

## Product Information

**Description**

GSK-5959 exhibits >100-fold selectivity for BRPF1 ( $IC_{50} = 80$  nM) over a panel of 35 other bromodomains, including BRPF2/3 and BET family bromodomains. Inhibits BRPF1 interaction with histone H3.3, which is involved in the epigenetic regulation of gene expression and have been implicated in human cancer.

**Molecular Formula:** $C_{22}H_{26}N_4O_3$ **Supplied As:**

Solid

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

394.5 Da

**Cas No.**

901245-65-6

## Chemical Structure

