

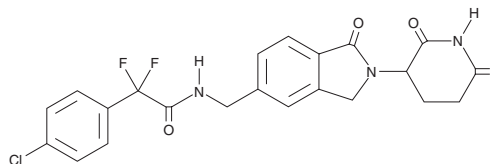
PRODUCT INFORMATION



CC-90009

Item No. 33692

CAS Registry No.: 1860875-51-9
Formal Name: 4-chloro-N-[[2-(2,6-dioxo-3-piperidiny)-2,3-dihydro-1-oxo-1H-isoindol-5-yl]methyl]- α,α -difluoro-benzeneacetamide
MF: C₂₂H₁₈ClF₂N₃O₄
FW: 461.9
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CC-90009 is supplied as a crystalline solid. A stock solution may be made by dissolving the CC-90009 in the solvent of choice, which should be purged with an inert gas. CC-90009 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CC-90009 in these solvents is approximately 10 and 20 mg/ml, respectively.

Description

CC-90009 is a modulator of the E3 ligase cereblon.¹ It binds to cereblon and enhances the protein-protein interaction between cereblon and the cell cycle regulator and translation termination factor G₁ to S phase transition 1 (GSPT1) to induce ubiquitination and proteasomal degradation of GSPT1. CC-90009 (0.01-10 μ M) inhibits the growth of U937 cells in a concentration-dependent manner, an effect that can be reversed by overexpression of the GSPT1 stabilization mutant GSPT1^{G575N}. It inhibits cell growth in a panel of acute myeloid leukemia (AML) cells, an effect that can be reversed by hyperactivation of the mammalian target of rapamycin (mTOR) pathway or knockout of the interleukin enhancer binding factor 2 (ILF2) and ILF3 heterodimeric complex to decrease production of full-length cereblon. CC-90009 induces intratumor GSPT1 degradation in AML mouse xenograft models and decreases leukemic engraftment of 35 patient-derived AML samples in NOD.SCID mice.

Reference

1. Surka, C., Jin, L., Mbong, M., *et al.* CC-90009, a novel cereblon E3 ligase modulator, targets acute myeloid leukemia blasts and leukemia stem cells. *Blood* **137**(5), 661-677 (2021).

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, 03/15/2021

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