

Binding kinetics of PHA-793887 on CDK9/CycK

