

Binding kinetics of LY2228820 on p38 (inactive)

inactive p38 LY2228820

