

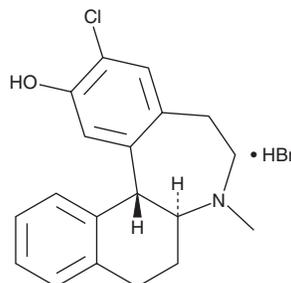
PRODUCT INFORMATION



SCH 39166 (hydrobromide)

Item No. 25331

CAS Registry No.: 1227675-51-5
Formal Name: 11-chloro-7-methyl-6,6 α S,7,8,9,13 β R-hexahydro-5H-benzo[d]naphtho[2,1- β]azepin-12-ol, monohydrobromide
MF: C₁₉H₂₀ClNO • HBr
FW: 394.7
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

SCH 39166 (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the SCH 39166 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. SCH 39166 (hydrobromide) is slightly soluble in DMSO and methanol.

SCH 39166 (hydrobromide) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

SCH 39166 is a dopamine D₁ receptor antagonist (K_i = 5 nM).¹ It is selective for D₁ over D₂₋₄ receptors (K_is = 3,751, >1,000, and 5,934 nM, respectively), however, it also binds to D₅ receptors (K_i = 4.4 nM). SCH 39166 inhibits food intake in a dose-dependent manner in rats (ED₅₀ = 0.84 mg/kg).² It reduces ethanol intake in Sardinian alcohol-preferring rats and sucrose intake in water-deprived and water-sated rats without affecting food or total fluid intake.³ SCH 39166 also suppresses cocaine-induced arrhythmias in anesthetized dogs.⁴

References

1. Tice, M.A., Hashemi, T., Taylor, L.A., *et al.* Characterization of the binding of SCH 39166 to the five cloned dopamine receptor subtypes. *Pharmacol. Biochem. Behav.* **49**(3), 567-571 (1994).
2. Terry, P. and Katz, J.L. A comparison of the effects of the D₁ receptor antagonists SCH 23390 and SCH 39166 on suppression of feeding behavior by the D₁ agonist SKF38393. *Psychopharmacology (Berl)*. **113**(3-4), 328-333 (1994).
3. Panocka, I., Ciccocioppo, R., Mosca, M., *et al.* Effects of the dopamine D₁ receptor antagonist SCH 39166 on the ingestive behaviour of alcohol-preferring rats. *Psychopharmacology (Berl)*. **120**(2), 227-235 (1995).
4. Kanani, P.M., Guse, P.A., Smith, W.M., *et al.* Acute deleterious effects of cocaine on cardiac conduction, hemodynamics, and ventricular fibrillation threshold: Effects of interaction with a selective dopamine D₁ antagonist SCH 39166. *J. Cardiovasc. Pharmacol.* **32**(1), 42-48 (1998).

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, 10/06/2022

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