

Binding kinetics of ibrutinib on BTK[T316A] non activated

BTK[T316A] non activated Ibrutinib

