

Product Information

Description

(S,R,S)-AHPC-PEG3-amine hydrochloride salt is comprised of VHL Ligand coupled to an amine via a PEG3 linker. The 1 VHL Ligand 1 is a synthetic compound used in PROTAC research based on E3 ligase ligand. NH₂ group is reactive with carboxylic acids in the presence of coupling reagent such as EDC, or directly with NHS esters, to form a stable amide bond.

Molecular Formula:

C₃₀H₄₅N₅O₇S

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

619.8 Da

Cas No.

2097971-11-2

Chemical Structure

