

Binding kinetics of ibrutinib on BTK[T474I]

BTK[T474I] Ibrutinib

