

Product Information

Description	(S,R,S)-AHPC-C4-NH2 is a E3 ligase ligand-linker conjugate that is comprised of the (S,R,S)-AHPC based VHL ligand and a four carbon alkane linker used for EED-Targeted PROTAC.
Molecular Formula:	C ₂₇ H ₃₉ N ₅ O ₄ S
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	529.7 Da
Cas No.	2138439-53-7

Chemical Structure

