

Binding kinetics of PHA-793887 on CDK9/CycK

CDK9/CycK

PHA-793887

RU

