

Binding kinetics of crizotinib on AXL

