

Binding kinetics of ibrutinib on BTK

