

Binding kinetics of crizotinib on AXL

AXL crizotinib
RU

