

Product Information

Description

THAL-SNS-032 is a selective CDK9 degrader PROTAC composed of a CDK-binding SNS-032 ligand linked to a thalidomide derivative that binds the E3 ubiquitin ligase Cereblon (CRBN).

Molecular Formula:

C₄₀H₅₂N₈O₁₀S₂

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

869.0 Da

Cas No.

2139287-33-3

Chemical Structure

