

Product Information

Description	N-(2-(2-Aminoethoxy)ethyl)-2-((2-(2,6-dioxopiperidin-3-yl)-1,3-dioxoisindolin-4-yl)oxy)acetamide is an E3 ligase ligand-linker conjugate used in PROTAC dTAG-13 as a degrader of FKBP12 F36V and BET.
Molecular Formula:	C ₁₉ H ₂₂ N ₄ O ₇
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	418.4 Da
Cas No.	2022182-59-6

Chemical Structure

