

## Binding kinetics of afatinib on EGFR [L858R]

