

Binding kinetics of staurosporine on JAK2

