

Binding kinetics of ibrutinib on BTK [T474I] non-activated

