

Binding kinetics of crizotinib on MET

MET Crizotinib
RU

