

# Binding kinetics of bosutinib on ABL

ABL bosutinib



