

## Product Information

<b>Description</b>	3-(5-Bromo-1-oxo-2-isoindolinyl)piperidine-2,6-dione is a lenalidomide analog that can be useful in PROTAC research because it is applicable to the recruitment of CRBN protein.
<b>Molecular Formula:</b>	C <sub>13</sub> H <sub>11</sub> BrN <sub>2</sub> O <sub>3</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>	323.1 Da
<b>Cas No.</b>	1010100-26-1

## Chemical Structure

