

Binding kinetics of dasatinib on FLT3

