

# PRODUCT INFORMATION



## CM11

Item No. 41885

**CAS Registry No.:** 2244684-49-7  
**Formal Name:** (4R,4'R)-1,1'-(1,20-dioxo-3,6,9,12,15,18-hexaoxaicosane-1,20-diy)bis[3-methyl-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide

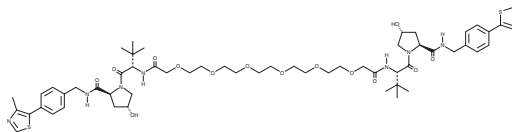
**MF:** C<sub>58</sub>H<sub>82</sub>N<sub>8</sub>O<sub>14</sub>S<sub>2</sub>  
**FW:** 1,178.5

**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

CM11 is supplied as a crystalline solid. A stock solution may be made by dissolving the CM11 in the solvent of choice, which should be purged with an inert gas. CM11 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

CM11 is a homodimeric proteolysis-targeting chimera (PROTAC) composed of two molecules of the von Hippel-Lindau (VHL) E3 ligase ligand VH 032 (Item No. 39982) linked by a PEG linker.<sup>1</sup> It selectively induces the degradation of the long isoform of VHL, pVHL30, over the short isoform, pVHL19, and a panel of 6,449 proteins in HeLa cells at 1 μM but does decrease protein levels of cullin 2 (Cul-2).

### Reference

1. Maniaci, C., Hughes, S.J., Testa, A., *et al.* Homo-PROTACs: Bivalent small-molecule dimerizers of the VHL E3 ubiquitin ligase to induce self-degradation. *Nat. Commun.* **8(1)**, 830 (2017).

#### 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, 05/06/2025

#### 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