

# Binding kinetics of staurosporine on CDC2 (CDK1)/CycA2

