

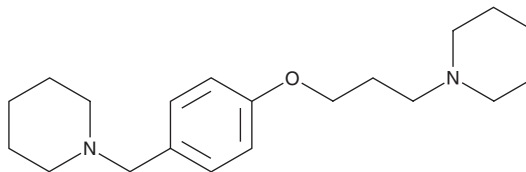
PRODUCT INFORMATION



JNJ-5207852

Item No. 11998

CAS Registry No.: 398473-34-2
Formal Name: 1-[3-[4-(1-piperidinylmethyl)phenoxy]propyl]-piperidine
MF: C₂₀H₃₂N₂O
FW: 316.5
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 275 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

JNJ-5207852 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of JNJ-5207852 in these solvents is approximately 5 and 20 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of JNJ-5207852 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of JNJ-5207852 in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histamine has been implicated in a variety of biological functions including thermo- and immunoregulation, food intake, hyperexcitability, pain transmission, arousal, reward, memory, and emotional responses. The histamine H₃ receptor is a G_i-coupled presynaptic auto- and hetero-receptor whose activation leads to a decreased release of histamine, glutamate, norepinephrine, and acetylcholine in the brain. H₃R antagonists are expected to increase the release of these neurotransmitters, which may contribute therapeutic benefit in diseases characterized by sleep/wake disturbances or cognitive disorders.^{1,2} JNJ-5207852 is a selective, non-imidazole histamine H₃R antagonist with high affinity at the rat (pK_i = 8.9) and human (pK_i = 9.24) H₃ receptors.³ In rodents, 1-10 mg/kg JNJ-5207852 administered subcutaneously increases time spent awake and decreases REM sleep and slow-wave sleep.³

References

1. Tiligada, E., Zampeli, E., Sander, K., *et al.* Histamine H₃ and H₄ receptors as novel drug targets. *Expert Opin. Investig. Drugs* **18**(10), 1519-1532 (2009).
2. Barbier, A.J., Dugovic, B.C., Laposky, A.D., *et al.* Acute wake-promoting actions of JNJ-5207852, a novel, diamine-based H₃ antagonist. *Br. J. Pharmacol.* **143**, 649-661 (2004).
3. Esbenshade, T.A., Browman, K.E., Bitner, R.S., *et al.* The histamine H₃ receptor: An attractive target for the treatment of cognitive disorders. *Br. J. Pharmacol.* **154**, 1166-1181 (2008).

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, 01/29/2019

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