

## Binding kinetics of ponatinib on TIE2

