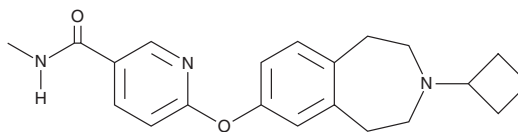


PRODUCT INFORMATION



GSK189254
Item No. 36705

CAS Registry No.: 720690-73-3
Formal Name: 6-[(3-cyclobutyl-2,3,4,5-tetrahydro-1H-3-benzazepin-7-yl)oxy]-N-methyl-3-pyridinecarboxamide
Synonym: GSK189254A
MF: C₂₁H₂₅N₃O₂
FW: 351.4
Purity: ≥98%
UV/Vis.: λ_{max}: 239 nm
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

GSK189254 is supplied as a solid. A stock solution may be made by dissolving the GSK189254 in the solvent of choice, which should be purged with an inert gas. GSK189254 is soluble in DMSO (warmed and sonicated).

Description

GSK189254 is a histamine H₃ receptor antagonist (K_is = 0.13, 0.68, and 1.74 nM for the recombinant human, rat, and mouse receptors, respectively).¹ It is selective for the histamine H₃ receptor over a panel of 50 receptors and ion channels at 1 μM. GSK189254 prevents imetit-induced decreases in forskolin-stimulated cAMP accumulation in HEK293-Gα_o cells expressing the human histamine H₃ receptor (pA₂ = 9.06). It inhibits increases in water intake induced by the histamine H₃ receptor agonist R-(-)-α-methylhistamine (Item No. 25601) and reverses scopolamine-induced memory deficits in the passive avoidance test in rats. GSK189254 increases hindlimb grip force in a rat model of monoiodoacetate-induced osteoarthritic pain and increases the paw withdrawal threshold in a rat model of spinal nerve ligation-induced neuropathic pain (ED₅₀s = 0.77 and 1.5 mg/kg, i.p., respectively).²

References

1. Medhurst, A.D., Atkins, A.R., Beresford, I.J., *et al.* GSK189254, a novel H₃ receptor antagonist that binds to histamine H₃ receptors in Alzheimer's disease brain and improves cognitive performance in preclinical models. *J. Pharmacol. Exp. Ther.* **321**(3), 1032-1045 (2007).
2. Hsieh, G.C., Honore, P., Pai, M., *et al.* Antinociceptive effects of histamine H₃ receptor antagonist in the preclinical models of pain in rats and the involvement of central noradrenergic systems. *Brain Res.* **1354**, 74-84 (2010).

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, 12/21/2022

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