

Product Information

Description	Pomalidomide 4'-PEG3-azide is an E3 ligase ligand-linker that incorporates the Pomalidomide-based cereblon ligand and a linker.
Molecular Formula:	C ₂₁ H ₂₆ N ₆ O ₇
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	474.5 Da
Cas No.	2271036-46-3
Purity:	95%

Chemical Structure

