

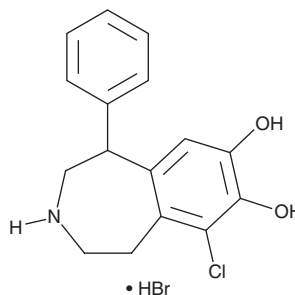
PRODUCT INFORMATION



(±)-SKF 81297 (hydrobromide)

Item No. 15067

CAS Registry No.: 67287-39-2
Formal Name: 6-chloro-2,3,4,5-tetrahydro-1-phenyl-1H-3-benzazepine-7,8-diol, monohydrobromide
Synonym: (±)-6-chloro-PB
MF: C₁₆H₁₆ClNO₂ • HBr
FW: 370.7
Purity: ≥98%
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

(±)-SKF 81297 (hydrobromide) is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-SKF 81297 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. (±)-SKF 81297 (hydrobromide) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)-SKF 81297 (hydrobromide) in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

(±)-SKF 81297 (hydrobromide) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (±)-SKF 81297 (hydrobromide) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (±)-SKF 81297 (hydrobromide) has a solubility of approximately 0.04 mg/ml in a 1:20 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-SKF 81297 is a selective agonist of the dopamine D₁-like receptor (K_i = 1.9 nM).¹ It demonstrates comparatively lower binding affinity for the dopamine D₂, dopamine D₃, serotonin 5-HT_{2A}, and adrenergic α₂ receptors (K_is = 1,272, >10,000, 955, and 509 nM, respectively).¹ Activation of dopamine D₁-like receptors by selective agonists such as (±)-SKF 81297 have been implicated in initiating a cascade of signaling events leading to reward-related incentive learning.²

References

1. Neumeyer, J.L., Kula, N.S., Bergman, J., *et al.* Receptor affinities of dopamine D₁ receptor-selective novel phenylbenzazepines. *Eur. J. Pharmacol.* **474(2-3)**, 137-140 (2003).
2. Beninger, R.J. and Miller, R. Dopamine D₁-like receptors and reward-related incentive learning. *Neurosci. Biobehav. Rev.* **22(2)**, 335-345 (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, 12/01/2022

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