

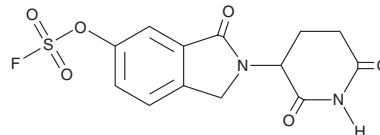
PRODUCT INFORMATION



EM-12-FS

Item No. 41983

CAS Registry No.: 2839138-44-0
Formal Name: 3-[6-[(fluorosulfonyl)oxy]-1,3-dihydro-1-oxo-2H-isoindol-2-yl]-2,6-piperidinedione
MF: C₁₃H₁₁FN₂O₆S
FW: 342.3
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

EM-12-FS is supplied as a solid. A stock solution may be made by dissolving the EM-12-FS in the solvent of choice, which should be purged with an inert gas. EM-12-FS is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

EM-12-FS is a degrader of protein N-terminal glutamine amidohydrolase (NTAQ1), an enzyme involved in N-end protein degradation, and a derivative of (±)-thalidomide (Item No. 14610).¹ It binds to cereblon (CRBN; IC₅₀ = 256 nM) and covalently labels the histidine in position 353 (His³⁵³). EM-12-FS (1 and 5 μM) reduces the levels of NTAQ1 in MOLT-4 cells in a NEDDylation-dependent manner.

Reference

1. Cruite, J.T., Dann, G.P., Che, J., et al. Cereblon covalent modulation through structure-based design of histidine targeting chemical probes. *RSC Chem. Biol.* **3(9)**, 1105-1110 (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, 10/31/2024

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