

# Binding kinetics of bosutinib on BMX

BMX Bosutinib

