

## Binding kinetics of bosutinib on TNIK

