

# Binding kinetics of ibrutinib on BTK[C418S]

