

Product Information

Description

(S,R,S)-AHPC-PEG2-acid is a synthetic PROTAC linker molecule that incorporates an E3 ligase ligand with a PEG linker to empower PROTAC drug research & discovery. The hydrophilic PEG spacer increases the solubility of a compound in aqueous media.

Molecular Formula:

C₃₀H₄₂N₄O₈S

Supplied As:

Solid

Solubility:

Water, DMSO, DMF, Methanol

Stability:

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

Storage:

-20°C

MW:

618.8 Da

Cas No.

2172820-09-4

Purity:

98%

Chemical Structure

