

Binding kinetics of fenebrutinib on BTK [C481S]

