

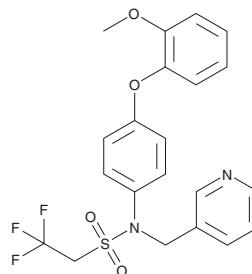
PRODUCT INFORMATION



LY487379

Item No. 45330

CAS Registry No.: 353231-17-1
Formal Name: 2,2,2-trifluoro-N-[4-(2-methoxyphenoxy)phenyl]-N-(3-pyridinylmethyl)ethanesulfonamide
MF: C₂₁H₁₉F₃N₂O₄S
FW: 452.4
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

LY487379 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY487379 in the solvent of choice, which should be purged with an inert gas. LY487379 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

LY487379 is a positive allosteric modulator of metabotropic glutamate receptor 2 (mGluR2).¹ It is selective for mGluR2 over mGluR3 in cells expressing recombinant mGluR2 or mGluR3 (IC₅₀s = 1.7 and >10 μM, respectively). LY487379 reduces MK-801-induced superoxide dismutase (SOD) activity in mouse frontal cortex and hippocampal homogenates.² *In vivo*, LY487379 (30 mg/kg) increases cognitive flexibility in the attentional set-shifting task and decreases impulsive behavior in the differential reinforcement of low rate 72s (DRL72) schedule of food reinforcement in rats.³ It also decreases the severity of dyskinesia and psychotic-like behaviors induced by L-DOPA (Item No. 13248) in a 6-MPTP-lesioned marmoset model of Parkinson's disease when administered at a dose of 1 or 10 mg/kg.⁴

References

1. Burnat, G., Santocki, M., Kalinowski, L., *et al.* The impact of muscarinic and mGlu receptors modulators on MK-801-induced impairments in NO-dependent processes both in vitro and in vivo. *Pharmacol. Rep.* **77(5)**, 1309-1322 (2025).
2. Nikiforuk, A., Popik, P., Drescher, K.U., *et al.* Effects of a positive allosteric modulator of group II metabotropic glutamate receptors, LY487379, on cognitive flexibility and impulsive-like responding in rats. *J. Pharmacol. Exp. Ther.* **335(3)**, 665-673 (2010).
3. Sid-Otmane, L., Hamadjida, A., Nuara, S.G., *et al.* Selective metabotropic glutamate receptor 2 positive allosteric modulation alleviates L-DOPA-induced psychosis-like behaviours and dyskinesia in the MPTP-lesioned marmoset. *Eur. J. Pharmacol.* **873**, 172957 (2020).
4. Schaffhauser, H., Rowe, B.A., Morales, S., *et al.* Pharmacological characterization and identification of amino acids involved in the positive modulation of metabotropic glutamate receptor subtype 2. *Mol. Pharmacol.* **64(4)**, 798-810 (2003).

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, 03/24/2026

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