

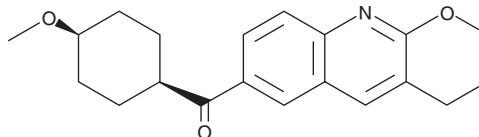
PRODUCT INFORMATION



JNJ-16259685

Item No. 35377

CAS Registry No.: 409345-29-5
Formal Name: (3,4-dihydro-2H-pyrano[2,3-b]quinolin-7-yl)(cis-4-methoxycyclohexyl)-methanone
MF: C₂₀H₂₃NO₃
FW: 325.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

JNJ-16259685 is supplied as a solid. A stock solution may be made by dissolving the JNJ-16259685 in the solvent of choice, which should be purged with an inert gas. JNJ-16259685 is slightly soluble in chloroform and DMSO.

Description

JNJ-16259685 is an antagonist of metabotropic glutamate receptor 1a (mGluR1a; $K_i = 0.34$ nM in CHO(dHFr) cell membranes expressing the rat receptor).¹ It is selective for mGluR1a over mGluR5, -2, -3, -4, and -6, as well as a panel of neurotransmitter receptors, ion channels, and transporters, at 10 μ M. JNJ-16259685 inhibits glutamate-induced inositol phosphate accumulation in primary rat cerebellar neurons with an IC₅₀ value of 1.73 nM. It increases drinking in the Vogel punished drinking task, indicating anxiolytic-like activity, when administered at doses of 2.5, 5, or 10 mg/kg.²

References

1. Lavreysen, H., Wouters, R., Bischoff, F., *et al.* JNJ16259685, a highly potent, selective and systemically active mGlu1 receptor antagonist. *Neuropharmacology* **47(7)**, 961-972 (2004).
2. Steckler, T., Lavreysen, H., Oliveira, A.M., *et al.* Effects of mGlu1 receptor blockade on anxiety-related behaviour in the rat lick suppression test. *Psychopharmacology (Berl.)* **179(1)**, 198-206 (2005).

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, 10/08/2021

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