

Binding kinetics of SNS-032 on CDK2/CycE1

