

Binding kinetics of PHA-793887 on CDK2/CycA2

