

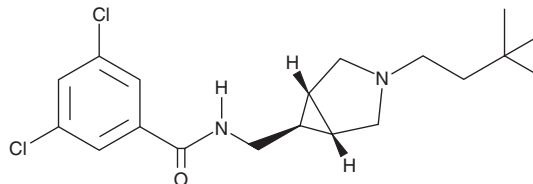
PRODUCT INFORMATION



ML-218

Item No. 43541

CAS Registry No.: 1346233-68-8
Formal Name: 3,5-dichloro-N-[[[(1 α ,5 α ,6 α)-3-(3,3-dimethylbutyl)-3-azabicyclo[3.1.0]hex-6-yl]methyl]-benzamide
MF: C₁₉H₂₆Cl₂N₂O
FW: 369.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ML-218 is supplied as a solid. A stock solution may be made by dissolving the ML-218 in the solvent of choice, which should be purged with an inert gas. ML-218 is sparingly soluble (1-10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

Description

ML-218 is an inhibitor of T-type voltage-gated calcium channel 3.2 (Ca_v3.2) and Ca_v3.3.¹ It inhibits currents in whole-cell patch clamp assays using HEK293 cells expressing Ca_v3.2 or Ca_v3.3 (IC₅₀s = 310 and 274 nM, respectively). ML-218 (0.1-30 mg/kg) decreases the latency to paw withdrawal in a rat model of catalepsy induced by haloperidol (Item No. 12014).

Reference

- Xiang, Z., Thompson, A.D., Brogan, J.T., *et al.* The discovery and characterization of ML218: A novel, centrally active T-type calcium channel inhibitor with robust effects in STN neurons and in a rodent model of Parkinson's Disease. *ACS Chem. Neurosci.* **2(12)**, 730-742 (2011).

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, 07/17/2025

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