

## Binding kinetics of AZ23 on TRKA (NTRK1)

