

Binding kinetics of staurosporine on CDK9/CycT2

