

## Binding kinetics of AZ23 on TRKB (NTRK2)

