

# Binding kinetics of sunitinib on CHK1

