

Binding kinetics of staurosporine on RET [V804M] non-activated

