

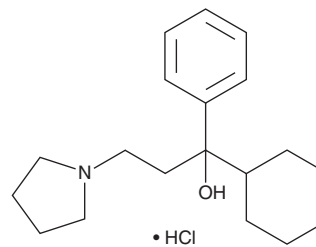
PRODUCT INFORMATION



Procyclidine (hydrochloride)

Item No. 33141

CAS Registry No.: 1508-76-5
Formal Name: α -cyclohexyl- α -phenyl-1-pyrrolidinepropanol, monohydrochloride
MF: $C_{19}H_{29}NO \cdot HCl$
FW: 323.9
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Procyclidine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the procyclidine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Procyclidine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of procyclidine (hydrochloride) in these solvents is approximately 1 mg/ml.

Description

Procyclidine is an antagonist of muscarinic acetylcholine receptors ($K_s = 4.47, 42.66, 7.24, 6.03,$ and 4.79 nM for M_{1-5} receptors, respectively).¹ It increases fine motor activity counts ($ED_{50} = 1.74$ mg/kg) without affecting ambulatory activity in rats in a home-cage activity test.² Formulations containing procyclidine have been used in the treatment of parkinsonism and extrapyramidal dysfunction.

References

1. Lazareno, S., Buckley, N.J., and Roberts, F.F. Characterization of muscarinic M_4 binding sites in rabbit lung, chicken heart, and NG108-15 cell. *Mol. Pharmacol.* **38(6)**, 805-815 (1990).
2. Sipos, M.L., Burchnell, V., and Galbicka, G. Dose-response curves and time-course effects of selected anticholinergics on locomotor activity in rats. *Psychopharmacology* **147(3)**, 250-256 (1999).

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, 10/24/2022

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