

# Binding kinetics of dinaciclib on CDK4/CycD1

