

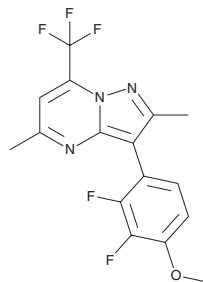
PRODUCT INFORMATION



VU6005649

Item No. 43138

CAS Registry No.: 2137047-43-7
Formal Name: 3-(2,3-difluoro-4-methoxyphenyl)-
2,5-dimethyl-7-(trifluoromethyl)-
pyrazolo[1,5-a]pyrimidine
MF: C₁₆H₁₂F₅N₃O
FW: 357.3
Purity: ≥98%
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

VU6005649 is supplied as a solid. A stock solution may be made by dissolving the VU6005649 in the solvent of choice, which should be purged with an inert gas. VU6005649 is soluble (≥10 mg/ml) in DMSO. VU6005649 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

VU6005649 is a positive allosteric modulator (PAM) of metabotropic glutamate receptor 7 (mGluR7) and mGluR8.¹ It enhances calcium mobilization induced by L-AP4 (Item No. 14538) in HEK cells expressing mGluR7 or mGluR8 (EC₅₀s = 0.65 and 2.6 μM, respectively). It is selective for mGluR7 and mGluR8 over mGluR1-6, as well as a panel of 67 GPCRs, ion channels, transports, and nuclear hormone receptors, at 10 μM but does inhibit the neurokinin-1 (NK₁) receptor (IC₅₀ = 3.4 μM). *In vivo*, VU6005649 (50 mg/kg, i.p) increases the length of freezing time induced by contextual fear conditioning in mice.

Reference

1. Abe, M., Seto, M., Gogliotti, R.G., *et al.* Discovery of VU6005649, a CNS penetrant mGlu_{7/8} receptor PAM derived from a series of pyrazolo[1,5- a]pyrimidines. *ACS Med. Chem. Lett.* **8**(10), 1110-1115 (2017).

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, 04/02/2025

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