

## (+)-JQ1

**Catalog #:** 27400

**Lot #:** 130118

**Size:** 10mg

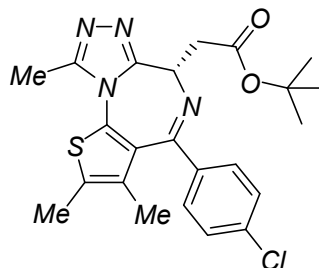
**Structure:**

**CAS Registry #:** N/A

**Purity:** ≥ 95%

**Chemical Formula:** C<sub>23</sub>H<sub>25</sub>ClN<sub>4</sub>O<sub>2</sub>S

**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<sub>d</sub> of 50 nM, and BRD4 (BD2) with a K<sub>d</sub> 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.