

PRODUCT INFORMATION



LSN2814617

Item No. 33607

CAS Registry No.: 1313498-08-6

Formal Name: 3-(1,1-dimethylethyl)-7-[3-(4-fluorophenyl)-1,2,4-oxadiazol-5-yl]-5,6,7,8-tetrahydro-1,2,4-triazolo[4,3-a]pyridine

MF: C₁₈H₂₀FN₅O

FW: 341.4

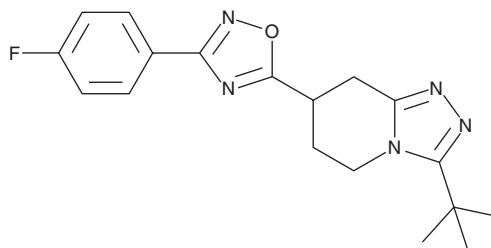
Purity: ≥98%

UV/Vis.: λ_{max}: 240 nm

Supplied as: A 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

LSN2814617 is supplied as a solid. A stock solution may be made by dissolving the LSN2814617 in the solvent of choice, which should be purged with an inert gas. LSN2814617 is slightly soluble in ethanol, DMSO, and dimethyl formamide.

Description

LSN2814617 is a positive allosteric modulator of metabotropic glutamate receptor 5 (mGluR5).¹ It potentiates glutamate-induced calcium mobilization in MCB3901 cells expressing human mGluR5 (EC₅₀ = 52 nM). LSN2814617 increases rat hippocampal mGluR5 receptor occupancy (ED₅₀ = 13 mg/kg). It increases wakefulness and decreases non-rapid eye movement (NREM) and REM sleep in rats when administered at doses of 0.3, 1, and 3 mg/kg.

Reference

1. 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).

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, 11/15/2022

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