

Product Information

Description

PFI-3 is a selective chemical probe for SMARCA bromodomains that inhibits SMARCA2, SMARCA4 and PB1(5) bromodomains. PFI-3 is a probe that binds avidly to the structurally-similar SMARCA4 bromodomain and PB1(bromodomain 5) with K_d values of 89 and 48 nM, respectively. PFI-3 displays 30-fold selectivity over other sub-family branches and accelerates FRAP recovery in cells at a concentration of 1 μ M.

Molecular Formula:

$C_{19}H_{19}N_3O_2$

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

321.4 Da

Cas No.

1819363-80-8

Chemical Structure

