

Binding kinetics of A-770041 on SRC

