

Product Information

Description

BAY-299 is a potent and selective inhibitor of BRD1 and the second bromodomain of TAF1 (Transcription initiation factor TFIID subunits. BAY-299 has IC₅₀ values of 6-67 and 8-13 nM for BRD1 and TAF1 inhibitor, respectively. BAY-299 shows selectivity over other bromodomains (>30-fold over other members of the BRPF family; BRD9 and ATAD2; >300-fold over BRD4).

Molecular Formula:C₂₅H₂₃N₃O₄**Supplied As:**

Solid

Stability:

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

Storage:

-20°C

MW:

429.5 Da

Cas No.

2080306-23-4

Chemical Structure

