

Product Information

Description

(S,R,S)-AHPC-PEG2-Tos is a combination von Hippel-Lindau (VHL)-recruiting ligand and a PEGylated crosslinker with a tosyl group. The tosyl group is a good leaving group for nucleophilic substitution reactions. (S,R,S)-AHPC-PEG2-Tos is useful for synthesis of PROTAC compounds and molecules for targeted protein degradation.

Molecular Formula:

C₃₄H₄₄N₄O₈S₂

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

700.2 Da

Purity:

98%

Chemical Structure

