

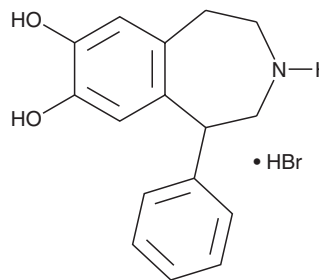
PRODUCT INFORMATION



SKF 38393 (hydrobromide)

Item No. 19150

CAS Registry No.: 20012-10-6
Formal Name: 2,3,4,5-tetrahydro-1-phenyl-1H-3-benzazepine-7,8-diol, monohydrobromide
MF: C₁₆H₁₇NO₂ • HBr
FW: 336.2
Purity: ≥98%
UV/Vis.: λ_{max}: 286 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

SKF 38393 (hydrobromide) is supplied as a crystalline solid. A stock solution may be made by dissolving the SKF 38393 (hydrobromide) in the solvent of choice. SKF 38393 (hydrobromide) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of SKF 38393 (hydrobromide) in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

SKF 38393 (hydrobromide) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SKF 38393 (hydrobromide) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SKF 38393 (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 38393 is a partial agonist of the dopamine D₁-like receptors D₁ and D₅ (K_is = 1 and ~0.5 nM, respectively).¹ It less potently binds D₂, D₃, and D₄ receptors (K_is = ~150, 5,000, and 1,000 nM, respectively).^{1,2}

References

- Seeman, P. and Van Tol, H.H.M. Dopamine receptor pharmacology. *Trends Pharmacol. Sci.* **15**(7), 264-270 (1994).
- 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).

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