

Binding kinetics of pictilisib on PIK3CA/PIK3R1

