

# 3-(5-(Aminomethyl)-1-oxoisindolin-2-yl) piperidine-2,6-dione hydrochloride

#82006  
Lot: 231106

## Product Information

<b>Description</b>	3-(5-(Aminomethyl)-1-oxoisindolin-2-yl)piperidine-2,6-dione hydrochloride is used in the recruitment of the cereblon protein and used as a PROTAC linker.
<b>Molecular Formula:</b>	C <sub>14</sub> H <sub>16</sub> ClN <sub>3</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>	309.8 Da
<b>Cas No.</b>	1158264-69-7

## Chemical Structure

