

Procainamide . hydrochloride

Na⁺ channel blocker

Na⁺ channel blocker and class I antiarrhythmic.

Citations: 4

[View Online »](#)

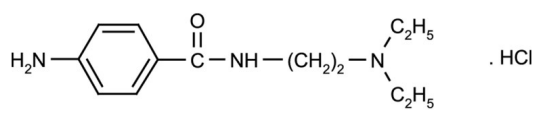
Ordering Information

[Order Online »](#)

ALX-550-287-G010	10g
------------------	-----

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 4-Amino-N-[2'-(diethylamino)ethyl]benzamide . HCl

Appearance White to off-white solid.

CAS 614-39-1

Couple Target Sodium channel

Couple Type Blocker

Formula $C_{13}H_{21}N_3O \cdot HCl$

MI 14: 7756

MW 235.3 . 36.5

Purity $\geq 98\%$ (HPLC)

Solubility Soluble in water; slightly soluble in 100% ethanol or methanol; insoluble in ethyl acetate.

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)