

# Binding kinetics of crizotinib on ALK

ALK cricotinib

