

Binding kinetics of staurosporine on GSK3 β

