

## Product Information

---

<b>Description</b>	4-Hydroxy-thalidomide is the Thalidomide-based Cereblon ligand used in the recruitment of CRBN protein. It can be connected to the ligand for protein by a linker to form PROTACs.
<b>Molecular Formula:</b>	C <sub>13</sub> H <sub>10</sub> N <sub>2</sub> O <sub>5</sub>
<b>Supplied As:</b>	Solid
<b>Stability:</b>	At least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>MW:</b>	274.2 Da
<b>Cas No.</b>	5054-59-1
<b>Purity:</b>	98%

## Chemical Structure

---

