

Product Information

Description

Thalidomide-O-Hexyl-Acid is an E3 ligase ligand-linker conjugate that incorporates the Thalidomide based cereblon ligand and an alkane linker used in PROTAC technology.

Molecular Formula:

C₂₀H₂₂N₂O₇

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

402.4 Da

Cas No.

2169266-69-5

Chemical Structure

