

Amiloride . hydrochloride . dihydrate

Ca²⁺ channel inhibitor

Selectively blocks low threshold (T-type) calcium channel ($K_D=30\mu\text{M}$) with no effect on L-type channels or tetrodotoxin Na⁺ channels. Inhibits the epithelial tetrodotoxin insensitive Na⁺ channel ($IC_{50}=1\mu\text{M}$). Clinically useful K⁺ sparing diuretic.

Citations: 7

[View Online »](#)

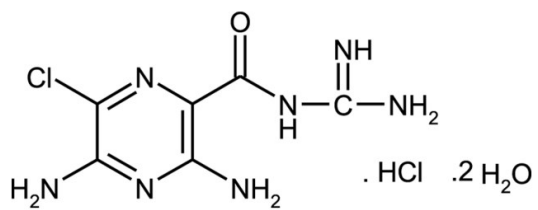
Ordering Information

[Order Online »](#)

ALX-550-212-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 Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3,5-Diamino-N-(aminoiminomethyl)-6-chloropyrazinecarboxamide . HCl . 2H₂O

Appearance Light yellow solid.

CAS 17440-83-4

Couple Target Calcium channel

Couple Type Inhibitor

Formula C₆H₈ClN₇O . HCl . 2H₂O

MW 229.6 . 36.5 . 36.0

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25mg/ml) or water (1mg/ml warm).

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)