

Product Information

Description	E3 Ligase ligand 9 targets ubiquitin-protein ligases in order to activate the E3 ubiquitin ligase. It can then treat metabolic disorders and psychiatric disorders.
Molecular Formula:	$C_{21}H_{32}N_2O_6$
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	408.5 Da
Cas No.	87304-15-2

Chemical Structure

