

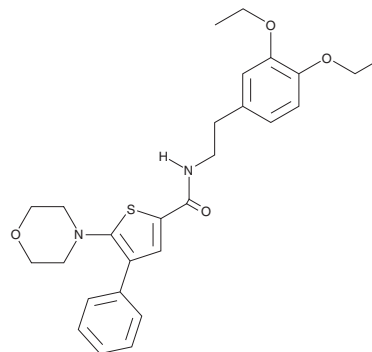
PRODUCT INFORMATION



ML-262

Item No. 21902

CAS Registry No.: 902502-82-3
Formal Name: N-[2-(3,4-diethoxyphenyl)ethyl]-5-(4-morpholinyl)-4-phenyl-2-thiophenecarboxamide
MF: C₂₇H₃₂N₂O₄S
FW: 480.6
Purity: ≥95%
UV/Vis.: λ_{max}: 228, 256, 328 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-262 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-262 in the solvent of choice. ML-262 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of ML-262 in these solvents is approximately 25 and 12.5 mg/ml, respectively.

ML-262 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-262 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ML-262 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ML-262 is an inhibitor of hepatic lipid droplet formation (IC₅₀ = 6.4 nM in murine AML-12 cells), which is associated with non-alcoholic fatty liver disease.¹ ML-262 does not induce cytotoxicity (up to 33 μM) or inhibit fatty acid uptake (up to 50 μM).

Reference

1. Zou, J., Ganji, S., Pass, I., *et al.* Potent inhibitors of lipid droplet formation. In Probe Reports from the NIH Molecular Libraries Program [Internet]. 1-17 (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, 09/29/2022

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