

Product Information

Description

CC-90009 is a modulator of the E3 ligase cereblon, binding to cereblon and leading to the induction of ubiquitination and proteasomal degradation of GSPT1.

Molecular Formula:

C₂₂H₁₈ClF₂N₃O₄

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

461.9 Da

Cas No.

1860875-51-9

Chemical Structure

