

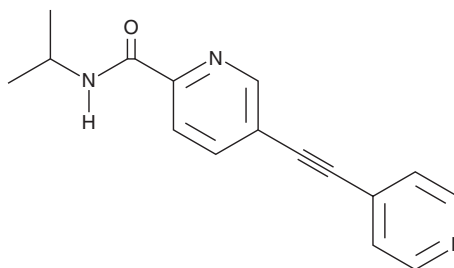
PRODUCT INFORMATION



LSN2463359

Item No. 22104

CAS Registry No.: 1401031-52-4
Formal Name: N-(1-methylethyl)-5-[2-(4-pyridinyl)ethynyl]-2-pyridinecarboxamide
MF: C₁₆H₁₅N₃O
FW: 265.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 247, 297 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

LSN2463359 is supplied as a crystalline solid. A stock solution may be made by dissolving the LSN2463359 in the solvent of choice. LSN2463359 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of LSN2463359 in these solvents is approximately 30 mg/ml.

Description

LSN2463359 is a positive allosteric modulator of the metabotropic glutamate receptor 5 (mGluR5; EC₅₀ = 24 nM).¹ It is without effect at other mGluRs. LSN2463359 is brain penetrant and reverses learning deficits in a rat model of schizophrenia.^{1,2} Through its effects on mGluR5, LSN2463359 attenuates deficits in performance in operant behavior induced by SDZ-220581, an NMDA (Item No. 14581) receptor antagonist, in rats.^{2,3} It also promotes wakefulness in animals.²

References

1. Gastambide, F., Cotel, M.-C., Gilmour, G., *et al.* Selective remediation of reversal learning deficits in the neurodevelopmental MAM model of schizophrenia by a novel mGlu5 positive allosteric modulator. *Neuropsychopharmacology* **37(4)**, 1057-1066 (2012).
2. Gilmour, G., Broad, L.M., Wafford, K.A., *et al.* *In vitro* characterisation of the novel positive allosteric modulators of the mGlu₅ receptor, LSN2463359 and LSN2814617, and their effects on sleep architecture and operant responding in the rat. *Neuropharmacology* **64**, 224-239 (2013).
3. Gastambide, F., Gilmour, G., Robbins, T.W., *et al.* The mGlu₅ positive allosteric modulator LSN2463359 differentially modulates motor, instrumental and cognitive effects of NMDA receptor antagonists in the rat. *Neuropharmacology* **64**, 240-247 (2013).

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, 12/02/2022

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