

Product Information

Description	Thalidomide-Acid is a linker of Thalidomide which is commonly used as a precursor to PROTAC that hijacks cereblon as the E3 ubiquitin ligase component.
Molecular Formula:	C ₁₅ H ₁₂ N ₂ O ₇
Supplied As:	Solid
Solubility:	DCM
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	332.3 Da
Cas No.	1061605-21-7
Purity:	98%

Chemical Structure

