

Product Information

Description

ARV-771 is a BET PROTAC based on E3 ligase von Hippel-Lindau with Kds of 34 nM, 4.7 nM, 8.3 nM, 7.6 nM, 9.6 nM, and 7.6 nM for BRD2(1), BRD2(2), BRD3(1), BRD3(2), BRD4(1) and BRD4(2), respectively.

Molecular Formula:

C₄₉H₆₀ClN₉O₇S₂

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

986.7 Da

Cas No.

1949837-12-0

Chemical Structure

