

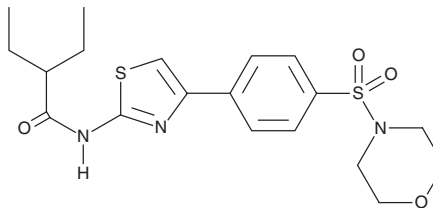
PRODUCT INFORMATION



ML-356

Item No. 20940

CAS Registry No.: 1808260-45-8
Formal Name: 2-ethyl-N-[4-[4-(4-morpholinylsulfonyl)phenyl]-2-thiazolyl]-butanamide
MF: C₁₉H₂₅N₃O₄S₂
FW: 423.5
Purity: ≥98%
UV/Vis.: λ_{max}: 240, 292 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-356 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-356 in the solvent of choice, which should be purged with an inert gas. ML-356 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ML-356 in these solvents is approximately 0.5, 20, and 10 mg/ml, respectively.

ML-356 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-356 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ML-356 has a solubility of approximately 0.09 mg/ml in a 1:10 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-356 is an inhibitor of the thioesterase domain of fatty acid synthase (FASN-TE; IC₅₀ = 334 nM).¹ It is selective for FASN over a number of other human thioesterases expressed in tumor cells, and selective for FASN-TE over its most closely related human homolog, ACOT4.¹ In cells, ML-356 has been shown to block the biosynthesis of palmitate, the end product of FASN.¹

Reference

1. Ardecky, R., Hedrick, M.P., Gosalia, P., *et al.* Selective inhibitors of FAS-TE. Probe Reports from the NIH Molecular Libraries Program, (2010).

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, 10/31/2022

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