

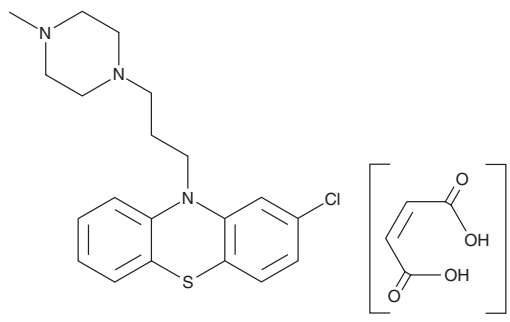
PRODUCT INFORMATION



Prochlorperazine (maleate)

Item No. 20742

CAS Registry No.: 84-02-6
Formal Name: 2-chloro-10-[3-(4-methyl-1-piperazinyl)propyl]-10H-phenothiazine, dimaleate
MF: C₂₀H₂₄ClN₃S • 2C₄H₄O₄
FW: 606.1
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 312 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Prochlorperazine (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the prochlorperazine (maleate) in the solvent of choice, which should be purged with an inert gas. Prochlorperazine (maleate) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Prochlorperazine (maleate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, prochlorperazine (maleate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Prochlorperazine (maleate) has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Prochlorperazine is a dopamine D₂ receptor antagonist with K_i values of 4.7 and 2.9 nM for rat recombinant D₂ receptors in CHO cells and rat striatal membranes, respectively.^{1,2} It also binds to rat recombinant D₃ receptors expressed in CHO cells (K_i = 35 nM) and to the serotonin (5-HT) receptor subtype 5-HT₃ in N1E-115 mouse neuroblastoma cell membranes (K_i = 1,200 nM).^{1,3} Prochlorperazine (2 mg/kg) increases the latency to paw licking in a hot plate test, indicating analgesia, an effect that is blocked by antisense oligonucleotides against the M₁ muscarinic receptor.⁴ It also inhibits emesis induced by apomorphine (Item No. 16094) in dogs (ED₅₀ = 0.34 mg/kg).⁵ Formulations containing prochlorperazine have been used in the treatment of psychotic disorders and as antiemetics.

References

1. Sokoloff, P., Giros, B., Martres, M.P., *et al.* Molecular cloning and characterization of a novel dopamine receptor (D₃) as a target for neuroleptics. *Nature* **137(6289)**, 146-151 (1990).
2. Tsuchihashi, H., Sasaki, T., Kojima, S., *et al.* Binding of [³H]haloperidol to dopamine D₂ receptors in the rat striatum. *J. Pharm. Pharmacol.* **44(11)**, 911-914 (1992).
3. Lummis, S.C. and Baker, J. Radioligand binding and photoaffinity labelling studies show a direct interaction of phenothiazines at 5-HT₃ receptors. *Neuropharmacology* **36(4-5)**, 665-670 (1997).
4. Ghelardini, C., Galeotti, N., Uslenghi, C., *et al.* Prochlorperazine induces central antinociception mediated by the muscarinic system. *Pharmacol. Res.* **50(3)**, 351-358 (2004).
5. Niemegeers, C.J.E. Antiemetic specificity of dopamine antagonists. *Psychopharmacology (Berl)*. **78(3)**, 210-213 (1982).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/01/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM