

Chlorpromazine . hydrochloride

Antipsychotic working on a variety of
receptors in the CNS

Inhibits calmodulin stimulation of cyclic nucleotide phosphodiesterase. Dopamine antagonist. D₂ antagonist. H₁ antagonist. Inhibits TNF-α production. Potent PLA₂ inhibitor. Inhibits nitric oxide synthase (NOS) in mouse brain and prevents lipopolysaccharide induction of NOS in mouse lung. Antipsychotic.

Citations: 13

[View Online »](#)

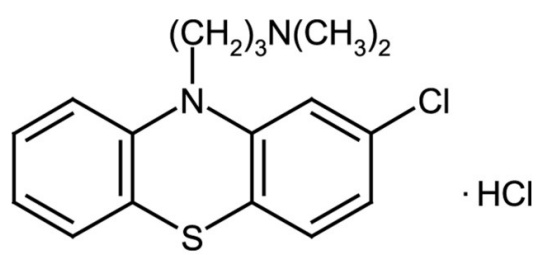
Ordering Information

[Order Online »](#)

ALX-270-171-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Not stable in aqueous solution. We strongly recommend always to use freshly prepared solutions.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature – Dangerous Good

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Chloro-10-(3-dimethylaminopropyl)phenothiazine . HCl
Appearance	White to off-white powder.
CAS	69-09-0
Couple Target	Acetylcholine receptor, Adrenergic receptor, Calmodulin, Dopamine receptor, Histamine receptor, PDE, PLA, Serotonin receptor
Couple Type	Inhibitor
Formula	$C_{17}H_{19}ClN_2S \cdot HCl$
MI	14: 2185
MW	355.3
Purity	≥98% (UHPLC)
RTECS	SO1750000
Solubility	Soluble in 100% ethanol, methanol or water (50mg/ml).



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com