

Product Information

Description

LC-2 is a von Hippel-Lindau-based PROTAC which can degrade endogenous KRAS G12C. It covalently binds KRAS G12C with a MARTX849 warhead and recruits the E3 ligase VHL which degrades the KRAS G12C and suppresses MAPK signaling in homozygous and heterozygous KRAS G12C cell lines.

Molecular Formula:

C₅₉H₇₁ClFN₁₁O₇S

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

1132.8 Da

Cas No.

2502156-03-6

Chemical Structure

