

Product Information

Description

GSK-6853 is a potent and selective BRPF1 inhibitor. GSK-6853 exhibits >1600-fold selectivity for BRPF1 over other bromodomains (pIC₅₀ values by TR-FRET are 8.1, 5.1/4.8 and 4.7/<4.3 for BRPF1, BRPF2/3 and BRD4 BD1/2, respectively). The structural analog of GSK-6853, GSK-9311, serves as a good inactive control (125- and 185-fold reduced potency in cell-free and cell-based assays, respectively).

Molecular Formula:C₂₂H₂₇N₅O₃**Supplied As:**

Solid

Stability:

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

Storage:

-20°C

MW:

409.5 Da

Cas No.

1910124-24-1

Chemical Structure

