

Binding kinetics of staurosporine on MAP3K5

MAP3K5 Staurosporine

