

# Binding kinetics of dasatinib on FLT3

FLT3 Dasatinib

