

## Binding kinetics of staurosporine on SYK non-activated

