

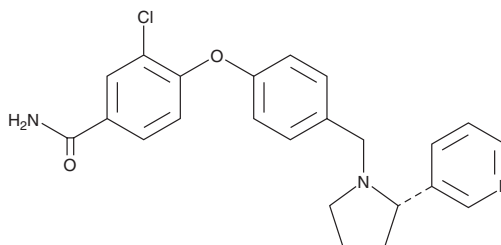
PRODUCT INFORMATION



LY2795050

Item No. 33604

CAS Registry No.: 1346133-08-1
Formal Name: 3-chloro-4-[4-[[[(2S)-2-(3-pyridinyl)-1-pyrrolidinyl]methyl]phenoxy]-benzamide
MF: C₂₃H₂₂ClN₃O₂
FW: 407.9
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

LY2795050 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2795050 in the solvent of choice, which should be purged with an inert gas. LY2795050 is soluble in the organic solvent DMSO.

Description

LY2795050 is a κ -opioid receptor antagonist ($K_i = 0.722$ nM).¹ It is selective for κ -opioid receptors over μ - and δ -opioid receptors ($K_{iS} = 25.8$ and 153 nM, respectively). LY2795050 (0.1 mg/kg, i.p.) reduces immobility time in the forced swim test, as well as reduces salvinorin A-induced decreases in grooming time in the splash test, indicating antidepressant-like activity, in mice.²

References

1. Mitch, C.H., Quimby, S.J., Diaz, N., *et al.* Discovery of aminobenzyloxyarylamides as κ opioid receptor selective antagonists: Application to preclinical development of a κ opioid receptor antagonist receptor occupancy tracer. *J. Med. Chem.* **54**(23), 8000-8012 (2011).
2. Butelman, E.R., McElroy, B.D., Prisinzano, T.E., *et al.* Impact of pharmacological manipulation of the κ -opioid receptor system on self-grooming and anhedonic-like behaviors in male mice. *J. Pharmacol. Exp. Ther.* **370**(1), 1-8 (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, 10/20/2022

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