

Binding kinetics of idelalisib on PIK3CA [H1047R]/PIK3R1

