

Product Information

Description	Homo-PROTAC Cereblon Degrader 1 is a highly potent and efficient Cereblon (CRBN) degrader with minimal effects on IKZF1 and IKZF3.
Molecular Formula:	$C_{32}H_{32}N_6O_{10}$
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	660.6 Da
Cas No.	2244520-98-5

Chemical Structure

