

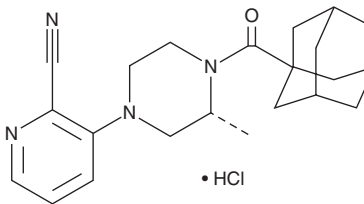
PRODUCT INFORMATION



VU0469650 (hydrochloride)

Item No. 32933

Formal Name: 3-[(3R)-3-methyl-4-(tricyclo[3.3.1.1^{3,7}]-dec-1-ylcarbonyl)-1-piperazinyl]-2-pyridinecarbonitrile, monohydrochloride
MF: C₂₂H₂₈N₄O • HCl
FW: 401.0
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

VU0469650 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the VU0469650 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. VU0469650 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of VU0469650 (hydrochloride) in these solvents is approximately 50 and 100 mM, respectively.

Description

VU0469650 is a brain-penetrant negative allosteric modulator of metabotropic glutamate receptor 1 (mGluR1).¹ It inhibits glutamate-induced calcium flux in cells expressing human mGluR₁ (IC₅₀ = 99 nM). VU0469650 is greater than 100-fold selective for mGluR1 over mGluR2-8 at 10 μM.

Reference

1. Lovell, K.M., Felts, A.S., Rodriguez, A.L., et al. N-Acyl-N'-arylpiperazines as negative allosteric modulators of mGlu₁: Identification of VU0469650, a potent and selective tool compound with CNS exposure in rats. *Bioorg. Med. Chem. Lett.* **23**(13), 3713-3718 (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, 07/27/2023

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