

RediSolution™ (+)-JQ1

Catalog #: 27403

Lot #: 121101

Size: 1 mg

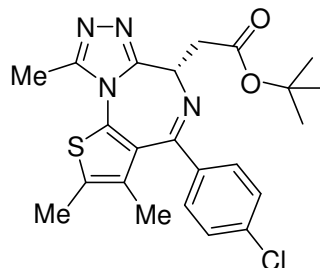
Structure:

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.

Formulation: Provided as a 10 mM solution in anhydrous DMSO

Appearance: Liquid.

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 -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/L/L43669, and any related U.S. and foreign patents and patent applications.