

Binding kinetics of staurosporine on PAK4

PAK4 Staurosporine
RU

