

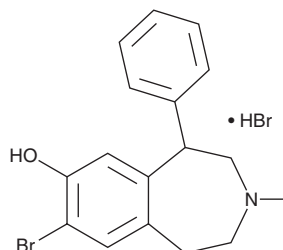
PRODUCT INFORMATION



SKF 83566 (hydrobromide)

Item No. 29480

CAS Registry No.: 108179-91-5
Formal Name: 8-bromo-2,3,4,5-tetrahydro-3-methyl-5-phenyl-1H-3-benzazepin-7-ol, monohydrobromide
MF: C₁₇H₁₈BrNO • HBr
FW: 413.2
Purity: ≥98%
Supplied as: A 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 83566 (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the SKF 83566 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. SKF 83566 (hydrobromide) is soluble in organic solvents such as ethanol and DMSO. It is also soluble in water. The solubility of SKF 83566 (hydrobromide) in ethanol, DMSO, and water is approximately 25, 100, and 5 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

SKF 83566 is a dopamine D₁ receptor antagonist.¹⁻³ *In vivo*, SKF 83566 (0.005, 0.01, and 0.03 mg/kg) reduces dexbenzetimide- and scopolamine-induced locomotor stereotypy in rats.¹ It decreases ghrelin-induced increases in novel object exploration and food intake in rats when administered at a dose of 100 µg/kg.² SKF 83566 also reduces schedule-induced polydipsia in rats.³

References

1. Fritts, M.E., Mueller, K., and Morris, L. Locomotor stereotypy produced by dexbenzetimide and scopolamine is reduced by SKF 83566, not sulpiride. *Pharmacol. Biochem. Behav.* **60(3)**, 639-644 (1998).
2. Jacoby, S.M. and Currie, P.J. SKF 83566 attenuates the effects of ghrelin on performance in the object location memory task. *Neurosci. Lett.* **504(3)**, 316-320 (2011).
3. Mittleman, G., Rosner, A.L., and Schaub, C.L. Polydipsia and dopamine: Behavioral effects of dopamine D1 and D2 receptor agonists and antagonists. *J. Pharmacol. Exp. Ther.* **271(2)**, 638-650 (1994).

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, 11/10/2022

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