

## Lenalidomide

Thalidomide analog. Immune modulatory drug (IMiD) and cereblon binding compound. Induces ubiquitination and degradation of neosubstrates such as IKZF1/3 and casein kinase (CK) 1 $\alpha$  by the E3 ubiquitin ligase CRL4 (CRBN).

### Materials Provided

<b>Catalog Number:</b>	<b>92-1361</b>
<b>Quantity Supplied:</b>	1 vial(s), 100 mg per vial

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

<b>Molecular Weight:</b>	259.26
<b>Source:</b>	Synthetic
<b>Purity:</b>	>98% (HPLC)
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	-20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	100 mM
<b>Reconstitution Volume:</b>	3.86 mL
<b>Reconstitution Solvent:</b>	DMSO

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)