

Product Information

Description

E3 Ligase ligand 1a is the (S,R,S)-AHPC-based VHL ligand used in the recruitment of the von Hippel-Lindau (VHL) protein. E3 Ligase ligand 1a can be used to synthesize ARV-771, a BET PROTAC degrader.

Molecular Formula:

C₂₃H₃₂N₄O₃S

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

444.6 Da

Cas No.

1948273-02-6

Chemical Structure

