

# Binding kinetics of fenebrutinib on BTK [T316A] non-activated

