

Product Information

Description	(S,R,S)-AHPC-Me-C10-NH2 is a E3 ligase ligand-linker conjugate that couples the VHL ligand to a 10 carbon alkane linker. (S,R,S)-AHPC-Me-C10-NH2 can be used in PROTAC MS432.
Molecular Formula:	C ₃₄ H ₅₃ N ₅ O ₄ S
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	627.9 Da
Cas No.	2376139-52-3

Chemical Structure

