

Binding kinetics of imatinib on ABL

