

Binding kinetics of ponatinib on BTK

