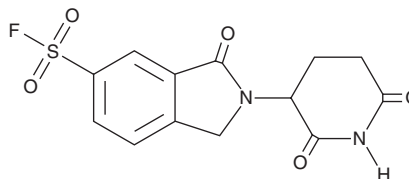


# PRODUCT INFORMATION



**EM12-SO<sub>2</sub>F**  
Item No. 41446

**CAS Registry No.:** 2803819-61-4  
**Formal Name:** 2-(2,6-dioxo-3-piperidiny)-2,3-dihydro-3-oxo-1H-isindole-5-sulfonyl fluoride  
**MF:** C<sub>13</sub>H<sub>11</sub>FN<sub>2</sub>O<sub>5</sub>S  
**FW:** 326.3  
**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

EM12-SO<sub>2</sub>F is supplied as a solid. A stock solution may be made by dissolving the EM12-SO<sub>2</sub>F in the solvent of choice, which should be purged with an inert gas. EM12-SO<sub>2</sub>F is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

EM12-SO<sub>2</sub>F is a covalent inhibitor of cereblon (CRBN; IC<sub>50</sub> = 0.88 nM).<sup>1</sup> It inhibits the degradation of Ikaros family zinc finger protein 1 (IKZF1) induced by the thalidomide derivative lenalidomide (Item No. 14643) in MOLT-4 lymphoblasts when used at a concentration of 1 μM.

## 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, 09/10/2024

## 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](http://WWW.CAYMANCHEM.COM)