

Binding kinetics of dinaciclib on CDK9/CycT1

CDK9/CycT1

dinaciclib

RU

