

5-(N-Ethyl-N-isopropyl)amiloride

Na⁺/H⁺ antiporter inhibitor

Selective inhibitor of the Na /H antiporter.

Citations: 11

[View Online »](#)

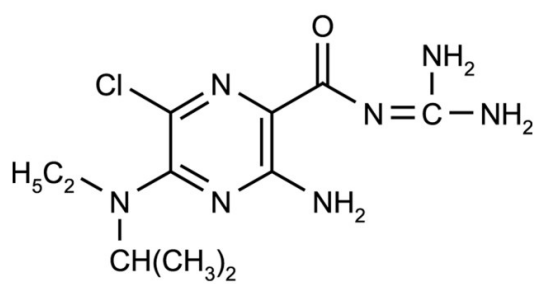
Ordering Information

[Order Online »](#)

ALX-550-266-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-(N-Ethyl-N-isopropyl)amiloride, EIPA
Appearance	Pale yellow to green solid.
CAS	1154-25-2
Couple Target	Na ⁺ /H ⁺ antiporter
Couple Type	Inhibitor
Formula	C ₁₁ H ₁₈ ClN ₇ O
Identity	Determined by NMR.
MW	299.8
Purity	≥95% (HPLC)
Solubility	Slightly soluble in Acetonitrile/DMSO.

Last modified: June 16, 2025



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)