

Lidocaine . hydrochloride . hydrate

Na⁺ channel blocker

Na⁺ channel blocker.

Citations: 2

[View Online »](#)

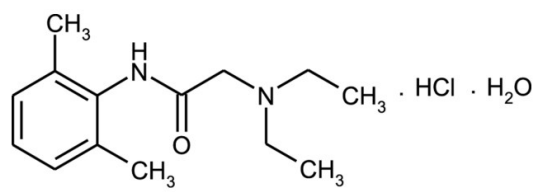
Ordering Information

[Order Online »](#)

ALX-550-251-G005	5g
------------------	----

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 2-Diethylamino-N-(2',6'-dimethylphenyl)acetamide . HCl . xH₂O

Appearance White to off-white crystalline solid.

CAS 6108-05-0

Couple Target Sodium channel

Couple Type Blocker

Formula C₁₄H₂₂N₂O . HCl . xH₂O

MI 14: 5482

MW 234.3 . 36.5 (anhydrous basis)

Purity ≥98% (Assay)

RTECS AN7700000

Solubility Soluble 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)