

PRODUCT INFORMATION



ML-402

Item No. 42962

CAS Registry No.: 298684-44-3

Formal Name: N-[2-(4-chloro-2-methylphenoxy)ethyl]-2-thiophenecarboxamide

MF: C₁₄H₁₄ClNO₂S

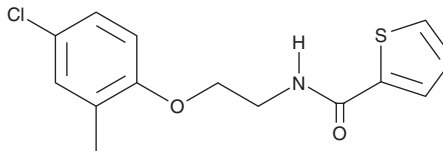
FW: 295.8

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

ML-402 is supplied as a solid. A stock solution may be made by dissolving the ML-402 in the solvent of choice, which should be purged with an inert gas. ML-402 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

ML-402 is an activator of the two-pore domain potassium channels K_{2p}2.1/TREK-1 and K_{2p}10.1/TREK-2.¹ It selectively activates K_{2p}2.1/TREK-1 and K_{2p}10.1/TREK-2 expressed in *Xenopus* oocytes (EC₅₀s = 13.7 and 5.9 μM, respectively) over K_{2p}4.1/TRAAC expressed in *Xenopus* oocytes at 100 μM

Reference

1. Lolicato, M., Arrigoni, C., Mori, T., *et al.* K_{2p}2.1(TREK-1): Activator complexes reveal a cryptic selectivity filter binding site. *Nature* **547(7663)**, 364-368 (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, 03/24/2025

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