

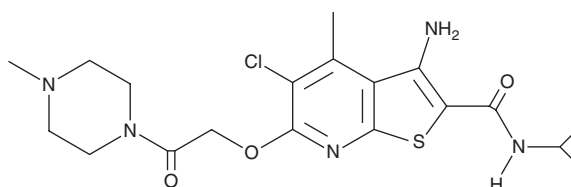
PRODUCT INFORMATION



LY2119620

Item No. 26373

CAS Registry No.: 886047-22-9
Formal Name: 3-amino-5-chloro-N-cyclopropyl-4-methyl-6-[2-(4-methyl-1-piperazinyl)-2-oxoethoxy]-thieno[2,3-b]pyridine-2-carboxamide
MF: C₁₉H₂₄ClN₅O₃S
FW: 437.9
Purity: ≥98%
UV/Vis.: λ_{max}: 291 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

LY2119620 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2119620 in the solvent of choice, which should be purged with an inert gas. LY2119620 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LY2119620 in ethanol is approximately 1 mg/ml and approximately 50 mg/ml in DMSO and DMF.

LY2119620 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY2119620 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LY2119620 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LY2119620 is a positive allosteric modulator of the M₂ and M₄ muscarinic acetylcholine receptors (mAChRs).¹ It binds to allosteric sites on M₂ and M₄ receptors with K_B values ranging from 1.8 to 3.4 μM. LY2119620 potentiates the activity of the AChR agonist acetylcholine (ACh; Item No. 23829) and the mAChR agonists iperoxo and oxotremorine M (Item No. 20847) at human recombinant M₂ and M₄ receptors expressed in CHO cell membranes. LY2119620 also exhibits allosteric agonism at M₂ and M₄ receptors (23.2 and 16.8%, respectively, of the maximal ACh response).

Reference

1. Croy, C.H., Schober, D.A., Xiao, H., *et al.* Characterization of the novel positive allosteric modulator, LY2119620, at the muscarinic M₂ and M₄ receptors. *Mol. Pharmacol.* **86(1)**, 106-115 (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, 11/28/2022

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