

(+)-JQ1

Catalog #: 27402

Lot #: 140627

Size: 5 mg

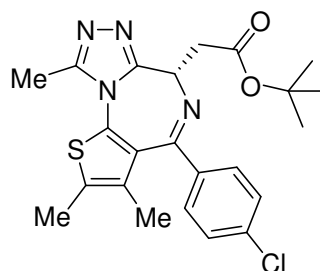
Structure:

CAS Registry #: N/A

Purity: ≥ 95%

Chemical Formula: C₂₃H₂₅ClN₄O₂S

Molecular Weight: 456.99



Description: (+)-JQ1 is a specific inhibitor of the BET family of acetyl-lysine recognition motifs, including the BRD2, BRD3, BRD4 and BRDT bromodomains.

Appearance: A crystalline solid.

Solubility: Soluble in DMSO or ethanol; insoluble in water.

Biological Activity: (+)-JQ1 has been shown to inhibit BRD4 (BD1) with a K_d of 50 nM, and BRD4 (BD2) with a K_d of 90 nM.

Storage/Stability: Store at or below +4 °C. For DMSO stock solutions, store at -20 °C and avoid multiple freeze-thaw cycles to prevent decomposition.

Quality Control: The purity was determined by HPLC.

Reference(s):

Filippakopoulos P., *et al. Nature*. **468(7327)**:1067-73 (2011).

Manufactured, marketed and sold with authorization from Tensha Therapeutics, Inc. Patent Pending relating to PCT Publ. No. WO/2011/143669, and any related U.S. and foreign patents and patent applications.