

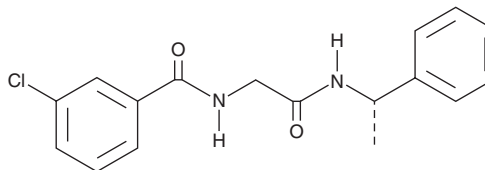
PRODUCT INFORMATION



JNJ-63533054

Item No. 23450

CAS Registry No.: 1802326-66-4
Formal Name: 3-chloro-N-[2-oxo-2-[[[(1S)-1-phenylethyl]amino]ethyl]-benzamide
MF: C₁₇H₁₇ClN₂O₂
FW: 316.8
Purity: ≥98%
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-63533054 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-63533054 in the solvent of choice, which should be purged with an inert gas. JNJ-63533054 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JNJ-63533054 in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

JNJ-63533054 is a potent agonist of the orphan G protein-coupled receptor GPR139 (EC₅₀ = 16 nM in HEK293 cells expressing human GPR139).¹ It is selective for GPR139 over 50 known G protein-coupled receptors, ion channels, and transporters as determined by an external selectivity panel. Tritium-labeled JNJ-63533054 has been used in displacement assays to identify physiological ligands of GPR139.²

References

1. Dvorak, C.A., Coate, H., Nepomuceno, D., *et al.* Identification and SAR of glycine benzamides as potent agonists for the GPR139 receptor. *ACS Med. Chem. Lett.* **6**(9), 1015-1018 (2015).
2. Liu, C., Bonaventure, P., Lee, G., *et al.* GPR139, an orphan receptor highly enriched in the habenula and septum, is activated by the essential amino acids L-Tryptophan and L-Phenylalanine. *Mol. Pharmacol.* **88**(5), 911-925 (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, 11/08/2022

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