

Binding kinetics of crizotinib on MET

