

Binding kinetics of dinaciclib on CDK7/CycH/MAT1

