

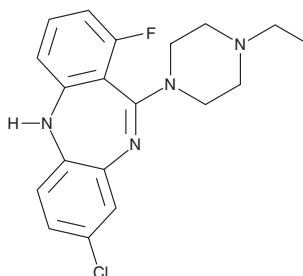
PRODUCT INFORMATION



JHU37152

Item No. 39532

CAS Registry No.: 2369979-67-7
Formal Name: 8-chloro-11-(4-ethyl-1-piperazinyl)-1-fluoro-5H-dibenzo[b,e][1,4]diazepine
MF: C₁₉H₂₀ClFN₄
FW: 358.8
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

JHU37152 is supplied as a solid. A stock solution may be made by dissolving the JHU37152 in the solvent of choice, which should be purged with an inert gas. JHU37152 is soluble in methanol and DMSO.

Description

JHU37152 is an activator of designer receptors exclusively activated by designer drugs (DREADDs) derived from human muscarinic acetylcholine receptors (mAChRs).¹ It binds to hM3Dq and hM4Di receptors (K_s = 1.8 and 8.7 nM, respectively) and induces G_{α_{o1}} activation in HEK293 cells expressing hM3Dq or hM4Di (EC₅₀s = 5 and 0.5 nM, respectively). *In vivo*, JHU37152 (1 mg/kg) inhibits locomotor activity in D1-hM3Dq and D1-hM4Di transgenic mice.

Reference

1. Bonaventura, J., Eldridge, M.A.G., Hu, F., *et al.* High-potency ligands for DREADD imaging and activation in rodents and monkeys. *Nat. Commun.* **10**(1), 4627 (2019).

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, 01/25/2024

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