

Product Information

Description	Pomalidomide-PEG3-NH-Boc is synthesized intermediate that incorporates the Pomalidomide based cereblon ligand and 4-unit PEG linker used in PROTAC technology.
Molecular Formula:	C ₂₆ H ₃₆ N ₄ O ₉
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	548.6 Da
Cas No.	2204246-03-5
Purity:	98%

Chemical Structure

