

Binding kinetics of LY2228820 on p38

