

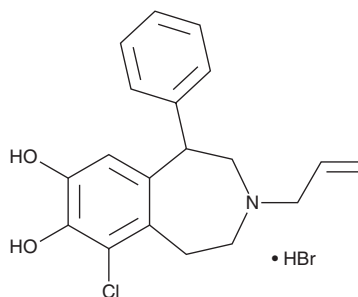
PRODUCT INFORMATION



SKF 82958 (hydrobromide)

Item No. 29033

CAS Registry No.: 74115-01-8
Formal Name: 6-chloro-2,3,4,5-tetrahydro-1-phenyl-3-(2-propen-1-yl)-1H-3-benzazepine-7,8-diol, monohydrobromide
Synonym: (±)-6-chloro-APB
MF: C₁₉H₂₀ClNO₂ • HBr
FW: 410.7
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 82958 (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the SKF 82958 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. SKF 82958 (hydrobromide) is soluble in the organic solvent DMSO and in water (warmed). The solubility of SKF 82958 (hydrobromide) in DMSO and water is approximately 100 mM and 5 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

SKF 82958 is a dopamine D₁ receptor agonist (K_{0.5} = 4 nM in a radioligand binding assay).¹ It is selective for dopamine D₁ over D₂ receptors (K_{0.5} = 73 nM). SKF 82958 induces dopamine D₁ receptor-dependent adenylate cyclase activity in rat striatal membranes (EC₅₀ = 491 nM). SKF 82958 (3 mg/kg) prevents and reverses fentanyl-induced respiratory depression, without affecting analgesia, in cats.² It also increases locomotion and rearing in rats.³

References

1. Mottola, D.M., Laiter, S., Watts, V.J., *et al.* Conformational analysis of D₁ dopamine receptor agonists: Pharmacophore assessment and receptor mapping. *J. Med. Chem.* **39**(1), 285-296 (1996).
2. Lalley, P.M. D₁-dopamine receptor agonists prevent and reverse opiate depression of breathing but not antinociception in the cat. *Am. J. Physiol. Regul. Integr. Comp. Physiol.* **289**(1), R45-R51 (2005).
3. Swanson, C.J., Heath, S., Stratford, T.R., *et al.* Differential behavioral responses to dopaminergic stimulation of nucleus accumbens subregions in the rat. *Pharmacol. Biochem. Behav.* **58**(4), 933-945 (1997).

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, 12/16/2022

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