

Binding kinetics of NTRC 0799-0 on IKK α (CHUK) inactive mutant

