

Binding kinetics of ponatinib on DDR2

DDR2 ponatinib

