

Binding kinetics of dasatinib on FYN

