

Product Information

Description	(S,R,S)-AHPC-amido-C6-acid is a VHL ligand with a 6 carbon alkyl linker and terminal acid for PROTAC compounds.
Molecular Formula:	C ₃₀ H ₄₂ N ₄ O ₆ S
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	586.8 Da
Cas No.	2172819-75-7

Chemical Structure

