

Binding kinetics of staurosporine on RET [V804M]

