

Binding kinetics of idelalisib on PIK3CD/PIK3R1

