

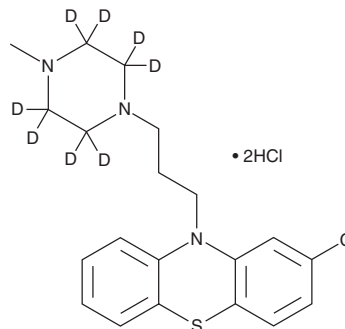
PRODUCT INFORMATION



Prochlorperazine-d₈ (hydrochloride)

Item No. 31916

CAS Registry No.: 2930627-64-6
Formal Name: 2-chloro-10-[3-(4-methyl-1-piperazinyl-d₈) propyl]-10H-phenothiazine, dihydrochloride
MF: C₂₀H₁₆ClD₈N₃S • 2HCl
FW: 454.9
Chemical Purity: ≥98% (Prochlorperazine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
Supplied as: A 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-d₈ (hydrochloride) is intended for use as an internal standard for the quantification of prochlorperazine (Item No. 20742) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Prochlorperazine-d₈ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the prochlorperazine-d₈ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Prochlorperazine-d₈ (hydrochloride) is soluble in DMSO.

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, 07/27/2023

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