

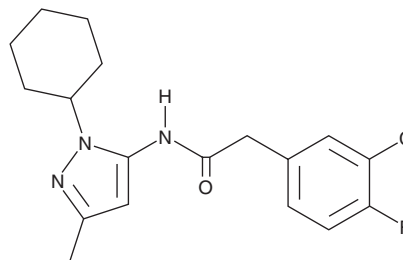
# PRODUCT INFORMATION



**VU0810464**

Item No. 43073

**CAS Registry No.:** 2126040-21-7  
**Formal Name:** 3-chloro-N-(1-cyclohexyl-3-methyl-1H-pyrazol-5-yl)-4-fluoro-benzeneacetamide  
**MF:** C<sub>18</sub>H<sub>21</sub>ClFN<sub>3</sub>O  
**FW:** 349.8  
**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

VU0810464 is supplied as a solid. A stock solution may be made by dissolving the VU0810464 in the solvent of choice, which should be purged with an inert gas. VU0810464 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

## Description

VU0810464 is an activator of G protein-activated inward rectifier potassium channels (GIRKs), also known as K<sub>ir</sub>3.<sup>1</sup> It induces GIRK1/2 and GIRK1/4 activity in a thallium flux assay using HEK293 cells expressing GIRK1 and GIRK2 or GIRK1 and GIRK4 (EC<sub>50</sub>s = 165 and 720 nM, respectively). VU0810464 (10 and 30 mg/kg, i.p.) reduces stress-induced hyperthermia in mice, but it does not affect the percentage of time spent or distance traveled in the open arms of the elevated plus maze in mice at 30 mg/kg.<sup>2</sup>

## References

1. Wieting, J.M., Vadukoot, A.K., Sharma, S., *et al.* Discovery and characterization of 1H-pyrazol-5-yl-2-phenylacetamides as novel, non-urea-containing GIRK1/2 potassium channel activators. *ACS Chem. Neurosci.* **8**(9), 1873-1879 (2017).
2. Vo, B.N., Abney, K.K., Anderson, A., *et al.* VU0810464, a non-urea G protein-gated inwardly rectifying K<sup>+</sup> (K<sub>ir</sub>3/GIRK) channel activator, exhibits enhanced selectivity for neuronal K<sub>ir</sub>3 channels and reduces stress-induced hyperthermia in mice. *Br. J. Pharmacol.* **176**(13), 2238-2249 (2019).

### 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, 05/08/2025

## 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