

Binding kinetics of dasatinib on FMS

