

Product Information

Description

I-BRD9 is a selective cellular chemical probe for bromodomain-containing protein 9 (BRD9). I-BRD9 has a pIC_{50} value of 7.3 for BRD9 inhibitor, with greater than 700-fold selectivity over the BET family and 200-fold over the highly homologous bromodomain-containing protein 7 (BRD7). I-BRD9 is also shows a greater than 70-fold selectivity against a panel of 34 bromodomains. I-BRD9 downregulates CLEC1, DUSP6, FES and SAMS1 genes in Kasumi-1 cells.

Molecular Formula:

$C_{22}H_{22}F_3N_3O_3S_2$

Supplied As:

Solid

Stability:

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

Storage:

$-20^{\circ}C$

MW:

497.6 Da

Cas No.

1714146-59-4

Chemical Structure

