

Binding kinetics of ibrutinib on BTK[T474S] non activated

BTK[T474S] non activated Ibrutinib

