

Binding kinetics of apitolisib on PIK3CG

