

Binding kinetics of GDC-0941 on PIK3CG

GDC-0941 on PIK3CG

