

Binding kinetics of idelalisib on PIK3CA [R88Q]/PIK3R1

