

## Product Information

**Description**

E3 ligase ligand-linker conjugate, Thalidomide-O-PEG6-Acid, is synthesized compound that incorporates the Thalidomide based cereblon ligand and the PEG linker used in PROTAC technology.

**Molecular Formula:**

$C_{28}H_{38}N_2O_{13}$

**Supplied As:**

Solid

**Solubility:**

DMSO, DCM, DMF

**Stability:**

At least 6 months at -20°C. Avoid freeze/thaw cycles.

**Storage:**

-20°C

**MW:**

610.6 Da

**Purity:**

98%

## Chemical Structure

