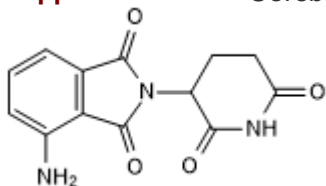


Pomalidomide

Cat. # PC-1005

Background: Potent TNF- α inhibitor (IC_{50} = 13 nM). Also potently inhibits IL-2 (EC_{50} = 8 nM). Thalidomide derivative. Binds cereblon and inhibits its ubiquitination. Exhibits antiproliferative effects in a Namalwa lymphoma cell line. Antiangiogenic.

Application: Cereblon ubiquitination inhibitor; also TNF- α inhibitor and antiangiogenic



Product Information

Purity:	>95%
MW:	273.24
Formula:	C₁₃H₁₁N₃O₄
CAS Number	19171-19-8
Physical State:	Lyophilized pale-yellow to yellow-brown solid
Quantity:	1g; 5g
Solubility:	25 mg/mL in DMSO
Storage:	Room Temperature

References

1. **Lohbeck & 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/)
2. **Nanthakumar et al** (2015) Dissecting fibrosis: therapeutic insights from the small-molecule toolbox. *Nat.Rev.Drug.Discov.* **14** 693 PMID: [26338155](https://pubmed.ncbi.nlm.nih.gov/26338155/)
3. **Lai et al** (2016) Modular PROTAC design for the degradation of oncogenic BCR-ABL. *Angew.Chem.Int.Ed.Engl.* **55** 807 PMID: [26593377](https://pubmed.ncbi.nlm.nih.gov/26593377/)
4. **Winter et al** (2015) Phthalimide conjugation as a strategy for *in vivo* target protein degradation. *Science.* **348**1376 PMID: [25999370](https://pubmed.ncbi.nlm.nih.gov/25999370/)

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