

Product Information

Description

E3 ligase ligand-linker conjugate, Thalidomide-O-PEG2-NHS ester, is synthesized compound that incorporates the Thalidomide based cereblon ligand and the PEG linker used in PROTAC technology.

Molecular Formula:

C₂₄H₂₅N₃O₁₁

Supplied As:

Solid

Solubility:

DMSO, DCM, DMF

Stability:

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

Storage:

-20°C

MW:

531.5 Da

Cas No.

2639395-34-7

Purity:

98%

Chemical Structure

