

## Product Information

<b>Description</b>	Thalidomide-5,6-F is used in the recruitment of the cereblon protein and also used as a PROTAC linker.
<b>Molecular Formula:</b>	C <sub>13</sub> H <sub>8</sub> F <sub>2</sub> N <sub>2</sub> O <sub>4</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>	294.2 Da
<b>Cas No.</b>	1496997-41-1

## Chemical Structure

