

Binding kinetics of crizotinib on ITK

ITK Crizotinib

RU

