

Binding kinetics of ibrutinib on BRK

