

Lorcainide . HCl

Na⁺ channel blocker

Possesses both local anesthetic and antiarrhythmic properties (class 1C agent). Elevates the ventricular fibrillation threshold and protects against ouabain-induced arrhythmias in animal models. Its mechanism of action is primarily inhibition of sodium channels in a voltage- and frequency-dependent fashion. Also inhibits the microsomal Na⁺,K⁺-ATPase in a concentration-dependent manner (IC₅₀=33.5μM) . Reduces neuronal death in cellular models of hypoxic injury when combined with a glutamate receptor antagonist and blocks aconitine-induced mortality in a rat model (ED₅₀=30mg/kg).

Citations: 4

[View Online »](#)

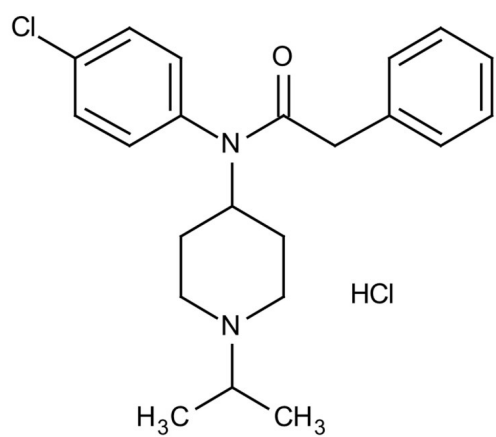
Ordering Information

[Order Online »](#)

BML-NA143-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-(4-Chlorophenyl)-N-[1-(1-isopropyl)-4-piperidiny]phenylacetamide . HCl
Appearance	White solid.
CAS	58934-46-6
Couple Target	Potassium channel, Sodium channel
Couple Type	Blocker
Formula	$C_{22}H_{27}ClN_2O \cdot HCl$
MW	370.9 . 36.5
Purity	≥98% (TLC)
Solubility	Soluble in water (>30mg/ml), DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).

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-
be@enzolifesciences.com](mailto:info-be@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)