

Binding kinetics of ibrutinib on BTK[T474S]

BTK[T474S] Ibrutinib
RU

