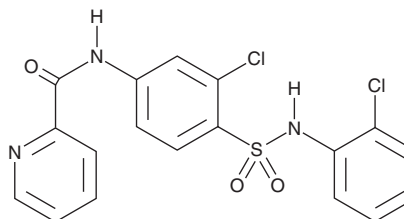


PRODUCT INFORMATION



VU0364439
Item No. 16296

CAS Registry No.: 1246086-78-1
Formal Name: N-[3-chloro-4-[[[(2-chlorophenyl)amino]sulfonyl]phenyl]-2-pyridinecarboxamide
MF: C₁₈H₁₃Cl₂N₃O₃S
FW: 422.3
Purity: ≥98%
UV/Vis.: λ_{max}: 280 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

VU0364439 is supplied as a crystalline solid. A stock solution may be made by dissolving the VU0364439 in the solvent of choice, which should be purged with an inert gas. VU0364439 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of VU0364439 in these solvents is approximately 3 mg/ml.

VU0364439 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VU0364439 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. VU0364439 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VU0364439 is a positive allosteric modulator of metabotropic glutamate receptor 4 (mGluR4) that potentiates glutamate responses in CHO cells expressing recombinant human mGluR4 (EC₅₀ = 19.8 nM).¹

Reference

- Engers, D.W., Gentry, P.R., Williams, R., *et al.* Synthesis and SAR of novel, 4-(phenylsulfamoyl) phenylacetamide mGlu₄ positive allosteric modulators (PAMs) identified by functional high-throughput screening (HTS). *Bioorg. Med. Chem. Lett.* **20(17)**, 5175-5178 (2010).

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, 12/14/2022

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