

Product Information

Description

NI-57 is a selective and potent inhibitor of BRPF (Bromodomain and PHD Finger) family of proteins (BRPF1/2/3). NI-57 shows K_d values of 31, 108 and 408 nM, for BRPF1B, 2 and 3, respectively. NI-57 also exhibits >32-fold selectivity for BRPFs over BRD9 and other non-Class IV bromodomains. Inhibits RANKL-induced differentiation of primary murine bone marrow cells and human primary monocytes into osteoclasts.

Molecular Formula:C₁₉H₁₇N₃O₄**Supplied As:**

Solid

Stability:

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

Storage:

-20°C

MW:

383.4 Da

Cas No.

1883548-89-7

Chemical Structure

