

## Binding kinetics of sorafenib on RET

