

Binding kinetics of fenebrutinib on BTK [C481S] non-activated

