

Product Information

Description

Thalidomide-O-PEG4-Propargyl is synthesized compound that incorporates the Thalidomide based cereblon ligand and 4-unit PEG linker used in PROTAC technology, Propargyl is reactive with azide molecule in the presence of Cu catalyst.

Molecular Formula:C₂₄H₂₈N₂O₉**Supplied As:**

Solid

Solubility:

Water, DMF, DCM, DMSO

Stability:

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

Storage:

-20°C

MW:

488.5 Da

Cas No.

2098799-77-8

Purity:

98%

Chemical Structure

