

Binding kinetics of staurosporine on SRPK1

