

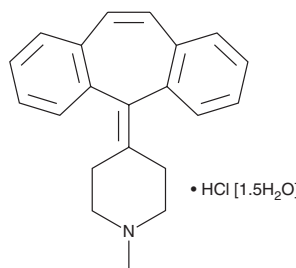
PRODUCT INFORMATION



Cyproheptadine (hydrochloride hydrate)

Item No. 19551

CAS Registry No.: 41354-29-4
Formal Name: 4-(5H-dibenzo[a,d]cyclohepten-5-ylidene)-1-methylpiperidine, monohydrochloride, sesquihydrate
MF: C₂₁H₂₁N • HCl [1.5H₂O]
FW: 350.9
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 287 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

Cyproheptadine (hydrochloride hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the cyproheptadine (hydrochloride hydrate) in the solvent of choice, which should be purged with an inert gas. Cyproheptadine (hydrochloride hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of cyproheptadine (hydrochloride hydrate) in these solvents is approximately 30 mg/ml.

Cyproheptadine (hydrochloride hydrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cyproheptadine (hydrochloride hydrate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Cyproheptadine (hydrochloride hydrate) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cyproheptadine is an antihistamine with antiserotonergic and anticholinergic activities.¹ It binds to histamine H₁, muscarinic, and serotonin 5-HT₂ receptors (K_is = 0.38, 1.26, and 0.83 nM, respectively, in radioligand binding assays).² Cyproheptadine reduces histamine-induced spasms in isolated guinea pig ileum (IC₇₅ = 0.0014 μg/ml).³ It protects against intravenous histamine diphosphate-induced death with a 50% protective dose (PD₅₀) value of 0.08 mg/kg and delays induction of aerosolized histamine diphosphate-induced coughing (ED_{100sec} = 0.29 mg/kg) in guinea pigs. Cyproheptadine also inhibits the lysine methyltransferase SET7/9 (IC₅₀ = 1 μM), decreasing the expression of estrogen receptor α (ERα) in MCF-7 cells.⁴ Formulations containing cyproheptadine have been used in the treatment of allergic reactions including rhinitis, conjunctivitis, and urticaria.

References

1. Greaves, M.W. Antihistamines. *Dermatol. Clin.* **19**(1), 53-62 (2001).
2. Kakiuchi, M., Ohashi, T., Musoh, K., *et al.* Studies on the novel antiallergic agent HSR-609: Its penetration into the central nervous system in mice and guinea pigs and its selectivity for the histamine H₁-receptor. *Jpn. J. Pharmacol.* **73**(4), 291-298 (1997).
3. Lish, P.M., Robbins, S.I., and Peters, E.L. Specificity of antihistamine drugs and involvement of the adrenergic system in histamine deaths in the guinea pig. *J. Pharmacol. Exp. Ther.* **153**(3), 538-543 (1966).
4. Takemoto, Y., Ito, A., Niwa, H., *et al.* Identification of cyproheptadine as an inhibitor of SET domain containing lysine methyltransferase 7/9 (Set7/9) that regulates estrogen-dependent transcription. *J. Med. Chem.* **59**(8), 3650-3660 (2016).

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, 11/29/2022

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