

## Product Information

## Description

BI-9564 inhibits BRD9 and BRD7 bromodomains and exhibits no significant off-target pharmacology against a panel of 324 kinases and 55 GPCRs at concentrations less than 5  $\mu\text{M}$ . BI-9564 is a selective inhibitor of BRD9 and BRD7 bromodomains ( $K_{\text{ds}}$  = 14.1 and 239 nM;  $\text{IC}_{50\text{S}}$  = 75 nM and 3.4  $\mu\text{M}$ , respectively) that demonstrates cellular activity by semi-quantitative FRAP assay with ~90% inhibition of BRD9 and BRD7 at 0.1  $\mu\text{M}$  and 1  $\mu\text{M}$ , respectively. Inhibits proliferation of AML cell lines in vitro and reduces AML tumor xenograft growth in vivo.

## Molecular Formula:

 $\text{C}_{20}\text{H}_{23}\text{N}_3\text{O}_3$ 

## Supplied As:

Solid

## Stability:

At least 6 months at  $-20^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

## Storage:

 $-20^{\circ}\text{C}$ 

## MW:

353.4 Da

## Cas No.

1883429-22-8

## Chemical Structure

