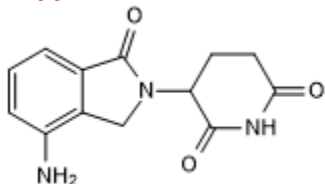


Lenalidomide

Cat. # PC-1004

Background: Thalidomide analog. Immune modulatory drug and cereblon binding compound. Induces ubiquitination and degradation of casein kinase (CK) 1 α by the E3 ubiquitin ligase CRL4^{CRBN}. Also TNF- α inhibitor and angiogenesis inhibitor.

Application: Cereblon binder; induces ubiquitination and degradation of CK1 α by E3 ubiquitin ligase



Product Information

Purity:	>97%
MW:	259.26
Formula:	C₁₃H₁₃N₃O₃
CAS Number	191732-72-6
Physical State:	Lyophilized white to yellow solid
Quantity:	1g; 5g
Solubility:	25 mg/mL in DMSO
Storage:	Room Temperature

References

1. **Petzold *et al*** (2016) Structural basis of lenalidomide-induced CK1 α degradation by the CRL4(CRBN) ubiquitin ligase. *Nature* **532** 127 PMID: [26909574](https://pubmed.ncbi.nlm.nih.gov/26909574/)
2. **Krönke *et al*** (2015) Lenalidomide induces ubiquitination and degradation of CK1 α in del(5q) MDS. *Nature* **523** 183 PMID: [26131937](https://pubmed.ncbi.nlm.nih.gov/26131937/)

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