

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:

$C_{19}H_{24}ClN_3O_5$

Supplied As:

Solid

Stability:

At least 6 months at $-20^{\circ}C$. Avoid freeze/thaw cycles.

Storage:

$-20^{\circ}C$

MW:

409.9 Da

Cas No.

2245697-88-3

Chemical Structure

