

Binding kinetics of infigratinib on FGFR1 non-activated

