

## Product Information

<b>Description</b>	4-Fluoro-thalidomide is the Thalidomide-based Cereblon ligand used in the recruitment of CRBN protein. It can be connected to the ligand for IRAK4 protein by a linker to form PROTAC IRAK4 degrader-1 (HY-129966).
<b>Molecular Formula:</b>	C <sub>13</sub> H <sub>9</sub> FN <sub>2</sub> O <sub>4</sub>
<b>Supplied As:</b>	Solid
<b>Solubility:</b>	Water, DMF, DCM, DMSO
<b>Stability:</b>	At least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>MW:</b>	276.2 Da
<b>Cas No.</b>	835616-60-9
<b>Purity:</b>	98%

## Chemical Structure

