

(-)-JQ1

Catalog #: 27335

Lot #: 230613

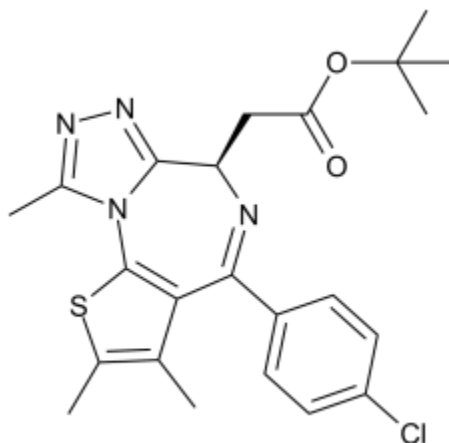
Size: 25 mg

Structure:

CAS Registry #: 1268524-71-5

Purity: ≥ 98%

Chemical Formula: C₂₃H₂₅ClN₄O₂S



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 IC₅₀ 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.