

# Binding kinetics of staurosporine on CDK8/CycC

