

PRODUCT INFORMATION



CRBN-6-5-5-VHL

Item No. 45748

CAS Registry No.: 2362575-45-7
Formal Name: N-[5-[[[5-[[[2-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4R-yl]amino]hexyl]oxy]pentyl]oxy]-1-oxopentyl]-3-methyl-L-valyl-4-hydroxy-N-[[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide

Synonyms: Cereblon-6-5-5-VHL,
Cereblon-6-5-5-von,
CRBN-6-5-5-von,
Hippel-Lindau

MF: C₅₁H₆₉N₇O₁₀S

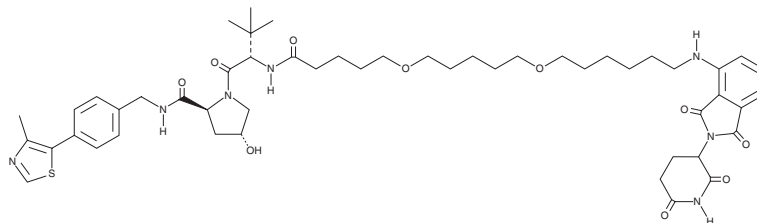
FW: 972.2

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

CRBN-6-5-5-VHL is supplied as a solid. A stock solution may be made by dissolving the CRBN-6-5-5-VHL in the solvent of choice, which should be purged with an inert gas. CRBN-6-5-5-VHL is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Description

CRBN-6-5-5-VHL is a proteolysis-targeting chimera (PROTAC) containing the cereblon (CRBN) inhibitor pomalidomide (Item No. 19877) and the von Hippel-Lindau (VHL) E3 ligase ligand VH 032 (Item No. 39982).¹ It selectively induces the degradation of CRBN over Ikaros family zinc finger protein 1 (IKZF1) in MM.1S cells at 10-500 nM. CRBN-6-5-5-VHL (100 nM) rescues pomalidomide-induced cytotoxicity in MM.1S cells.

Reference

1. Powell, C.E., Du, G., Bushman, J.W., *et al.* Selective degradation-inducing probes for studying cereblon (CRBN) biology. *RSC Med. Chem.* **12**(8), 1381-1391 (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, 06/17/2026

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