

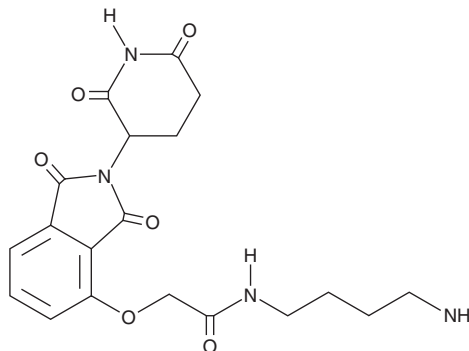
# PRODUCT INFORMATION



## E3 Ligand-Linker Conjugate 3

Item No. 36666

**CAS Registry No.:** 1799711-24-2  
**Formal Name:** N-(4-aminobutyl)-2-[[2-(2,6-dioxo-3-piperidiny)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]oxy]-acetamide  
**MF:** C<sub>19</sub>H<sub>22</sub>N<sub>4</sub>O<sub>6</sub>  
**FW:** 402.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 218 nm  
**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

E3 ligand-linker conjugate 3 is supplied as a solid. Aqueous solutions of E3 ligand-linker conjugate 3 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of E3 ligand-linker conjugate 3 in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

E3 ligand-linker conjugate 3 is a building block in the synthesis of proteolysis-targeting chimera (PROTAC) technologies containing thalidomide (Item No. 14610), which is a ligand for the E3 ubiquitin ligase cereblon.<sup>1,2</sup> It has been used in the synthesis of PROTACs for the degradation of several proteins, including anaplastic lymphoma kinase (ALK) and the bromodomain and extra terminal domain (BET) family of proteins, with anticancer activity.

### References

1. Yan, G., Zhong, X., Yue, L., *et al.* Discovery of a PROTAC targeting ALK with *in vivo* activity. *Eur. J. Med. Chem.* **212(1)**, 113150 (2021).
2. Zhou, B., Hu, J., Xu, F., *et al.* Discovery of a small-molecule degrader of bromodomain and extra-terminal (BET) proteins with picomolar cellular potencies and capable of achieving tumor regression. *J. Med. Chem.* **61(2)**, 462-481 (2018).

#### 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, 11/14/2022

#### 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