

Binding kinetics of idelalisib on PIK3CA [E545K]/PIK3R1

