

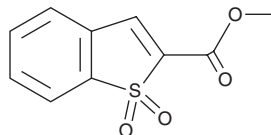
PRODUCT INFORMATION



FX12

Item No. 42089

CAS Registry No.: 1807639-36-6
Formal Name: 1,1-dioxide, benzo[b]thiophene-2-carboxylic acid, methyl ester
MF: C₁₀H₈O₄S
FW: 224.2
Purity: ≥95%
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

FX12 is supplied as a solid. A stock solution may be made by dissolving the FX12 in the solvent of choice, which should be purged with an inert gas. FX12 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

FX12 is an inhibitor and degrader of the E3 ubiquitin ligase RNF5.¹ It binds to RNF5 and inhibits its E3 ligase activity in cell-free assays. FX12 (1.25-20 μM) induces proteasomal degradation of RNF5 via the endoplasmic reticulum-associated proteasomal degradation (ERAD) pathway without modifying mRNA levels in HeLa cells.

Reference

1. Ruan, J., Liang, D., Yan, W., *et al.* A small-molecule inhibitor and degrader of the RNF5 ubiquitin ligase. *Mol. Biol. Cell* **33**(13), ar120 (2022).

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, 11/05/2024

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