

## Binding kinetics of ponatinib on LCK

