

Binding kinetics of staurosporine on SYK [non activated]

SYK - staurosporine

