

Binding kinetics of idelalisib on PIK3CA [E542K]/PIK3R1

