

Binding kinetics of dinaciclib on CDK9/CycT1

