

Product Information

Description	(S,R,S)-AHPC-BOC is a ligand used to recruit the von Hippel-Lindau protein.
Molecular Formula:	C ₂₇ H ₃₈ N ₄ O ₅ S
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	530.7 Da
Cas No.	1448189-98-7

Chemical Structure

