

Binding kinetics of dinaciclib on CDK6/CycD1

