

## Product Information

**Description**

CPI-203 is a primary amide analog of (+)-JQ1, which has shown superior bioavailability with oral or i.p. administration. CPI-203 downregulates Myc expression, causes G1 cell cycle arrest and attenuates cell proliferation in human pancreatic neuroendocrine tumors. CPI-203 also arrests the growth of T cell acute lymphoblastic leukemia cells in vitro ( $EC_{50} = 91.2$  nM).

**Molecular Formula:**C<sub>19</sub>H<sub>18</sub>ClN<sub>5</sub>OS**Supplied As:**

Solid

**Stability:**

At least 6 months at -20°C. Avoid freeze/thaw cycles.

**Storage:**

-20°C

**MW:**

399.9 Da

**Cas No.**

1446144-04-2

## Chemical Structure

