

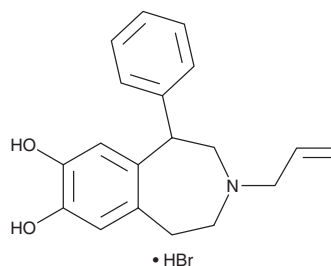
# PRODUCT INFORMATION



## N-Allyl-(±)-SKF 38393 (hydrobromide)

Item No. 32776

**CAS Registry No.:** 300561-58-4  
**Formal Name:** 2,3,4,5-tetrahydro-1-phenyl-3-(2-propen-1-yl)-1H-3-benzazepine-7,8-diol, monohydrobromide  
**MF:** C<sub>19</sub>H<sub>21</sub>NO<sub>2</sub> • HBr  
**FW:** 376.3  
**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

N-Allyl-(±)-SKF 38393 (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the N-allyl-(±)-SKF 38393 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. N-Allyl-(±)-SKF 38393 (hydrobromide) is soluble in organic solvents such as ethanol and DMSO. It is also soluble in water. The solubility of N-allyl-(±)-SKF 38393 (hydrobromide) in these solvents is approximately 10 mM in ethanol and water and 100 mM in DMSO. We do not recommend storing the aqueous solution for more than one day.

### Description

N-Allyl-(±)-SKF 38393 is a dopamine D<sub>1</sub> receptor partial agonist.<sup>1</sup> It is selective for dopamine D<sub>1</sub> over D<sub>2</sub> receptors in radioligand binding assays (K<sub>i</sub>s = 302 and 5,250 nM, respectively). N-Allyl-(±)-SKF 38393 reduces food intake in rats (ED<sub>50</sub> = 3.6 mg/kg).<sup>2</sup> It reduces cocaine and food self-administration in squirrel monkeys when administered at doses ranging from 0.1 to 10 mg/kg.<sup>3</sup>

### References

1. Daly, S.A. and Waddington, J.L. D-1 dopamine receptors and the topography of unconditioned motor behaviour: studies with the selective, 'full efficacy' benzazepine D-1 agonist SKF 83189. *J. Psychopharmacol.* **6**(1), 50-60 (1992).
2. Terry, P. and Katz, J.L. Differential antagonism of the effects of dopamine D1-receptor agonists on feeding behavior in the rat. *Psychopharmacology (Berl)* **109**(4), 403-409 (1992).
3. Platt, D.M., Rowlett, J.K., and Spealman, R.D. Modulation of cocaine and food self-administration by low- and high-efficacy D1 agonists in squirrel monkeys. *Psychopharmacology (Berl)* **157**(2), 208-216 (2001).

#### 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, 11/03/2022

#### 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](http://WWW.CAYMANCHEM.COM)