

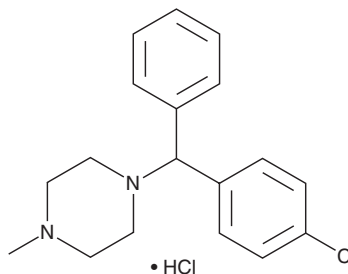
PRODUCT INFORMATION



Chlorcyclizine (hydrochloride)

Item No. 19239

CAS Registry No.: 14362-31-3
Formal Name: 1-[[4-chlorophenyl]phenylmethyl]-4-methyl-piperazine, monohydrochloride
Synonym: NSC 169496
MF: C₁₈H₂₁ClN₂ • HCl
FW: 337.3
Purity: ≥98%
UV/Vis.: λ_{max}: 230 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

Chlorcyclizine (hydrochloride) is supplied as a crystalline solid. Chlorcyclizine (hydrochloride) is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of chlorcyclizine (hydrochloride) be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of chlorcyclizine (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Chlorcyclizine is a phenylpiperazine that acts as a histamine H₁ receptor antagonist (K_i = 9 nM).¹ It has also been shown to be effective against hepatitis C virus (HCV; EC₅₀ = 44 nM *in vitro*), preventing viral entry into host cells.² In chimeric mice engrafted with primary human hepatocytes, 10-50 mg/kg chlorcyclizine significantly inhibited infection of HCV genotypes 1b and 2a.²

References

1. Tran, V.T., Chang, R.S.L., and Snyder, S.H. Histamine H₁ receptors identified in mammalian brain membranes with [³H]mepyramine. *Proc. Natl. Acad. Sci. USA* **75**(12), 6290-6294 (1978).
2. He, S., Xiao, J., Dulcey, A.E., *et al.* Discovery, optimization, and characterization of novel chlorcyclizine derivatives for the treatment of hepatitis C virus infection. *J. Med. Chem.* **59**(3), 841-853 (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, 12/19/2022

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