

Binding kinetics of GDC0941 on PIK3CA PIK3R1

PIK3CA/PIK3R1 GDC0941

