

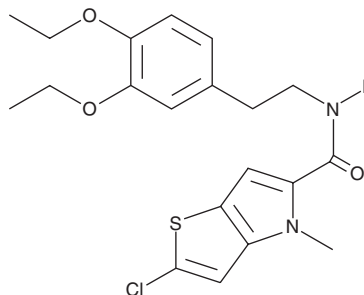
PRODUCT INFORMATION



ML-261

Item No. 15192

CAS Registry No.: 902523-58-4
Formal Name: 2-chloro-N-[2-(3,4-diethoxyphenyl)ethyl]-4-methyl-4H-thieno[3,2-b]pyrrole-5-carboxamide
Synonym: CID-9550710
MF: C₂₀H₂₃ClN₂O₃S
FW: 406.9
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 297 nm
Supplied as: A crystalline 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-261 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-261 in the solvent of choice, which should be purged with an inert gas. ML-261 is soluble in the organic solvent DMSO, at a concentration of approximately 5 mg/ml.

Description

ML-261 is a thienopyrrole-5-carboxamide that inhibits lipid droplet formation in murine AML12 hepatocytes (IC₅₀ = 69.7 nM) but not in human Huh7 hepatocellular carcinoma or primary hepatocyte cells.¹

Reference

1. Zou, J., Ganji, S., Pass, I., et al. Potent inhibitors of lipid droplet formation. *Probe Reports from the NIH Molecular Libraries Program*. (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, 12/13/2022

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