

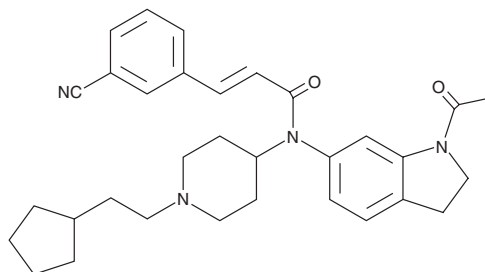
PRODUCT INFORMATION



JNJ-5207787

Item No. 13947

CAS Registry No.: 683746-68-1
Formal Name: (2E)-N-(1-acetyl-2,3-dihydro-1H-indol-6-yl)-3-(3-cyanophenyl)-N-[1-(2-cyclopentylethyl)-4-piperidinyl]-2-propenamide
MF: C₃₂H₃₈N₄O₂
FW: 510.7
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 262 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-5207787 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-5207787 in the solvent of choice. JNJ-5207787 is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of JNJ-5207787 in these solvents is approximately 5 mg/ml, with gentle warming required.

Description

JNJ-5207787 is an antagonist of the neuropeptide Y (NPY) receptor Y₂ that blocks peptide YY binding (IC₅₀ = 100 nM).^{1,2} It displays 100-fold selectivity for Y₂ over other NPY receptors and has minimal effect on a panel of 50 receptors, ion channels, and transporters, except for sodium channel 2.¹ JNJ-5207787 is intraperitoneally bioavailable and brain penetrant in rats, occupying maximally 50% of Y₂ receptor binding sites in the hypothalamic area.¹

References

1. Bonaventure, P., Nepomuceno, D., Mazur, C., *et al.* Characterization of N-(1-Acetyl-2,3-dihydro-1H-indol-6-yl)-3-(3-cyano-phenyl)-N-[1-(2-cyclopentyl-ethyl)-piperidin-4yl]acrylamide (JNJ-5207787), a small molecule antagonist of the neuropeptide Y Y₂ receptor. *J. Pharm. Exp. Ther.* **308**(3), 1130-1137 (2004).
2. Mittapalli, G.K. and Roberts, E. Ligands of the neuropeptide Y Y₂ receptor. *Bioorg. Med. Chem. Lett.* **24**(2), 430-441 (2014).

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, 11/17/2022

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