

Pinacidil . monohydrate

K⁺ channel activator

K⁺ channel activator. Antihypertensive.

Citations: 1

[View Online »](#)

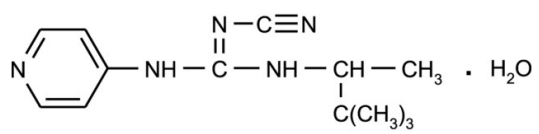
Ordering Information

[Order Online »](#)

ALX-550-285-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Handling Keep under inert gas.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (±)-N-Cyano-N'-4-pyridinyl-N''-(1,2,2-trimethylpropyl)guanidine monohydrate

Appearance White to off-white powder.

CAS 85371-64-8

Couple Target Potassium channel

Couple Type Activator

Formula $C_{13}H_{19}N_5 \cdot H_2O$

MI 14: 7438

MW 245.3 . 18.0

Purity ≥98% (HPLC)

Solubility Soluble in 100% ethanol (14mg/ml), DMSO (28mg/ml) or dilute aqueous acid; 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-
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)