

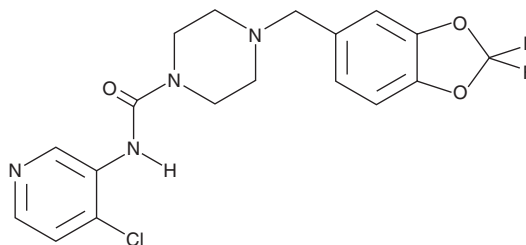
PRODUCT INFORMATION



JNJ-42165279

Item No. 19987

CAS Registry No.: 1346528-50-4
Formal Name: N-(4-chloro-3-pyridinyl)-4-[(2,2-difluoro-1,3-benzodioxol-5-yl)methyl]-1-piperazinecarboxamide
MF: C₁₈H₁₇ClF₂N₄O₃
FW: 410.8
Purity: ≥98%
UV/Vis.: λ_{max}: 241, 272 nm
Supplied as: A crystalline 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-42165279 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-42165279 in the solvent of choice, which should be purged with an inert gas. JNJ-42165279 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of JNJ-42165279 in these solvents is approximately 30 mg/ml.

JNJ-42165279 is sparingly soluble 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-42165279 is a potent, irreversible inhibitor of fatty acid amide hydrolase (FAAH; IC₅₀s = 70 and 313 nM for human and rat forms, respectively).¹ It displays selectivity for FAAH over a panel of other enzymes, receptors, transporters, and ion channels. JNJ-42165279 is active *in vivo*, blocking FAAH activity in brain and periphery of rats and raising concentrations of anandamide (Item No. 90050), oleoyl ethanolamide (Item No. 90265), and palmitoyl ethanolamide (Item No. 90350).¹ It is also efficacious in the spinal nerve ligation model of neuropathic pain.¹

Reference

1. Keith, J. M., Jones, W. M., Tichenor, M. *et al.* Preclinical characterization of the FAAH inhibitor JNJ-42165279. *ACS Med.Chem.Lett.* **6**(12), 1204-1208 (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, 11/21/2022

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