

# Binding kinetics of sunitinib on CHK1

CHK1 Sunitinib

