

# Binding kinetics of ponatinib on BTK

