

# PRODUCT INFORMATION



**CC-92480**

Item No. 31065

**CAS Registry No.:** 2259648-80-9  
**Formal Name:** 4-[4-[[[2-[(3S)-2,6-dioxo-3-piperidiny]-2,3-dihydro-1-oxo-1H-isoindol-4-yl]oxy)methyl]phenyl]methyl-1-piperaziny]-3-fluoro-benzonitrile

**MF:** C<sub>32</sub>H<sub>30</sub>FN<sub>5</sub>O<sub>4</sub>

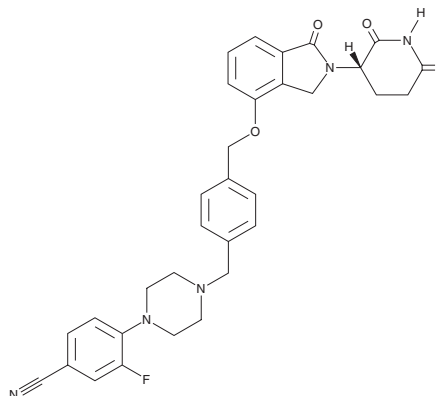
**FW:** 567.6

**Purity:** ≥98%

**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

CC-92480 is supplied as a solid. A stock solution may be made by dissolving the CC-92480 in the solvent of choice, which should be purged with an inert gas. CC-92480 is soluble in the organic solvent DMSO.

## Description

CC-92480 is a modulator of the E3 ligase cereblon.<sup>1</sup> It binds to cereblon (IC<sub>50</sub> = 0.03 μM in a FRET assay) and induces the degradation of Ikaros and Aiolos in lenalidomide-resistant H929 R10-1 cells when used at concentrations of 1, 10, and 100 nM. CC-92480 induces apoptosis in the same cells. It reduces tumor growth in an H929 R10-1 mouse xenograft model when administered at doses of 1, 3, and 10 mg/kg.

## Reference

1. Hansen, J.D., Correa, M., Nagy, M.A., *et al.* Discovery of CRBN E3 ligase modulator CC-92480 for the treatment of relapsed and refractory multiple myeloma. *J. Med. Chem.* **63**(13), 6648-6676 (2020).

### 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, 11/04/2022

## 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