

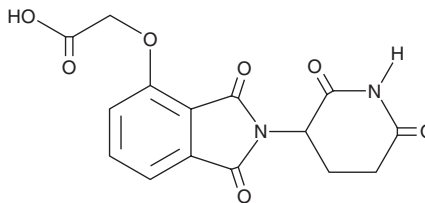
PRODUCT INFORMATION



Cereblon Ligand 1

Item No. 21589

CAS Registry No.: 1061605-21-7
Formal Name: 2-[[2-(2,6-dioxo-3-piperidinyloxy)-2,3-dihydro-1,3-dioxo-1H-indol-4-yl]oxy]-acetic acid
MF: C₁₅H₁₂N₂O₇
FW: 332.3
Purity: ≥95%
UV/Vis.: λ_{max}: 220, 330 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

Cereblon ligand 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the cereblon ligand 1 in the solvent of choice, which should be purged with an inert gas. Cereblon ligand 1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of cereblon ligand 1 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of cereblon ligand 1 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cereblon ligand 1 in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cereblon ligand 1 is a building block in the synthesis of proteolysis-targeting chimera technologies (PROTACs).¹

Reference

1. Lohbeck, J. and Miller, A.K. Practical synthesis of a phthalimide-based Cereblon ligand to enable PROTAC development. *Bioorg. Med. Chem. Lett.* **26(21)**, 5260-5262 (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, 12/22/2022

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