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-piperidinyl)-2,3-dihydro-1-oxo-1H-isoindol-

5-yl]methyl]-urea

MF: C22H21CIN4O4

FW: 440.9 **Purity:** ≥98%

λ_{max}: 244 nm UV/Vis.: A crystalline solid Supplied as:

-20°C Storage: Stability: ≥2 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. CC-885 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. 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. 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.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the material can be found on our website.

Copyright Cayman Chemical Company, 09/19/2017

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