

Binding kinetics of imatinib on DDR2

