

Binding kinetics of dasatinib on BMX

