

Binding kinetics of ibrutinib on BTK[T474I] non activated

BTK[T474I] non activated Ibrutinib
RU

