

## Product Information

<b>Description</b>	(S,R,S)-AHPC-C6-NH2 is a E3 ligase ligand-linker conjugate that is comprised of the (S,R,S)-AHPC based VHL ligand and a six carbon alkane linker used for EED-Targeted PROTAC.
<b>Molecular Formula:</b>	C <sub>29</sub> H <sub>43</sub> N <sub>5</sub> O <sub>4</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>	557.8 Da
<b>Cas No.</b>	2360522-76-3

## Chemical Structure

