

# PRODUCT INFORMATION



## PROTAC CRBN Degrader-1

Item No. 41878

CAS Registry No.: 2358775-70-7

Formal Name: N1-(2-((2-(2,6-dioxopiperidin-3-yl)-1,3-dioxoisindolin-4-yl)amino)ethyl)-N20-((S)-1-((2S,4R)-4-hydroxy-2-((4-(4-methylthiazol-5-yl)benzyl)carbamoyl)pyrrolidin-1-yl)-3,3-dimethyl-1-oxobutan-2-yl)-3,9,12,18-tetraoxaicosanediamide

Synonyms: PROTAC Cereblon Degrader-1, Proteolysis-targeting Chimera Cereblon Degrader-1, Proteolysis-targeting Chimera CRBN Degrader-1

MF: C<sub>53</sub>H<sub>72</sub>N<sub>8</sub>O<sub>13</sub>S

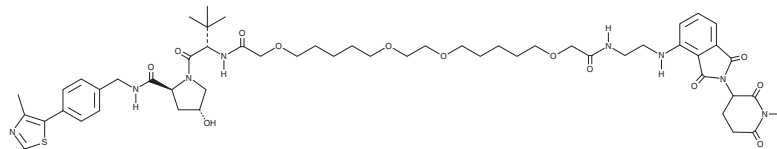
FW: 1,061.3

Purity: ≥95%

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

PROTAC CRBN Degrader-1 is supplied as a solid. A stock solution may be made by dissolving the PROTAC CRBN degrader-1 in the solvent of choice, which should be purged with an inert gas. PROTAC CRBN Degrader-1 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

PROTAC CRBN degrader-1 is a proteolysis-targeting chimera (PROTAC) that contains VHL ligand 1 (Item No. 21591) conjugated to the cereblon (CRBN) inhibitor pomalidomide (Item No. 19877).<sup>1</sup> It induces degradation of CRBN in HeLa cells with a half-maximal degradation concentration (DC<sub>50</sub>) value of 200 nM.

### Reference

1. Girardini, M., Maniaci, C., Hughes, S.J., *et al.* Cereblon versus VHL: Hijacking E3 ligases against each other using PROTACs. *Bioorg. Med. Chem.* **27**(12), 2466-2479 (2019).

#### 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, 09/23/2024

#### 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