

PRODUCT INFORMATION



CC-885

Item No. 19966

CAS Registry No.: 1010100-07-8

Formal Name: N-(3-chloro-4-methylphenyl)-
N'-[[2-(2,6-dioxo-3-piperidiny)-
2,3-dihydro-1-oxo-1H-isoindol-
5-yl)methyl]-urea

MF: C₂₂H₂₁ClN₄O₄

FW: 440.9

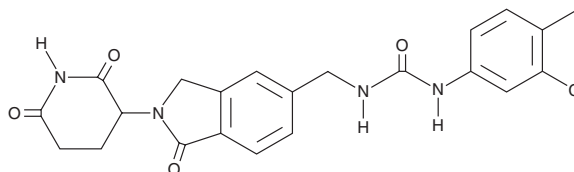
Purity: ≥98%

UV/Vis.: λ_{max}: 244 nm

Supplied as: A crystalline 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-885 is supplied as a crystalline solid. A stock solution may be made by dissolving the CC-885 in the solvent of choice, which should be purged with an inert gas. CC-885 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CC-885 in these solvents is approximately 12.5 mg/ml.

CC-885 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CC-885 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CC-885 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CC-885 is a modulator of the E3 ligase protein cereblon with antiproliferative effects on human myeloid leukemia cell lines.¹ It forms a complex with cereblon and the cell cycle regulator and translation termination factor GSPT1.¹ The antitumour activity of CC-885 relies on cereblon-dependent ubiquitination and degradation of the GSPT1.¹

Reference

1. Matyskiela, M.E., Lu, G., Ito, T., *et al.* A novel cereblon modulator recruits GSPT1 to the CRL4^{CRBN} ubiquitin ligase. *Nature* **535(7611)**, 252-257 (2016).

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, 02/06/2025

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