

## Product Information

<b>Description</b>	LCL-161 is a small molecule Smac mimetic that inhibits multiple inhibitors of apoptosis (IAP) family proteins.
<b>Molecular Formula:</b>	C <sub>26</sub> H <sub>33</sub> FN <sub>4</sub> O <sub>3</sub> S
<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>	500.9 Da
<b>Cas No.</b>	1005342-46-0

## Chemical Structure

