

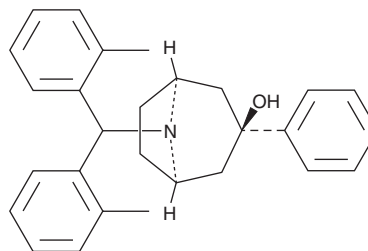
PRODUCT INFORMATION



SCH 221510

Item No. 43181

CAS Registry No.: 322473-89-2
Formal Name: (3-*endo*)-8-[bis(2-methylphenyl)methyl]-3-phenyl-8-azabicyclo[3.2.1]octan-3-ol
MF: C₂₈H₃₁NO
FW: 397.6
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

SCH 221510 is supplied as a solid. A stock solution may be made by dissolving the SCH 221510 in the solvent of choice, which should be purged with an inert gas. SCH 221510 is soluble (≥10 mg/ml) in ethanol. SCH 221510 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SCH 221510 is an agonist of nociceptin opioid peptide (NOP) receptor.¹ It induces [³⁵S]GTPγS binding to CHO membranes expressing recombinant human NOP receptors (EC₅₀ = 12 nM). SCH 221510 (10 mg/kg) increases time spent in the open arms and number of open arm entries in an elevated plus maze in rats. It increases drinking in the Vogel punished drinking task in rats when administered at doses of 3, 10, or 30 mg/kg. SCH 221510 (0.3, 1, or 3 mg/kg) reduces the number of vocalizations in a guinea pig pup model of separation-induced anxiety. It decreases the number of coughs in a guinea pig model of capsaicin-induced coughing (ED₅₀ = 0.07 mg/kg).²

References

1. Varty, G.B., Lu, S.X., Morgan, C.A., *et al.* The anxiolytic-like effects of the novel, orally active nociceptin opioid receptor agonist 8-[bis(2-methylphenyl)methyl]-3-phenyl-8-azabicyclo[3.2.1]octan-3-ol (SCH 221510). *J. Pharmacol. Exp. Ther.* **326**(2), 672-682 (2008).
2. Ho, G.D., Anthes, J., Bercovici, A., *et al.* The discovery of tropane derivatives as nociceptin receptor ligands for the management of cough and anxiety. *Bioorg. Med. Chem. Lett.* **19**(9), 2519-2523 (2009).

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, 04/09/2025

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