

Binding kinetics of staurosporine on YES (YES1)

