

Binding kinetics of idelalisib on PIK3CA [P539R]/PIK3R1

