

Binding kinetics of bosutinib on HPK1 (MAP4K1)

