

Binding kinetics of infigratinib on FGFR4

