

Product Information

Description

(S,R,S)-AHPC-PEG6-NHS ester is comprised of VHL Ligand coupled to an NHS ester via a PEG6 linker. The 1 VHL Ligand 1 is a synthetic compound used in PROTAC research based on E3 ligase ligand. The NHS ester can be used to label the primary amines (-NH₂) of proteins, amine-modified oligonucleotides, and other amine-containing molecules.

Molecular Formula:

C₄₂H₆₁N₅O₁₄S

Supplied As:

Solid

Solubility:

DMSO, DCM, DMF

Stability:

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

Storage:

-20°C

MW:

892.0 Da

Chemical Structure

