

Bepriidil . hydrochloride

Ca²⁺ channel inhibitor

Non-selective Ca²⁺ channel blocker. Stimulates the binding of 1,4-dihydropyridine-based drugs to Ca²⁺ channels. Inhibits Na⁺/Ca²⁺ exchange.

Citations: 5

[View Online »](#)

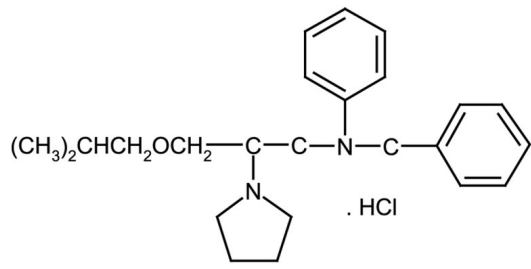
Ordering Information

[Order Online »](#)

ALX-550-237-M010	10mg
ALX-550-237-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Store tightly sealed.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	1-Isobutoxy-2-pyrrolidino-3-(N-benzylanilino)propane . HCl
Appearance	White to off-white solid.
CAS	74764-40-2
Couple Target	Calcium channel
Couple Type	Inhibitor
Formula	$C_{24}H_{34}N_2O \cdot HCl$
MI	14: 1150
MW	366.6 . 36.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol; insoluble in water.

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