

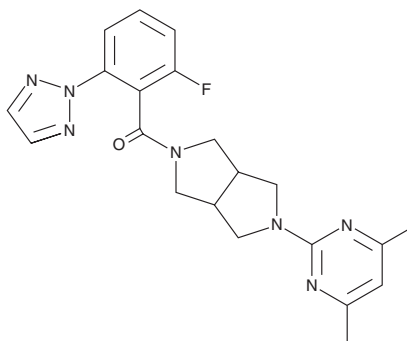
PRODUCT INFORMATION



JNJ-42847922

Item No. 41854

CAS Registry No.: 1293281-49-8
Formal Name: [5-(4,6-dimethyl-2-pyrimidinyl)hexahydropyrrolo[3,4-c]pyrrol-2(1H)-yl][2-fluoro-6-(2H-1,2,3-triazol-2-yl)phenyl]-methanone
Synonym: MIN-202
MF: C₂₁H₂₂FN₇O
FW: 407.5
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

JNJ-42847922 is supplied as a solid. A stock solution may be made by dissolving the JNJ-42847922 in the solvent of choice, which should be purged with an inert gas. JNJ-42847922 is slightly soluble (0.1-1 mg/ml) in DMSO.

JNJ-42847922 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

JNJ-42847922 is an orexin receptor 2 (OX2R) antagonist ($K_i = 10$ nM for the human receptor).¹ It is selective for OX2R over OX1R ($K_i = 794$ nM for the human receptor) and a panel of 50 receptors, ion channels, and transporters at 1 μ M. JNJ-42847922 (3, 10, and 30 mg/kg) reduces the latency to non-rapid eye movement (NREM) sleep onset and increases the duration of NREM sleep in rats.

Reference

1. Bonaventure, P., Shelton, J., Yun, S., *et al.* Characterization of JNJ-42847922, a selective orexin-2 receptor antagonist, as a clinical candidate for the treatment of insomnia. *J. Pharmacol. Exp. Ther.* **354**(3), 471-482 (2015).

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, 10/29/2024

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