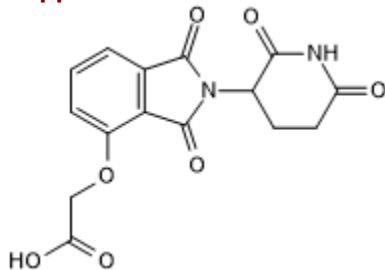


LS E3 1017
Cat. # PC-1008

Background: Derivative of Thalidomide (Cat.No. PC-1006), commonly used as a precursor to PROTACs that hijack cereblon as the E3 ubiquitin ligase component. Supplied with a carboxylic acid functional handle at a position known not to significantly affect binding to cereblon, for ready conjugation to a linker/target protein ligand.

Application: Functionalized cereblon ligand for PROTACs

**Product Information**

Purity:	>98%
MW:	332.26
Formula:	$C_{15}H_{12}N_2O_7$
CAS Number	1061605-21-7
Physical State:	Lyophilized off white solid
Quantity:	5mg; 10mg; 25mg
Solubility:	30 mg/mL in DMSO
Storage:	Store at room temperature.

References

1. **Lohbeck and Miller** (2016) Practical synthesis of a phthalimide-based cereblon ligand to enable PROTAC development. *Bioorg.Med.Chem.Lett.* **26** 5260 PMID: [27687673](https://pubmed.ncbi.nlm.nih.gov/27687673/)

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved