**

MW 404.5 . 72.9

Purity ≥99% (HPLC)

Solubility Soluble in DMSO (25mg/ml), 100% ethanol (5mg/ml) or chloroform:methanol (1:1; 50mg/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-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)