

# Binding kinetics of GDC 0853-0 on BTK[C418S] non activated

BTK[C418S] non activated GDC 0853-0

