

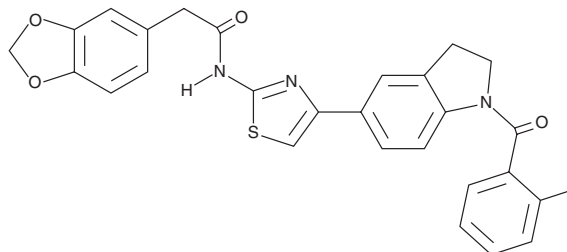
PRODUCT INFORMATION



ML-385

Item No. 21114

CAS Registry No.: 846557-71-9
Formal Name: N-[4-[2,3-dihydro-1-(2-methylbenzoyl)-1H-indol-5-yl]-5-methyl-2-thiazolyl]-1,3-benzodioxole-5-acetamide
MF: C₂₉H₂₅N₃O₄S
FW: 511.6
Purity: ≥98%
UV/Vis.: λ_{max}: 288 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-385 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-385 in the solvent of choice, which should be purged with an inert gas. ML-385 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ML-385 in these solvents is approximately 25 and 30 mg/ml, respectively.

ML-385 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-385 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ML-385 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ML-385 is an inhibitor of nuclear factor erythroid 2-related factor 2 (Nrf2; IC₅₀ = 1.9 μM) that blocks downstream target gene expression.¹ It substantially enhances cytotoxicity in non-small cell lung cancer (NSCLC) cells treated with platinum-based drugs.¹ ML-385 shows specificity and selectivity for NSCLC cells with KEAP1 mutation, which leads to gain of Nrf2 function.¹

Reference

1. Singh, A., Venkannagari, S., Oh, K.H., *et al.* Small molecule inhibitor of NRF2 selectively intervenes therapeutic resistance in KEAP1-deficient NSCLC tumors. *ACS Chem Biol.* **11**(11), 3214-3225 (2016).

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, 11/21/2022

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