

Binding kinetics of NTRC 0799-0 on IKKa

IKKalpha NTRC 0799-0

