

PRODUCT INFORMATION



VL285

Item No. 41538

CAS Registry No.: 1448188-57-5

Formal Name: 1-[(2S)-2-(1,3-dihydro-1-oxo-2H-isoindol-2S-yl)-3-methyl-1-oxobutyl]-4-hydroxy-N-[[4-(4R-methyl-5-thiazolyl)phenyl]methyl]-2-pyrrolidinecarboxamide

MF: C₂₉H₃₂N₄O₄S

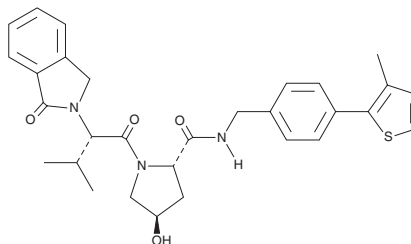
FW: 532.7

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

VL285 is supplied as a solid. A stock solution may be made by dissolving the VL285 in the solvent of choice, which should be purged with an inert gas. VL285 is soluble (≥10 mg/ml) in DMSO. VL285 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

VL285 is a starting material that has been used in the synthesis of proteolysis-targeting chimeras (PROTACs) that drive the degradation of tagged proteins.¹

Reference

1. Buckley, D.L., Raina, K., Darricarrere, N., *et al.* HaloPROTACS: Use of small molecule PROTACs to induce degradation of HaloTag fusion proteins. *ACS Chem. Biol.* **10**(8), 1831-1837 (2015).

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, 10/09/2024

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