

# PRODUCT INFORMATION



## PROTAC Bcl-2 Degradator-1

Item No. 45385

**CAS Registry No.:** 2378801-85-3  
**Formal Name:** N<sup>1</sup>-[2-[6-[(4-bromophenyl)thio]-1,3-dioxo-1H-benz[de]isoquinolin-2(3H)-yl]ethyl]-N<sup>6</sup>-[2-[2-[2-[[2-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]ethoxy]ethoxy]ethyl]-hexanediamide

**Synonyms.:** PROTAC BCL2 Degradator-1, Proteolysis-targeting Chimera Bcl-2 Degradator-1, Proteolysis-targeting Chimera BCL2 Degradator-1

**MF:** C<sub>45</sub>H<sub>45</sub>BrN<sub>6</sub>O<sub>10</sub>S

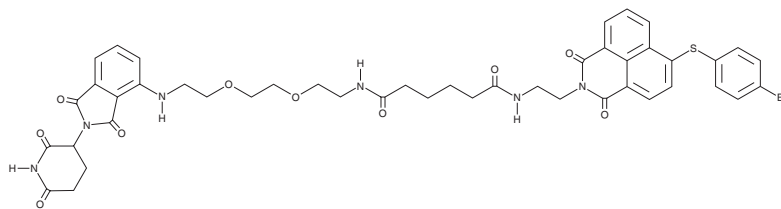
**FW:** 941.9

**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

PROTAC Bcl-2 degrader-1 is supplied as a solid. A stock solution may be made by dissolving the PROTAC Bcl-2 degrader-1 in the solvent of choice, which should be purged with an inert gas. PROTAC Bcl-2 degrader-1 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

PROTAC Bcl-2 degrader-1 is a proteolysis-targeting chimera (PROTAC) composed of a Bcl-2 inhibitor conjugated to the cereblon (CRBN) inhibitor pomalidomide (Item No. 19877) via a PEG linker.<sup>1</sup> It induces the degradation of Bcl-2 in HeLa cells with a half-maximal degradation concentration (DC<sub>50</sub>) value of 3 μM. PROTAC Bcl-2 degrader-1 induces selective degradation of Bcl-2 over myeloid cell leukemia-1 (Mcl-1), Bcl-xL, murine double minute 2 (MDM2), and tubulin in HeLa cells when used at a concentration of 10 μM.

### Reference

1. Wang, Z., He, N., Guo, Z., *et al.* Proteolysis targeting chimeras for the selective degradation of Mcl-1/Bcl-2 derived from nonselective target binding ligands. *J. Med. Chem.* **62**(17), 8152-8163 (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, 04/17/2026

#### 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