

## (-)-JQ1

**Catalog #:** 27334

**Lot #:** 230613

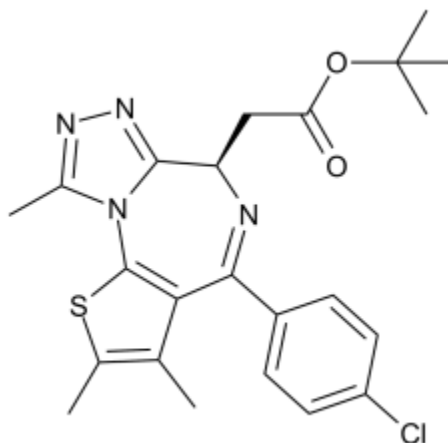
**Size:** 5 mg

**Structure:**

**CAS Registry #:** 1268524-71-5

**Purity:** ≥ 98%

**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 the inactive stereoisomer of the potent, cell-permeable, small molecule bromodomain inhibitor (+)-JQ1 that competitively binds to acetyl-lysine recognition motifs. (+)-JQ1 competitively binds to the bromodomain displacing the BRD4 fusion oncoprotein from chromatin, which induces squamous differentiation and specific anti-proliferative effect in BRD4-dependent cell lines and patient-derived xenograft models. However, study results have shown that (-)-JQ1 fails to significantly interact with any bromodomain tested and exhibits inhibition against BRD4(1) with an IC<sub>50</sub> of 10,000 nM.

**Appearance:** Yellowish, white crystalline powder.

**Solubility:** Soluble in DMSO or ethanol at 100 mM

**Biological Activity:** (-)-JQ1 showed no inhibition of BRD4-dependent NMC cells.

**Storage/Stability:** Store at or below -80°C. Stable for less than 6 months when stored as directed. Avoid freeze/thaw cycles.

**Quality Control:** The purity was determined by HPLC.

### References:

1. Filippakopoulos, P., *et al. Nature*. 2010; **468(7327)**: 1067-1073.