

Binding kinetics of ponatinib on LCK

