

# Binding kinetics of crizotinib on FAK

