

Binding kinetics of sunitinib on IRAK4

