

Product Information

Description

MZ1 is a PROTAC compound that is selective for proteasomal degradation of bromodomain-containing protein 4 (BRD4) in HeLa cells. MZ1 contains JQ-1, which binds bromo- and extra-terminal (BET) proteins, linked to a ligand for the E3 ubiquitin ligase VHL.

Molecular Formula:

$C_{49}H_{60}ClN_9O_8S_2$

Supplied As:

Solid

Stability:

At least 6 months at $-20^{\circ}C$. Avoid freeze/thaw cycles.

Storage:

$-20^{\circ}C$

MW:

1002.7 Da

Cas No.

1797406-69-9

Chemical Structure

