

Product Information

Description

VHL Ligand 1 HCl salt is a synthetic E3 ligase ligand that is useful for PROTAC drug research & discovery. VHL Ligand 1 works by binding the Hippel-Lindau (VHL) E3 ligase portion of VHL-based PROTACs.

Molecular Formula:

C₂₂H₃₀N₄O₃S

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

430.6 Da

Cas No.

1448189-80-7

Purity:

98%

Chemical Structure

