

Binding kinetics of staurosporine on BTN-GSK3B

