

Product Information

Description	Thalidomide-O-amido-PEG4-propargyl is a compound that incorporates the Thalidomide based cereblon ligand and PEG4 linker used in PROTAC technology.
Molecular Formula:	C ₂₆ H ₃₀ N ₂ O ₁₁
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	546.5 Da
Cas No.	2496687-00-2
Purity:	98%

Chemical Structure

