

Binding kinetics of ibrutinib on BTK [T474S] non-activated

