

## Product Information

## Description

BI-2536 is a potent, ATP-competitive, well tolerated and effects Polo-like kinases (Plks) as well as a pan-BET (bromodomain and extra-terminal motif) family inhibitor. BI-2536 has an IC<sub>50</sub> of 0.83 and 25 nM with PLK1 and BRD4 inhibitor, respectively. BI-2536 also blocks the activities of Plk2 and Plk3 to a slightly lesser extent with IC<sub>50</sub> of 3.5 nM and 9.0 nM, respectively. BI-2536 suppresses IFNB (encoding IFN-β) gene transcription. It also shows good blood-brain barrier permeability and causes mitotic arrest in glioblastoma-derived neural stem cells but not normal neural stem cells.

## Molecular Formula:

C<sub>28</sub>H<sub>39</sub>N<sub>7</sub>O<sub>3</sub>

## Supplied As:

Solid

## Stability:

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

## Storage:

-20°C

## MW:

521.7 Da

## Cas No.

755038-02-9

## Purity:

95%

## Chemical Structure

